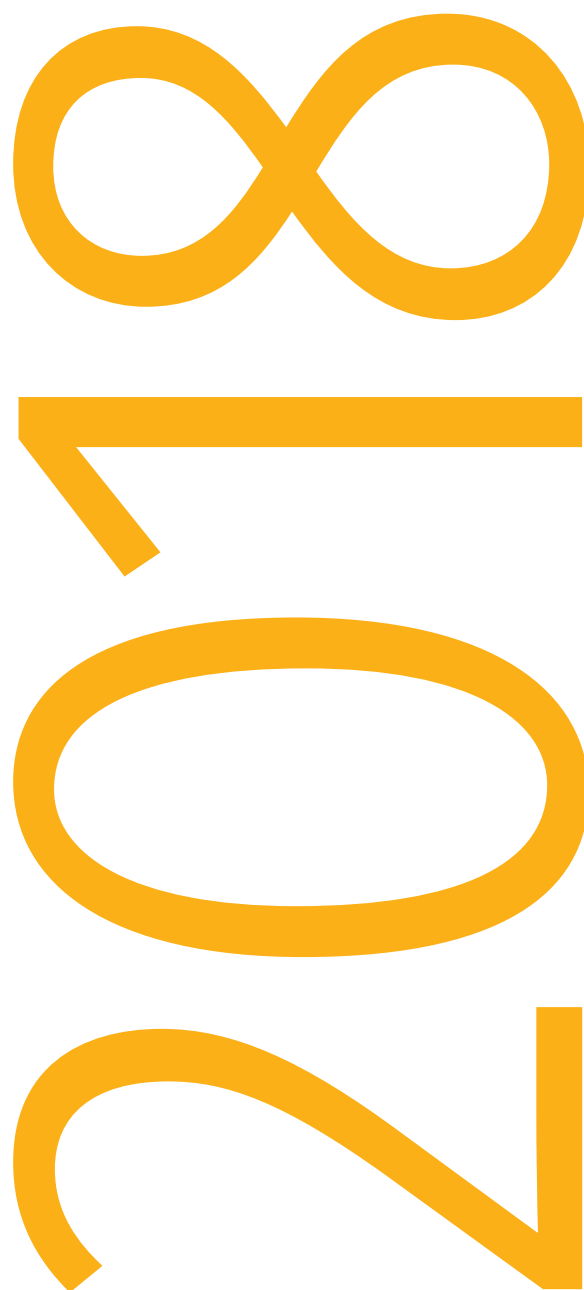




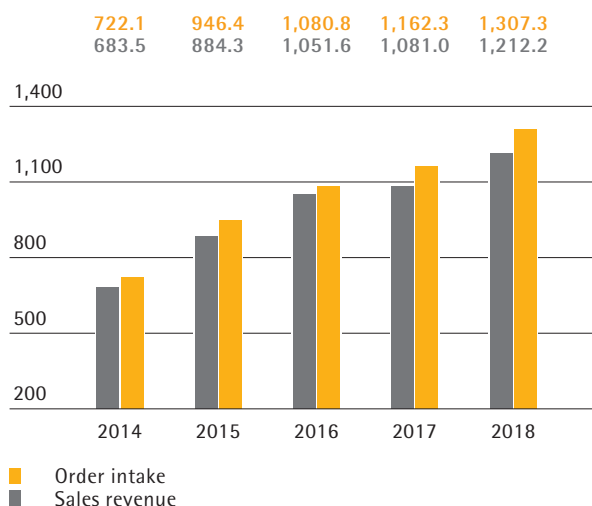
sartorius stedim
biotech

Sartorius Stedim Biotech
Reference Document 2018
Including the Annual Financial Report

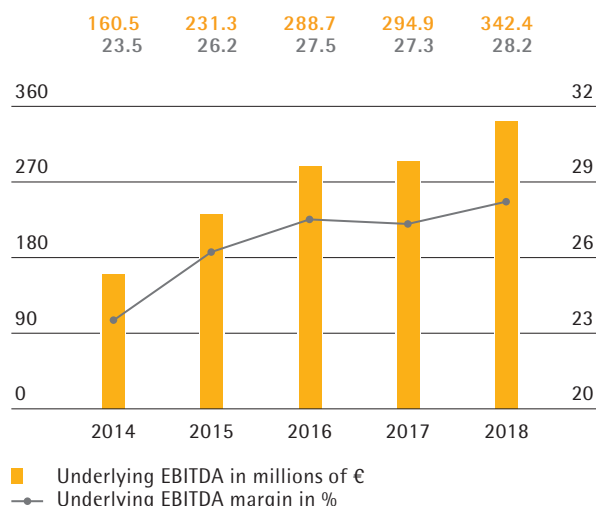


Order Intake and Sales Revenue

€ in millions



Underlying EBITDA and Margin¹⁾



Key Figures

All figures are given in millions of € according to IFRS, unless otherwise specified

| | 2018 | 2017 | 2016 | 2015 | 2014 |
|--|--------------------|---------|---------|---------|--------------------|
| Order intake, sales revenue and earnings | | | | | |
| Order intake | 1,307.3 | 1,162.3 | 1,080.8 | 946.4 | 722.1 |
| Sales revenue | 1,212.2 | 1,081.0 | 1,051.6 | 884.3 | 683.5 |
| Underlying EBITDA ^{1,2)} | 342.4 | 294.9 | 288.7 | 231.3 | 160.5 |
| Underlying EBITDA ^{1,2)} as % of sales revenue | 28.2 | 27.3 | 27.5 | 26.2 | 23.5 |
| Net profit after non-controlling interest | 208.1 | 161.1 | 153.7 | 118.0 | 72.4 |
| Underlying net profit ¹⁾ after non-controlling interest ²⁾ | 219.3 | 180.4 | 176.6 | 139.3 | 87.2 |
| Research and development costs | 60.6 | 53.2 | 47.5 | 41.5 | 34.1 ³⁾ |
| Financial data per share⁴⁾ | | | | | |
| Earnings per share | 2.26 | 1.75 | 1.67 | 1.28 | 0.79 |
| Earnings per share (in €) ^{1,5)} | 2.38 | 1.96 | 1.92 | 1.51 | 0.95 |
| Dividend per share (in €) | 0.57 ⁶⁾ | 0.46 | 0.42 | 0.33 | 0.22 |
| Balance sheet | | | | | |
| Balance sheet total | 1,571.5 | 1,403.9 | 1,195.8 | 1,066.1 | 907.3 |
| Equity | 1,044.9 | 879.5 | 763.6 | 647.2 | 539.1 |
| Equity ratio (in %) | 66.5 | 62.6 | 63.9 | 60.7 | 59.4 |
| Financials | | | | | |
| Capital expenditures | 177.0 | 136.7 | 80.2 | 54.5 | 44.2 |
| Capital expenditures as % of sales revenue | 14.6 | 12.6 | 7.6 | 6.2 | 6.5 |
| Depreciation and amortization | 60.9 | 50.6 | 44.7 | 39.4 | 35.6 |
| Net cash flow from operating activities | 227.3 | 174.7 | 156.7 | 142.8 | 111.3 |
| Net debt ⁷⁾ | 125.7 | 127.1 | 67.6 | 86.4 | 87.4 |
| Ratio of net debt to underlying EBITDA ^{1,2)} | 0.4 | 0.4 | 0.2 | 0.4 | 0.5 |
| Total number of employees as of December 31 | 5,637 | 5,092 | 4,725 | 4,202 | 3,697 |

¹⁾ Adjusted for extraordinary items

²⁾ For more information on EBITDA, net profit and the underlying presentation, please refer to the Group Business Development chapter and to the Glossary.

³⁾ Restated

⁴⁾ 2014 to 2015 adjusted for stock split; rounded values

⁵⁾ Adjusted for extraordinary items, non-cash amortization acc. to IFRS 3 and fair value adjustments of hedging instruments, as well as the corresponding tax effects for each of these items.

⁶⁾ Amount suggested by the Board of Directors (Conseil d'administration) and subject to approval by the Annual General Shareholders' Meeting.

⁷⁾ Net debt excludes the liability for the remaining purchase price for acquisitions; 2018: 8.7 million euros, 2017: 46.5 million euros, 2016: 49.6 million euros, 2015: 47.5 million euros, 2014: 42.8 million euros



sartorius stedim
biotech

Mission & Vision

Mission

We empower engineers to simplify and accelerate progress in bioprocessing. In this way, we enable new and better pharmaceuticals to be manufactured and help keep medications affordable.

Vision

As pioneers, we are a magnet and a dynamic platform for the leading experts in our field. We bring creative minds together for a common goal: technological breakthroughs that lead to better health for more people.



sartorius stedim
biotech

Reference Document 2018



The present original French "Document de Référence" of this translated Reference Document was filed with the Autorité des Marchés Financiers on 18 February, 2019, in accordance with Article 212 - 13 of its "règlement général". It may be used in connection with an offering of securities if it is supplemented by a prospectus ("note d'opération") for which the Autorité des Marchés Financiers has issued an endorsement. This Reference Document has been made out by the issuer and engages the responsibility of his signatory.

This Reference Document incorporates by reference the preceding Reference Documents D.18-0041 filed on February 20, 2018 and D.17-0102 filed on 22 February 2017.

The following information is included by reference in the present Reference Document:

- The year 2017 consolidated financial statements of Sartorius Stedim Biotech prepared using international accounting standards and the report of the statutory auditors relating to these statements, and the Group 2017 management report appearing on pages 114 to 161 and 18 to 78 respectively, of the Reference Document filed with the Autorité des Marchés Financiers on 20 February 2018, under the number D.18-0041.
- The year 2016 consolidated financial statements of Sartorius Stedim Biotech prepared using international accounting standards and the report of the statutory auditors relating to these statements, and the Group 2016 management report appearing on pages 113 to 160 and 18 to 63 respectively, of the Reference Document filed with the Autorité des Marchés Financiers on 22 February 2017, under the number D.17 - 0102.

The sections of these documents not included are not of interest to an investor, and are covered in another part of this Reference Document.

Copies of the present Reference Document can be obtained from the following:

- Sartorius Stedim Biotech S.A.
Z.I. Les Paluds - Avenue de Jouques
CS 91051 - 13781 Aubagne Cedex
- Group website: www.sartorius-stedim.com
- Autorité des Marchés Financiers website:
www.amf-france.org

01 To Our Shareholders

| | |
|----|---------------------------------|
| 8 | Chairman's Message |
| 10 | Board of Directors |
| 12 | Sartorius Stedim Biotech Shares |

02 Management Report

| | |
|----|--|
| 18 | Structure and Management of the Group |
| 19 | Business Model, Strategy and Goals |
| 22 | Sector Conditions |
| 24 | Group Business Development |
| 29 | Net Worth and Financial Position |
| 34 | Sustainability |
| 35 | Opportunity and Risk Report |
| 47 | Forecast Report |
| 50 | Financial Statements of the Parent Company Sartorius Stedim Biotech S.A. as of December 31, 2018 |

This Reference Document contains statements concerning the future performance of Sartorius Stedim Biotech S.A. These statements are based on assumptions and estimates. Although we are convinced that these forward-looking statements are realistic, we cannot guarantee that they will actually apply. This is because our assumptions harbor risks and uncertainties that could lead to actual results diverging substantially from the expected ones. It is not planned to update our forward-looking statements.

This is a translation of the original French-language Reference Document "Document de Référence 2018". Sartorius shall not assume any liability for the correctness of this translation. The original French Reference Document is the legally binding version. Furthermore, Sartorius Stedim Biotech S.A. reserves the right not to be responsible for the topicality, correctness, completeness or quality of the information provided. Liability claims regarding damage caused by the use of any information provided, including any kind of information which is incomplete or incorrect, will therefore be rejected.

Throughout the Reference Document, differences may be apparent as a result of rounding during addition.

Contents

03 Corporate Governance Report

| | |
|----|---|
| 58 | The Board of Directors and Its Committees |
| 71 | Report on Corporate Governance |
| 72 | Corporate Governance Code/ AFEP MEDEF |
| 76 | Shareholders' Meetings |
| 78 | Delegation granted for increase in capital by the Shareholders' meeting to the Board of Directors |
| 79 | Remuneration of the Executive Director and Non-executive Members of the Board |
| 86 | Independent Auditors' Fees |

04 Consolidated Financial Statements and Notes

| | |
|-----|---|
| 90 | Statement of Profit or Loss and Other Comprehensive Income |
| 91 | Statement of Financial Position |
| 92 | Statement of Cash Flows |
| 93 | Statement of Changes in Equity |
| 94 | Notes to the Financial Statements |
| 102 | Notes to the Statement of Profit or Loss |
| 107 | Notes to the Individual Balance Sheet Items |
| 132 | Other Disclosures |
| 134 | Statutory Auditors' Report on the Consolidated Financial Statements |

05 Annual Financial Statements of Sartorius Stedim Biotech S.A. and Notes

| | |
|-----|--|
| 140 | Annual Financial Statements |
| 150 | Statutory Auditors' Report on the Financial Statements |

06 Supplementary Information

| | |
|-----|--|
| 154 | Other Information of a Legal Nature |
| 161 | Other Information on the Assets, Financial Position and Results for the Group |
| 164 | Special Report of the Statutory Auditors on Related Party Agreements and Commitments |
| 166 | Resolutions Submitted to the Annual General Shareholders' Meeting on March 26, 2019 |
| 173 | Report of the Board of Directors |
| 178 | Information on the Reference Document and the Annual Financial Report |
| 182 | Glossary |
| 184 | A Local Presence Worldwide |
| 187 | Financial Schedule |

To Our Shareholders

01

Chairman's Message

Dear Shareholders,

The publication of our Annual Report is always a good opportunity to stop for a moment and reflect upon the past year, not only from a financial perspective. In my letter to you last year, I explained a couple of effects that temporarily muted our growth in 2017, while our strong finish to the year already signaled that these effects would subside. Within this context, in 2018, our strong underlying growth trends became fully apparent again, and we even performed somewhat ahead of initial expectations. This led us to lift our targets quite substantially at mid-year. We have also been looking beyond our mid-term targets for 2020 and extended our time horizon to 2025 – more on that later.

Let us first turn to what we achieved in 2018: I am pleased to report that we closed the year with revenue growth in constant currencies of 13.7% and an increase in order intake of 12.5%. Operating profit rose by 16.1% to €342.4 million, and was mainly driven by economies of scale; the respective margin stood at 28.2%, 0.9 percentage points over the prior year. Underlying earnings per share were up, to €2.38 a gain of 21.5%. This dynamic performance was broad-based, encompassing both single-use products and equipment. Almost all of the increase in our sales revenue was achieved organically, while the acquisition of the software company Umetrics contributed around 0.5 percentage points of non-organic growth.

Regionally, the Americas led growth, reporting a gain of 16.9% to sales revenue of €508.2 million relative to a moderate previous-year base (reported +14.6%). Asia, especially China, again performed very strongly: Even relative to the high growth in the year-earlier period, sales revenue in Asia|Pacific again increased by double digits, 14.7%, to €281.3 million (reported +11.8%). That underscores the enormous growth potential of this region, where many previously underserved patients now begin to gain access to state-of-the-art medical drugs. The EMEA region recorded a solid gain of 10.7% to €508.2 million of sales revenue (reported +10.3%), which is excellent given the more mature markets and our already relatively high market share.

In light of these positive results, the Board of Directors will submit a proposal to the Annual General Shareholders' Meeting to raise dividends by 23.9% to

€0.57 per share. After development had been rather flat during 2017, the company's stock showed some volatility during 2018 and closed the year with a gain of 44.9%, strongly outperforming relevant indices such as the CAC 40, SBF 120 or NASDAQ Biotechnology.

Besides our strong financial performance, fiscal 2018 was also marked by important operational achievements, such as further development of our portfolio. We are pleased with how well our Umetrics data analytics business has been developing under our umbrella – as a stand-alone business, and regarding the progress we have made by integrating its data analytics technology into our equipment. Our customers are working intensely to better monitor, automate and predict their biomanufacturing processes and our Umetrics tools are helping them to make considerable strides in reaching these objectives.

Towards the end of 2018, we have modified our partnership in the area of cell culture media with our partner Lonza, the Swiss life science company, so that we both have relinquished our mutual exclusivity arrangement. While this modification will affect 2019 sales growth to some extent, it will create additional strategic leeway over the mid- and long term for building an integrated cell culture media franchise.

We significantly progressed in our multi-year investment program designed to nearly double manufacturing capacity of single-use bags and filters at our plant in Puerto Rico, create additional membrane casting capacity at our Göttingen site in Germany, and prepare to extend our aseptic bag site in Aubagne, France, and our cell line development site near Ulm, Germany. These major expansion projects are now nearing completion, and we anticipate lower investment requirements for the coming years. The Capex ratio is already set to decrease in 2019.

Ultimately, we have further developed our organization this year, gained 545 new employees, progressed with our digital agenda and are making strides in establishing our strong Sartorius brand even more firmly in the bioprocessing and life science tools community.



As we advance in 2019, we are striving to raise our sales revenue by about 7% to 11%. Without the effect from modification of the contract with Lonza, we estimate that sales revenue would be approximately 3 percentage points higher. Regarding profitability, we forecast that our operating profit margin will increase by slightly more than one percentage point over the prior-year figure of 28.2%. Both figures are stated in constant currencies. The projected increase in underlying EBITDA is expected to include an operational gain of approximately half a percentage point, whereas the remainder of this growth will result from changes to the IFRS accounting rules. The Capex ratio is projected to be at around 11% after 14.6% a year earlier.

Expectations for 2019 are well in line with our 2020 mid-term plan, which we still adhere to and which forecasts sales revenue to reach around €1.5 billion to €1.6 billion and our operating margin to attain about 29% to 30%. As 2020 approaches, we have extended our time horizon and defined our goals for 2025: we are again aiming to almost double revenue in this 5-year span to €2.8 billion and further expand our operating margin to around 30%.

Our targets remain ambitious because we have positioned ourselves well for the future and address a market with sustainable, fundamental growth drivers. We assume that demand for biopharmaceutical drugs will continuously increase due to a growing and aging population, improved access to medicines in emerging countries, an expanding market for biosimilars and due to completely new, emerging treatment options, such as personalized cell-based therapies. Regionally, the Asian market will play an increasingly important role, primarily China.

At the same time, cost-effective bioprocessing technologies will become even more essential as the biopharma market matures, become increasingly differentiated and also more competitive. Our customers are looking for partners that help them develop and produce innovative medicines faster and more effectively. In this context, we anticipate that the adoption of single-use systems will continue, encompassing a growing number of steps in our customers' value chain, and that these systems will progressively move from pre-commercial manufacture to commercial scale.

Sartorius Stedim Biotech will remain a prime vendor for this market and will continue to create innovative products and services, constantly enriching its portfolio through acquisitions, alliances and its own R&D. We will also further invest into our processes and systems to make doing business with us as convenient and easy as possible.

We could not have accomplished all this without our employees – now numbering 5.637 worldwide – who again did a remarkable job with their customer-first mentality and their fighting spirit. They drive our success, setting us apart in the bioprocessing world, and were key to making 2018 another successful year for Sartorius Stedim Biotech.

I also appreciate the continued trust of our customers, partners and shareholders, and cordially invite you to continue with us on the road to further achievements.

Sincerely

Joachim Kreuzburg

Chairman of the Board and CEO

Board of Directors



Board of Directors

The Board of Sartorius Stedim Biotech is the central management and supervisory entity of the company, and is composed of seven members. The directors are appointed for a three-year term.



Joachim Kreuzburg
Chairman | CEO



Susan Dexter



Anne-Marie Graffin



Lothar Kappich



Liliane de Lassus



Bernard Lemaître



Henri Riey

Sartorius Stedim Biotech Shares

Facts about the Share¹⁾

| | |
|----------------------|---|
| ISIN | FR0013154002 |
| Liquidity provider | Gilbert Dupont |
| Stock exchange | Euronext Paris |
| Market segment | Local Securities - Compartment A (Large Caps) |
| Indexes | SBF 120; CAC All-Tradable; CAC All SHARES; CAC MID & SMALL 190; CAC HEALTHCARE; STOXX Europe 600; MSCI France |
| Number of shares | 92,180,190 |
| thereof Sartorius AG | 74.3% |
| thereof free float | 25.7% |
| Voting rights | 161,024,614 |
| thereof Sartorius AG | 85.0% |
| thereof free float | 15.0% |

¹⁾ As of December 31, 2018

Persistent and High Volatility in Stock Markets

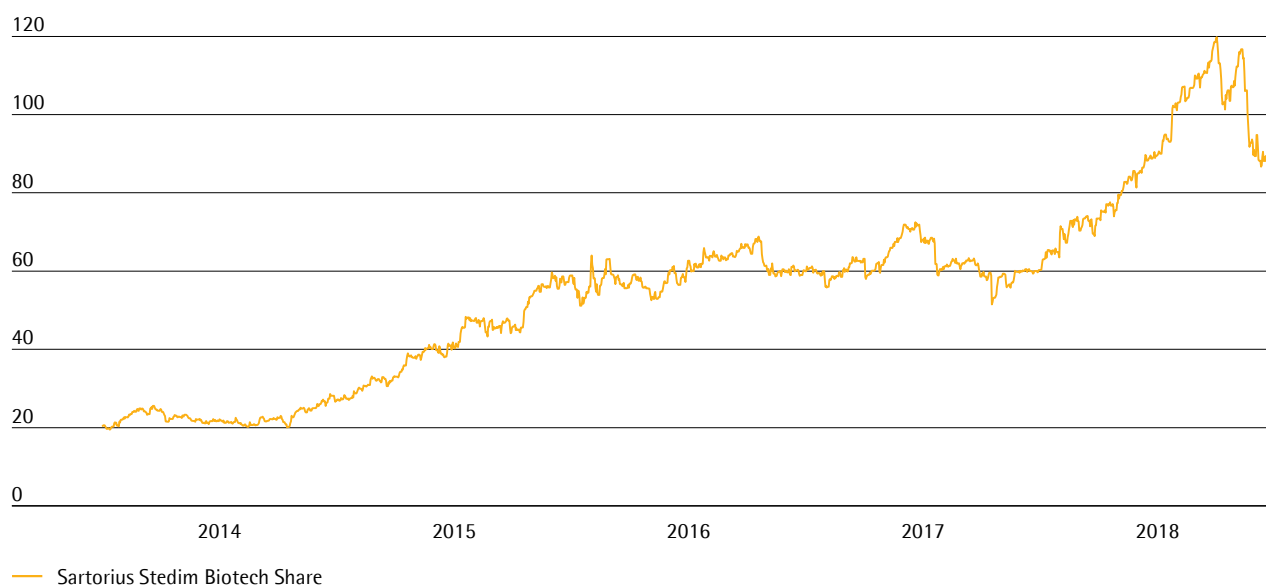
Gains and relatively steady growth among the international stock exchanges marked the year 2017. In contrast, the year under review was characterized by substantial fluctuations and falling share prices. Political uncertainties, international trade conflicts and the interest-rate reversal in the United States caused high volatility that also affected the leading French stock index. The CAC 40 reached an all-time high of more than 5,600 points in May 2018, then fell considerably, finishing the 2018 stock-market year down 11.0% at 4,730 points. The SBF 120, home to Sartorius Stedim Biotech shares, declined 11.7%. The index relevant to the biotech industry, NASDAQ Biotechnology, registered a drop of 9.3%.

Sartorius Shares Rise Substantially

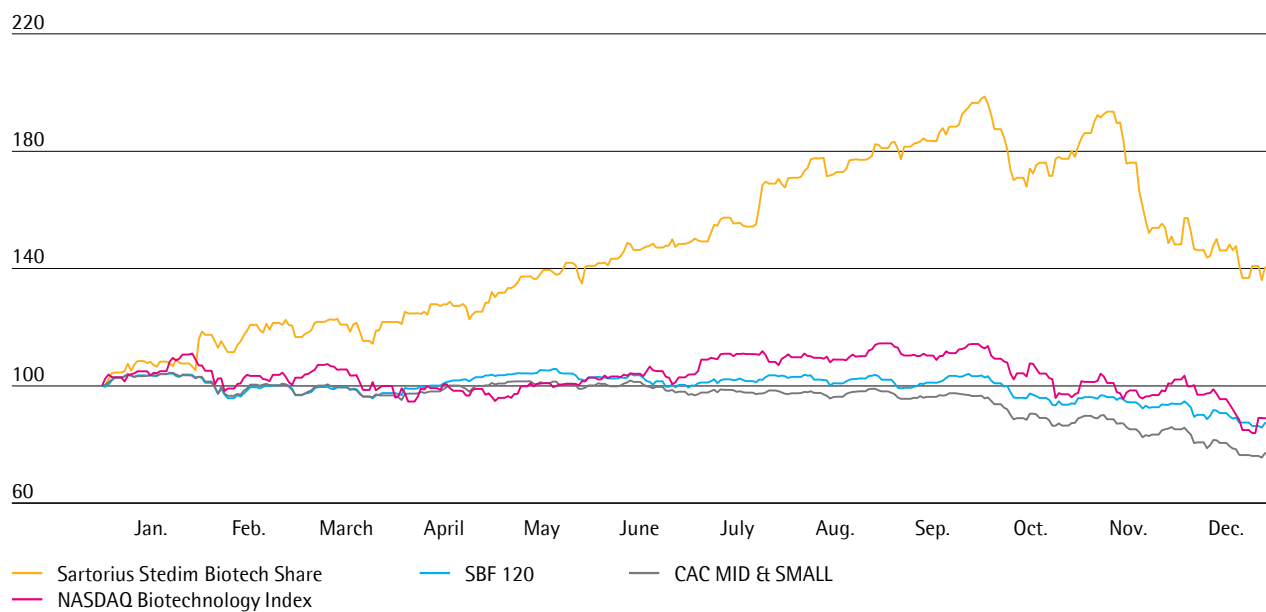
The Sartorius Stedim Biotech share price developed positively in spite of the volatile stock exchanges. Contributing factors included the announcement of medium-term financial targets for 2025 and better-than-expected business performance halfway through 2018. In the first half of the year the share recorded significant gains and reached an all-time high of €119.80 on October 02, 2018. In the following months the share lost in value, but finished the 2018 stock-market year at a closing price of €87.35 – up 44.9% year on year.

Sartorius Stedim Biotech Share in €¹⁾

January 1, 2014, to December 31, 2018

¹⁾ January 1, 2014, to May 9, 2017, adjusted for stock split**Sartorius Stedim Biotech Share¹⁾ in Comparison to the SBF 120, CAC MID & SMALL and NASDAQ Biotechnology Index (indexed)**

January 1, 2018, to December 31, 2018



Investor Relations Activities

Sartorius Stedim Biotech's investor relations activities follow the objective of making the current and future development of the company transparent for its stakeholders. To achieve this objective, Sartorius maintains an ongoing, open dialog with shareholders, potential investors and financial analysts.

Besides providing quarterly, first-half and annual reports, we inform the capital market and the interested public at quarterly teleconferences and in regularly published press releases about the current development of our business and other material events at the company. Moreover, Group management and our IR team were present for talks and factory tours with capital-market participants at our sites in Aubagne, France, and in Göttingen, Germany. Management and IR specialists also took part at conferences and roadshows in various financial centers. In February 2018, we hosted a capital market day, with a focus on discussing our medium-term financial targets for 2025.

All information and publications relating to our company and its shares may be found on our website at www.sartorius-stedim.com.

Analysts

The recommendations of financial analysts serve as a foundation for the decisions of private and institutional investors when investing in shares. Currently, seven institutions regularly prepare reports and updates on Sartorius Stedim Biotech shares.

Research Coverage

| Date | Company | Recommendation | Target price in € |
|------------------|---------------------|----------------|-------------------|
| January 30, 2019 | AlphaValue | Sell | 88.40 |
| January 30, 2019 | Société Générale | Hold | 105.00 |
| January 30, 2019 | Janney | Hold | 121.00 |
| January 29, 2019 | Oddo BHF | Hold | 90.00 |
| January 24, 2019 | New Street Research | Buy | 100.00 |
| October 23, 2018 | Gilbert Dupont | Sell | 107.00 |
| August 07, 2018 | UBS | Buy | 115.00 |

Key Figures for Sartorius Stedim Biotech Share¹⁾

| | | 2018 | 2017 | 2016 | 2015 | 2014 |
|---|----------------|---------|---------|---------|---------|---------|
| Share price ²⁾ in € | Reporting date | 87.35 | 60.29 | 59.97 | 58.90 | 26.88 |
| | High | 119.80 | 72.49 | 68.84 | 59.67 | 28.61 |
| | Low | 60.35 | 51.50 | 51.17 | 26.89 | 19.50 |
| Dividends ³⁾ in € | | 0.57 | 0.46 | 0.42 | 0.33 | 0.22 |
| Total dividends paid ³⁾ in millions of € | | 52.5 | 42.4 | 38.7 | 30.7 | 20.0 |
| Payout ratio ^{3,4)} in % | | 24.0 | 23.5 | 21.9 | 22.1 | 22.9 |
| Dividend yield ⁵⁾ in % | | 0.7 | 0.8 | 0.7 | 0.6 | 0.8 |
| Market capitalization in millions of € | | 8,051.9 | 5,557.5 | 5,528.0 | 5,430.8 | 2,477.4 |
| Average daily trading number of shares | | 80,140 | 52,753 | 46,752 | 44,115 | 42,084 |
| Trading volume of shares in millions of € | | 1,874.9 | 818.2 | 714.2 | 485.2 | 243.8 |
| CAC MID & SMALL (closing prices of the year) | | 11,337 | 14,456 | 11,848 | 11,054 | 9,354 |
| SBF 120 (closing prices of the year) | | 3,756 | 4,251 | 3,836 | 3,664 | 3,360 |

¹⁾ For 2014 to 2015, share prices, dividends and average daily trading number of shares adjusted for stock split; rounded values

²⁾ Daily closing price

³⁾ For 2018, amounts suggested by the Board of Directors and subject to approval by the Annual General Shareholders' Meeting

⁴⁾ Based on the underlying net result

⁵⁾ Dividends in relation to the corresponding closing prices of the year

Sources: Euronext; Bloomberg

Dividends

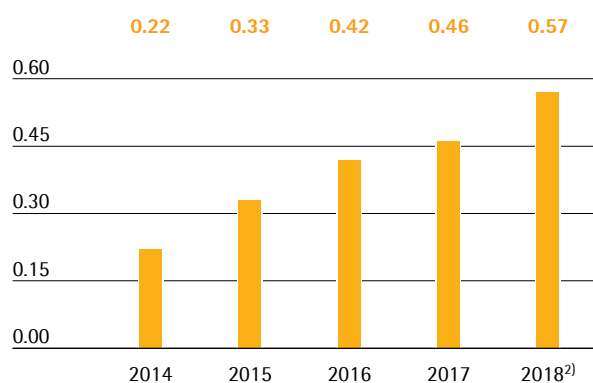
Sartorius Stedim Biotech strives to enable its shareholders to participate adequately in the company's success and has continuously increased its dividend in recent years. In line with this objective, we basically follow the policy of paying out a relatively stable share of relevant net profit to our shareholders.

Relevant net profit

The Board of Directors will submit a proposal to the Annual General Shareholders' Meeting on March 26, 2019, to pay a dividend from the underlying net profit of €219.3 million for fiscal 2018 of €0.57 per share compared to previous year's figure of €0.46. If approved, the dividend would increase for the ninth consecutive year and the total profit distributed would rise by a considerable 23.9%, from €42.4 million last year to €52.5 million. The corresponding dividend payout ratio would be 24.0% compared to 23.5% in the previous year. In relation to the shares' closing price of €87.35 on December 31, 2018, the dividend yield would be 0.7% (previous year: 0.8%).

Dividends¹⁾

in €



¹⁾ 2014 to 2015 adjusted for stock split; rounded values

²⁾ Amount suggested by the Board of Directors and subject to approval by the Annual General Shareholders' Meeting

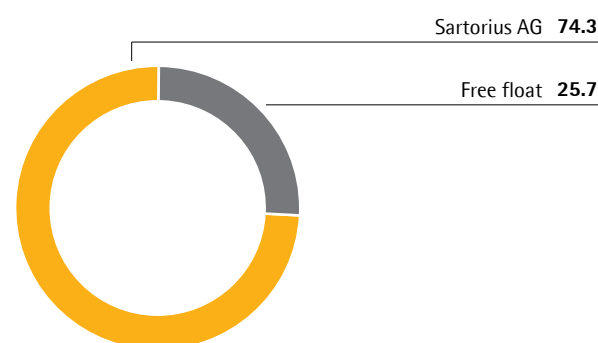
Shareholder Structure

As a consequence of the stock split and the increase of the individual par value of the company's shares, Stedim Biotech S.A.'s issued capital amounted to €18.4 million as of December 31, 2018, and was divided into 92,180,190 shares, each with a calculated par value of €0.20. As some of the shares convey double voting rights, there were a total of 161,024,614 voting rights as of the reporting date.

As of December 31, 2018, Sartorius AG holds 74.3% of the share capital and 85.0% of the outstanding voting rights. The remaining 25.7% of the shares are in free float; corresponding to 15.0% of the outstanding voting rights.

Shareholding Structure

in % of share capital



Management Report

02

Structure and Management of the Group

Group Legal Structure

Sartorius Stedim Biotech is a globally operating company with subsidiaries in more than 20 countries and more than 5,600 employees worldwide. The parent company of the Sartorius Stedim Biotech Group is Sartorius Stedim Biotech S.A., headquartered in Aubagne, France.

Sartorius Stedim Biotech S.A. is listed on the Euronext stock exchange in Paris. Approximately 74% of the share capital and around 85% of the voting rights of Sartorius Stedim Biotech S.A. are held by Sartorius AG.

Sartorius AG is a leading international bioprocess and laboratory equipment and technology provider headquartered in Göttingen, Germany. It is listed on the German Stock Exchange and operates two divisions: the bioprocess business as a subgroup under its parent corporation Sartorius Stedim Biotech S.A. and the laboratory business as a further subgroup.

The consolidated financial statements of the Sartorius Stedim Biotech Group include Sartorius Stedim Biotech S.A. and all affiliates in which Sartorius Stedim Biotech S.A. has a controlling interest pursuant to IFRS 10.

Organization and Management of the Group

The Sartorius Stedim Biotech Group is largely organized by function on a worldwide basis. Accordingly, the respective management responsibilities are performed along the company's core functions across all sites and regions.

This global functional organization forms an effective platform for central strategic control and for fast, efficient collaboration and execution within the Group. It enables the company to realize its total solutions provider strategy and position itself effectively in respect of global customers.

The Board of Directors of Sartorius Stedim Biotech S.A. is composed of seven members, one executive director and six non-executive directors.

Implementing the Group's various strategies and initiatives at the local level is the responsibility of the national affiliates. The management bodies of the local companies run their organizations in accordance with applicable statutory provisions, articles of association and rules of procedure and in keeping with the principles of corporate governance that apply throughout the Sartorius Stedim Biotech Group worldwide. Please see details of the Board of Directors in the section "Corporate Governance."

Financial Controlling and Key Performance Indicators

The Sartorius Stedim Biotech Group is managed using a number of key performance indicators, which are also decisive for determination of the variable remuneration component for the Executive Committee and managers.

The key management parameter that Sartorius Stedim Biotech uses to measure the development of its size is currency-adjusted growth of sales revenue.

The key performance measure for profitability is EBITDA adjusted for extraordinary items, i.e. underlying EBITDA, and the corresponding margin. For a definition of this term and more information on its presentation, see the Glossary on page 182.

Regarding the debt capacity of the Sartorius Stedim Biotech Group, a further key indicator is the ratio of net debt to underlying EBITDA for the last twelve months.

Moreover, the capex ratio, i.e. capital expenditures relative to sales revenue, represents a key control parameter.

The following financial and non-financial indicators are also reported on a regular basis:

- Order intake
- Underlying net profit | Earnings per share
- Net profit | Earnings per share
- Equity ratio
- Net working capital
- Net cash flow from operating activities
- Number of employees

The annual financial forecast published at the beginning of a fiscal year for the Group generally refers to the development of sales revenue and of underlying EBITDA margin. The expected capex ratio, as well as a directional forecast for the ratio of net debt to underlying EBITDA, is also indicated for the Group.

Business Model, Strategy and Goals

Market and Strategic Positioning

As a leading partner of the biopharmaceutical industry, we help our customers to develop their production processes and manufacture medications more efficiently. Our objective is to propel science forward and contribute toward enabling more people to have access to healthcare.

We are a leading player on this attractive market, which is characterized by strong growth momentum and long-term trends. Primary growth drivers are a growing world population and an increase in age-related diseases in industrialized countries. In addition, rising incomes in emerging countries are leading to improved access to healthcare and rising demand for medications. Medical advances are also driving the ongoing development and approval of new biopharmaceuticals as well as of emerging biosimilars, which are highly similar to established biologics already licensed, and account for a share of the biopharma market that is currently still small, but fast-growing. As a result of these factors, the volumes of biotech medications are steadily increasing and so is the demand for the appropriate production technologies. This market is largely independent of business cycles.

The maturity and intensity of competition in this still comparably young biopharmaceutical industry are successively increasing. In addition to achieving scientific success, our customers will find it more important, in view of mounting cost pressure on healthcare systems, to increase the efficiency of their research, development and manufacturing processes. We help them meet this challenge by further developing our product portfolio. One of the decisive success factors of Sartorius Stedim Biotech is to use technology in order to differentiate ourselves from our competitors. Our innovative power rests on three pillars: our own specialized product development, the integration of innovations via acquisitions, and alliances with partners in complementary fields.

Another competitive advantage of Sartorius Stedim Biotech is its broad understanding of applications, which is based on its clear focus on the sector. We are thoroughly familiar with our customers' entire added-value chains, especially the interactivity of the systems in these chains. All this makes us a strategic partner of these customers who drive forward innovations in bioprocess technology.

Products & Services

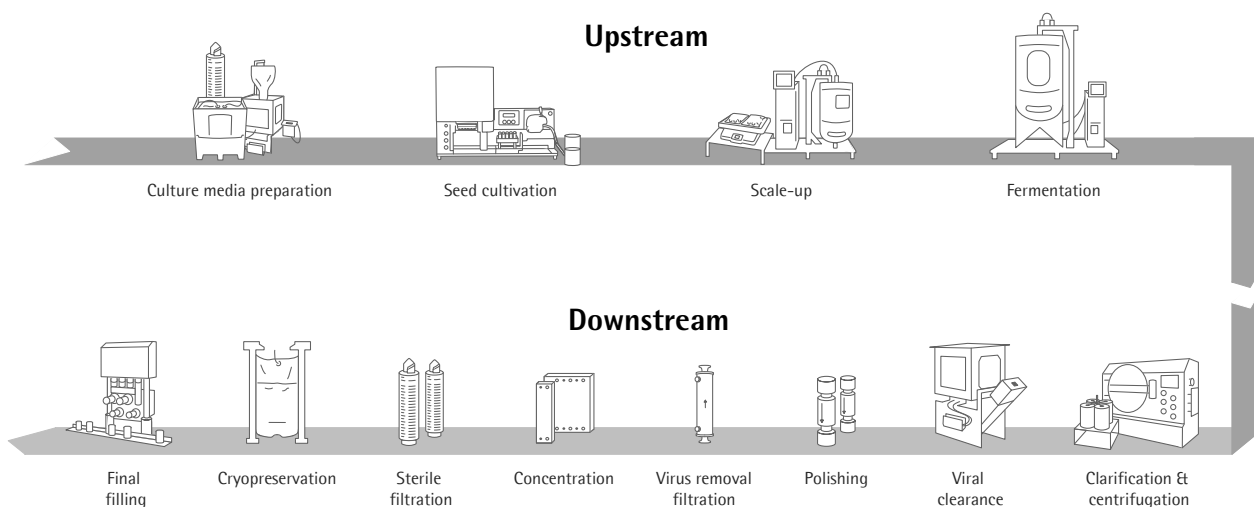
We are offering a broad portfolio of products that focuses on all major steps in the manufacture of a biopharmaceutical, as well as in process development as prerequisite procedures. Our technologies cover, inter alia, cell line technologies, cell culture media, bioreactors, and a wide range of products for separation, purification and concentration of biological intermediates and finished products, as well as solutions for their storage and transportation. Sartorius Stedim Biotech also offers data analytics software for modeling and optimizing processes of biopharmaceutical development and production. In its core technologies, the company has leading market positions with high double-digit market shares.

The breadth of our product portfolio sets us apart from our competitors. We provide customers with an entire production unit from a single source, as well as assist with preceding project planning, process integration and subsequent validation.

As an innovation leader, Sartorius Stedim Biotech was moreover one of the first biopharmaceutical suppliers to specialize in single-use technologies, which account for approximately three-quarters of the Group's sales revenue. Due to their cost advantages and their greater flexibility and safety compared with reusable technologies, the pharmaceutical industry is increasingly relying on single-use products. Particularly in pre-commercial production processes, single-use products have almost completely

supplanted classic stainless steel components. Industry observers believe that market penetration is likely to continue as commercial production also increasingly moves toward single-use products. As a result, we are generating a large share of sales from repeat business. The high approval requirements placed on our customers' products are also contributing to this growth. Because our customers' production processes must be validated by the health authorities responsible, the technological components initially used can be replaced only at considerable expense once they have been approved. The manufacturers of medications are therefore closely tied to the suppliers for the life cycle of a medication. Beyond this, our broad and stable customer base that we address through our specialized sales force directly for the most part also contributes to this favorable risk profile.

The strong strategic positioning of Sartorius Stedim Biotech and the above-average expansion of the sector are a good foundation for profitable growth in the future as well. Beyond realizing our organic development potential, we also aim to further expand the portfolio through complementary acquisitions and alliances.



Sartorius Stedim Biotech 2020 and 2025 Strategies

In 2011, Sartorius Stedim Biotech presented its strategy and targets for profitable growth up to 2020 according to which sales revenue is projected to increase in a range of around €1.5 to €1.6 billion with an underlying EBITDA margin of about 29% to 30%. At a Capital Markets Day in February of the reporting year, management extended its time horizon, introducing its strategy and long-term targets up to 2025.

While the targets for 2020 still continue to apply unchanged, Sartorius Stedim Biotech plans to increase its revenue in the period of 2020 to 2025 to around €2.8 billion, given the high market dynamics and the company's strong strategic positioning. The underlying EBITDA margin is forecasted to increase to around 30%.

In these projections, management takes into account that any future acquisitions in the Group would initially be margin dilutive and that no significant changes in key exchange rates would occur.

These targets are being implemented by various growth initiatives with the following focal points:

Expansion of the Product Portfolio

Sartorius Stedim Biotech offers a broad product portfolio that is continuously expanded in line with the value-added chain of the biopharmaceutical industry. Aside from our own research and development activities and strategic partnerships, acquisitions that are complementary to or extend our strengths appropriately will remain part of our strategy. We see opportunities in digital networking of products, for example, in the integration of software solutions for bioprocess production control, among others. Expansion into adjacent applications, such as regenerative medicine, is also conceivable. At the focus of our efforts will be products that offer solutions to the challenges our customers face and that make our offering even more attractive from the customers' perspective.

Regional Growth Initiatives

North America and Asia are the key focal areas of our regional growth strategy.

North America is the world's largest market for bioprocess equipment. Yet because it is home to our main competitors, Sartorius Stedim Biotech has lower market share in this region than in Europe and Asia. Accordingly, the company is striving to gain additional market share, primarily by strengthening its sales and service capacities.

A further strategic focus is on China. This market offers sizable growth potential owing to rising private and public healthcare expenditures and the rapid development of regional biopharmaceutical plants. To benefit from the dynamic development of this market, Sartorius Stedim Biotech has already been investing heavily in its sales infrastructure and plans to expand production capacity levels there over the medium term.

Optimization of Work Processes

Sufficient production capacity and a powerful supply chain are an essential foundation of future growth. For this reason, in recent years Sartorius Stedim Biotech has substantially expanded its capacities for membranes, filters and single-use bags at various Group sites. Beyond this, a new ERP system based on Group-wide standardized business processes was introduced, the implementation of which is due to be completed in 2019.

Following these significant infrastructural expansions, our focus is increasingly shifting to optimization of our processes. Thus, we are driving forward digitalization and process automation in all parts of the company to further enhance the performance power of our supply chain and our customer contact interfaces. This also includes extending our activities in the areas of e-commerce, digital marketing and analytics.

Sector Conditions

Sartorius Stedim Biotech serves customers mainly in the biopharmaceutical industry, which makes its business particularly sensitive to the development of this industry.

Strong Growth in the Biopharmaceutical Market

According to estimates from several market observers, the global pharmaceutical market showed a positive development once again in 2018, with an increase of approximately 4% to 5%. Within the pharmaceutical market, the segment for medications and vaccinations manufactured using biotech methods has grown faster than the rest of the market for many years now. In 2018, the biopharmaceutical market was estimated at a volume of €217 billion, an increase of approximately 8% to 9% over the previous year. The steadily growing significance and acceptance of biopharmaceutical drugs is reflected in its increasing share of the sales revenue in the global pharmaceutical market and the development activities of the pharmaceutical industry. For example, biopharmaceutical compounds account for more than 40% of the R&D pipeline.

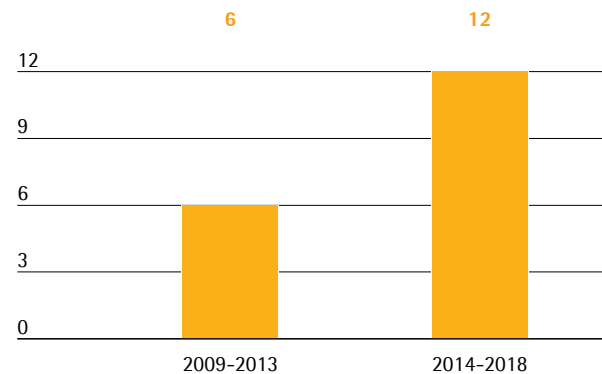
It is estimated that nearly €13 billion in sales were generated in 2018 with products and technologies for the manufacture of biopharmaceuticals. In contrast to its moderate growth in 2017, the supplier industry considerably increased sales in the year under review. There was once again a dynamic rise in demand in emerging countries, especially in China and India, where many commercial biopharmaceutical production facilities have opened in recent years to meet domestic demand. But there was also considerable investment in the United States and Europe, where the highest bioreactor capacities are located.

Market growth fundamentally depends more on medium- to long-term trends than on short-term economic developments. The major growth driver is the increasing demand worldwide for medications. In addition, the approval and market launch of new biopharmaceuticals boosts growth. In the year under review, the U.S. Food and Drug Administration (FDA), approved a record 17 biopharmaceuticals. A growing number of active pharmaceutical ingredients (APIs) manufactured using biotech production methods is being approved for the treatment of rare illnesses that have been untreatable so far. There has been recent progress in cell and gene therapies: the United States and Europe have now granted market approval to three therapies. A growing number of approved biopharmaceuticals as well as an increasing variety of

therapy types and API classes coupled with growing demand for medications are the main drivers for the worldwide increase in production capacities for biopharmaceuticals.

Biosimilars, or generic versions of biologics similar to originally patented medications, are playing an increasingly important role in the biotechnology market. Although sales volume was comparatively low at an estimated €6 billion in 2018, the market is expected to generate an average annual growth rate of around 30% until 2022. Through the approval and market launch of new biosimilars, further progress was made in the reporting period. The biosimilars market is projected to continue growing in coming years due to the expiration of patents for a number of high-margin biopharmaceuticals. In addition, the regulatory, patent law-related and marketing challenges and hurdles that have hindered faster market penetration of biosimilars to date are likely to decrease gradually.

Average Number of New Approvals of Biotech Medications in the USA per Year



Single-Use Systems for Biopharmaceutical Production Continue to Gain Importance

Biotech production methods are much more complex and cost-intensive than traditional methods for producing medications. Consequently, manufacturers and suppliers are continually looking to develop more efficient technologies. Single-use products play a decisive role in this effort: They require significantly less capital expenditure, reduce costs for cleaning and validation and minimize downtime. They also offer greater flexibility in production and help accelerate time to market.

Owing to these advantages, single-use technologies are already an established part of many process steps. Single-use systems are primarily employed in pre-commercial development activities and production phases and in small-batch manufacturing. It can be expected that single-use technologies will become increasingly popular for the production of high-volume commercial quantities. This is particularly relevant to the production of biotech drugs, whose clinical development takes place in single-use systems. Sartorius offers the sector's most extensive portfolio of single-use technologies, with scalable products for every step of manufacturing. The company is also actively helping the biopharmaceutical industry convert to these technologies in the production of medications.

Moderate Growth in the Global Laboratory Market

According to estimates by the market research firm Frost & Sullivan, the global laboratory market grew by approximately 3.6% to €39 billion in the reporting year. Following a lackluster first quarter, budget increases for academic and public research institutes during 2018 led to rising demand. In addition, investing activities in the private sector remained stable due to the positive economic situation.

While Europe saw an increase of 3.1%, the United States, the largest market for laboratory products, recorded growth of 3.3%. Significant growth was once again reported in Asian countries, such as China and India, in which the laboratory market enjoyed an above-average expansion of 7.5% (China) and 8.5% (India).

Competition

The primary means by which companies in the biotechnology market differentiate themselves from competitors are innovative process and the quality and performance of their products. The biotechnology sector is constantly discovering new areas of application and expects suppliers to be equally fast-moving and creative in developing new equipment for the manufacture of biotech products. New suppliers, in particular, seek to capitalize on the opportunities inherent in this environment to gain a foothold in the market with carefully targeted niche products. The more established suppliers, meanwhile, are expanding their product range continuously.

We generate around 90% of our sales revenue from validated processes in which replacing products during the production cycle is very expensive, so we receive a high proportion of follow-up and repeat business. The

particular strength of Sartorius Stedim Biotech lies in its integrated process solutions: we offer the broadest range in the industry, from drug discovery and development in the lab to commercial manufacture of the final product. Our strategic focus on single-use products gives us a further edge over the competition. Sartorius Stedim Biotech has leading positions on the global market for bioprocess filtration, fermentation, cell cultivation, fluid management and membrane chromatography.

Most of our competitors are multinationals based in the USA. Certain business units of Merck KGaA, Danaher Corp., General Electric Company and Thermo Fisher Scientific Inc. are among our main rivals in the process area; Thermo Fisher and Merck are key players in the laboratory field. We also face competition from smaller companies in individual segments.

Sources: IQVIA Institute: 2018 and Beyond: Outlook and Turning Points, March 2018; Evaluate Pharma: World Preview 2018, Outlook to 2024, June 2018; BioPlan: 15th Annual Report and Survey of Biopharmaceutical Manufacturing Capacity and Production, April 2018; Frost & Sullivan: 2018 Mid-year Report: Forecast and Analysis of the Global Market for Laboratory Products, May 2018; BCC Research: Biosimilars: Global Markets, March 2018; Daedal Research: Global Biologics Market: Size, Trends & Forecasts, February 2018; www.fda.gov

Group Business Development

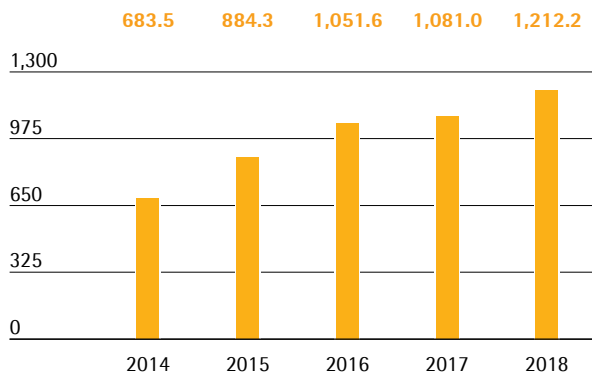
Sales Revenue and Order Intake

In fiscal 2018, Sartorius Stedim Biotech continued on the growth track, with double-digit gains in sales revenue and order intake. The dynamic growth was based on its very competitive product portfolio and fueled by strong demand across all product categories and geographies. Hence, after the previous year's comparably moderate performance, momentum considerably picked up in the reporting period. Group's sales revenue rose in constant currencies by 13.7% to €1,212.2 million (reported: +12.1%). Sartorius Stedim Biotech therefore reached the upper end of its guidance which had been revised upwards at mid-year 2018 to 11% to 14%. Most of this increase was organic, whereas the acquisition of the software company Umetrics contributed around 0.5 percentage points of non-organic growth.

Order intake outperformed sales, posting an increase of 14.2% in constant currencies to €1,307.3 million. On a reported basis, this gain was 12.5%.

Sales Revenue 2014 to 2018

in € million



Sales Revenue and Order Intake

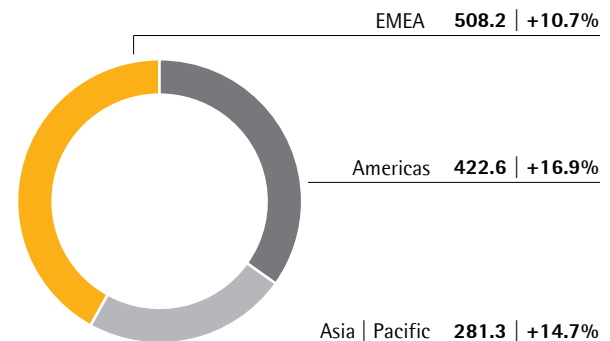
| in € millions | 2018 | 2017 | in % reported | in % const. fx |
|---------------|---------|---------|------------------|-------------------|
| Sales Revenue | 1,212.2 | 1,081.0 | 12.1 | 13.7 |
| Order Intake | 1,307.3 | 1,162.3 | 12.5 | 14.2 |

Geographically, all regions contributed to consolidated growth. EMEA, the region generating the highest revenue for the company with around 42% of its total sales, recorded a gain of 10.7% to €508.2 million against a moderate prior-year base. In the Americas region, sales increased dynamically by 16.9% to

422.6 million after a decline in the previous year, now representing around 35% of revenue. The Asia | Pacific region, which accounted for around 23% of the Group's sales, also grew significantly after an already strong performance in 2017: Sales rose in the clear double-digits by 14.7% to €281.3 million. All growth rates are in constant currencies unless otherwise stated.

Sales Revenue and Growth¹⁾ by Region²⁾

in € millions unless otherwise specified



¹⁾ In constant currencies

²⁾ Acc. to customers' location

Development of Costs and Earnings

In the reporting year, the cost of sales stood at €582.6million. In comparison with sales revenue growth of 12.1%, the cost of sales increased underproportionately by 10.7%, which was due to product mix effects and economies of scale. The cost of sales ratio was 48.1% relative to 48.7% a year ago.

Selling and distribution costs rose by 10.2% to €215.2million so the ratio of these costs to sales revenue decreased from 18.1% in the previous year to 17.8%.

Expenses for research and development rose year over year by 13.9% to €60.6million. The ratio of R&D expenses to sales revenue was 5.0%, slightly above the prior-year level of 4.9%.

Concerning general administrative expenses, Sartorius Stedim Biotech reported an increase of 8.6% to €67.0million. In relation to sales revenue, general administrative expenses decreased from 5.7% in the previous year to 5.5% in the reporting year.

In fiscal 2018, the balance of other operating income and expenses significantly improved to €13.5million relative to -€23.0million a year earlier. This year-over-year change was mainly driven by significant increases in extraordinary items. In the previous year, this balance totaled -€22.6million and was due to various corporate projects, expenses related to the most recent acquisitions and to the consequences of Hurricane Maria. In the reporting year, modification of the contract with the life science company Lonza on the exclusive distribution and marketing rights for specific cell culture media and buffers resulted in income of €35.2million so that extraordinary items totaled €12.7million.

In the year under review, the Group's EBIT increased strongly by 35.4% to €300.2million, especially due to higher sales revenue and to the improvement mentioned above in the other operating income and expenses. As a consequence and despite slightly higher depreciation, the Group's EBIT margin rose from 20.5% a year earlier to 24.8%.

The financial result was -€15.7million in 2018 relative to -€1.1million in 2017. This was essentially attributable to valuation effects related to foreign currency liabilities and hedging instruments.

In the reporting year, tax expenses of €74.6million were higher than the prior-year total of €56.8million. The company's tax rate was 26.2% compared to 25.8% in the year before.

In the reporting year, net profit attributable to shareholders of Sartorius Stedim Biotech S.A. amounted to €208.1million relative to €161.1million a year earlier.

Statement of Profit or Loss

| € in millions | 2018 | 2017 | in % |
|---|---------|---------|-------|
| Sales revenue | 1,212.2 | 1,081.0 | 12.1 |
| Cost of sales | -582.6 | -526.2 | -10.7 |
| Gross profit on sales | 629.6 | 554.8 | 13.5 |
| Selling and distribution costs | -215.2 | -195.2 | -10.2 |
| Research and development costs | -60.6 | -53.2 | -13.9 |
| General administrative expenses | -67.0 | -61.7 | -8.6 |
| Other operating income and expenses | 13.5 | -23.0 | 158.7 |
| Earnings before interest and taxes (EBIT) | 300.2 | 221.7 | 35.4 |
| Financial income | 5.3 | 9.5 | -43.7 |
| Financial expenses | -21.0 | -10.6 | -98.6 |
| Financial result | -15.7 | -1.1 | n.m. |
| Profit before tax | 284.5 | 220.6 | 29.0 |
| Income taxes | -74.6 | -56.8 | -31.2 |
| Net result | 210.0 | 163.8 | 28.2 |
| Attributable to: | | | |
| Equity holders of SSB S.A. | 208.1 | 161.1 | 29.2 |
| Non-controlling interest | 1.9 | 2.7 | -28.7 |

Earnings

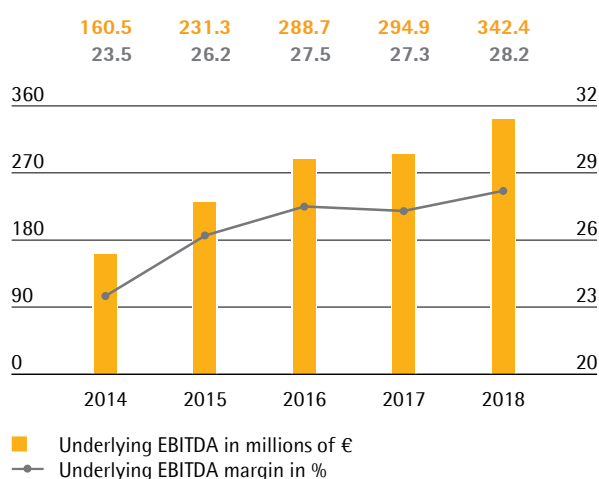
At the Sartorius Stedim Biotech Group, earnings before interest, taxes, depreciation and amortization (EBITDA) are used as the key profitability measure. To provide a complete and transparent picture of the Group's profitability, also in an international comparison, we report earnings adjusted for extraordinary items (underlying EBITDA). For more information about definitions, please refer to the Glossary on page 182. The underlying presentation is reconciled with the EBITDA key indicator (see Glossary) as follows:

Reconciliation between EBIT and underlying EBITDA

| € in millions | 2018 | 2017 |
|-------------------------------|-------|-------|
| EBIT | 300.2 | 221.7 |
| Extraordinary items | -12.7 | 22.6 |
| Depreciation and amortization | 54.9 | 50.6 |
| Underlying EBITDA | 342.4 | 294.9 |

In fiscal 2018, Sartorius Stedim Biotech increased its earnings significantly. Underlying EBITDA thus rose by 16.1% to €342.4 million. The Group's respective underlying EBITDA margin improved to 28.2% (2017: 27.3%) due to economies of scale, thus exceeding our forecast which had been revised upwards at mid-year.

Underlying EBITDA and margin¹⁾

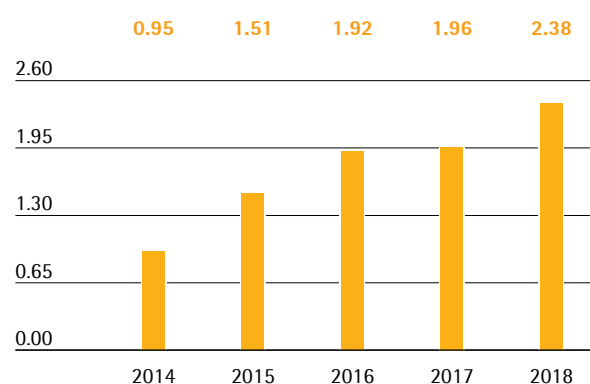


¹⁾ Adjusted for extraordinary items

The underlying net result after non-controlling interest for the Group rose strongly from €180.4 million a year ago to €219.3 million in fiscal 2018. This figure is the basis for calculating the profit to be appropriated and is calculated by adjusting for extraordinary items, eliminating non-cash amortization of €16.8 million (previous year: €16.6 million), and is based on the normalized financial result (see Glossary), as well as the corresponding tax effects for each of these items. Underlying earnings per share surged by 21.5% from €1.96 a year earlier to €2.38.

Underlying Earnings per Share¹⁾²⁾

in €



¹⁾ Excluding extraordinary items

²⁾ 2014 to 2015 adjusted for stock split; rounded values

| € in millions | 2018 | 2017 |
|---|--------------|--------------|
| EBIT (operating result) | 300.2 | 221.7 |
| Extraordinary items | -12.7 | 22.6 |
| Amortization IFRS 3 | 16.8 | 16.6 |
| Normalized financial result¹⁾ | -5.3 | -6.6 |
| Normalized income tax (2018: 26%, 2017: 28%) ²⁾ | -77.7 | -71.2 |
| Underlying net result | 221.2 | 183.1 |
| Non-controlling interest | -1.9 | -2.7 |
| Underlying net result after non-controlling interest | 219.3 | 180.4 |
| Underlying earnings per share (in €) | 2.38 | 1.96 |

¹⁾ Financial result excluding fair value adjustments of hedging instruments, as well as currency effects from foreign currency loans

²⁾ Underlying income tax, based on the underlying profit before taxes and non-cash amortization

See Glossary for the definitions of the totals listed above.

Research and Development

Activities in product development at Sartorius Stedim Biotech encompass both new and advanced in-house product developments in our own core technologies as well as the integration of new products through collaborations and acquisitions.

Development activities focus on technology areas such as membranes, which are the core component of all types of filter products; various base technologies, such as single-use containers and sensors; and control technologies for processes such as fermentation. Additional focal areas include developments in materials and components such as plastic, elastomers and intelligent polymers; expanded data analysis; and cell-line development.

In the reporting year, one emphasis was on developing the next generation of integrity-testing applications and a system for virus filtration that can be integrated as a single-use solution, even in stainless-steel units.

Another key task in 2018 was the ongoing integration into our systems of software from Umetrics, which Sartorius Stedim Biotech acquired in 2017. It should become possible in 2019 to use the programs in the downstream process, as well.

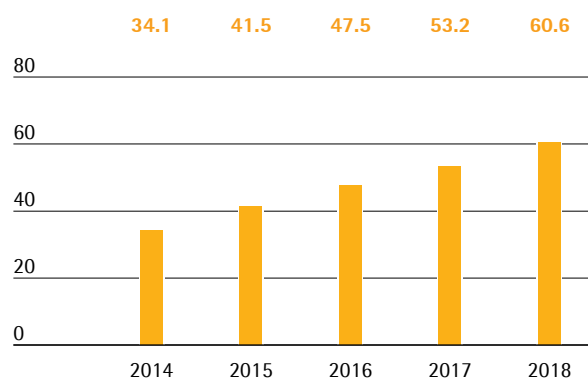
In addition, we are investing in the development of applications for regenerative medicine and plan to launch an analytical solution for lentiviruses, which include human immunodeficiency (HI) viruses. We also intend to launch an innovative single-use centrifuge and a membrane-based chromatography system on the market.

Our largest site for product development is Göttingen; other key sites are Aubagne (France), Guxhagen (Germany), Bangalore (India), Bohemia, New York (United States), Royston (United Kingdom) and Umeå (Sweden).

The Sartorius Stedim Biotech Group stepped up its research and development activities in the reporting year, increasing spending in this area by 13.9% to €60.6 million (previous year: €53.2 million). The ratio of R&D costs to sales revenue slightly increased to 5.0% compared to 4.9% a year earlier.

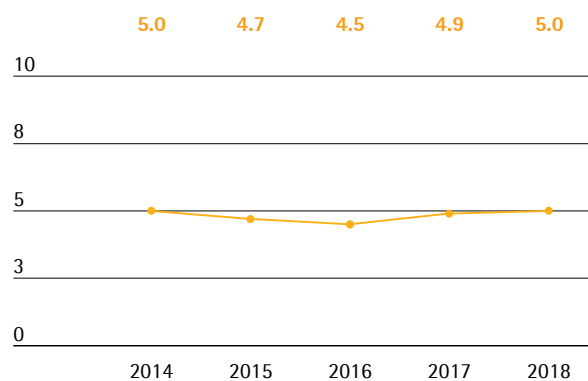
Research & Development Costs

€ in millions



Research & Development Ratio

In % of sales revenue



IFRS require that certain development costs be capitalized on the balance sheet and then amortized over subsequent years. In the reporting year, these development investments amounted to €22.8 million compared to €20.8 million the year before. This amounts to a share of 27.4% (2017: 28.1%) of the Group's total R&D expenses. Regular depreciation related to capitalized development costs totaled €4.3 million in the reporting period (2017: €4.7 million). These expenses are disclosed in the cost of sales.

To protect our know-how, we pursue a targeted intellectual and industrial property rights policy. We systematically monitor compliance with these rights and review from a cost|benefit viewpoint whether it is necessary to continue to maintain individual rights.

The number of applications for intellectual property rights filed in 2018 amounted to 125 compared to 135 in the previous year. As a result of the applications submitted in the past years, we were issued 154 patents and trademarks (previous year: 197). As of the balance sheet date, we had a total of 2,245 patents and trademarks in our portfolio (previous year: 2,073).

| | 2018 | 2017 |
|---|------|------|
| Number of patent and trademark applications | 125 | 135 |
| Registered patents and trademarks | 154 | 197 |

Capital Expenditures

The Sartorius Stedim Biotech Group increased capital expenditures considerably from €136.7 million in 2017 to €177.0 million in the reporting year. The ratio of capital expenditures to sales revenue was 14.6% (previous year: 12.6%), within the range of our forecast.

Owing to its strong organic growth, the company made significant investments during the reporting year in its production capacities. For instance, we build significant additional capacity for filters and bags at our facility in Yauco (Puerto Rico). The company is also making investments in additional capacity at our headquarter in Aubagne (France).

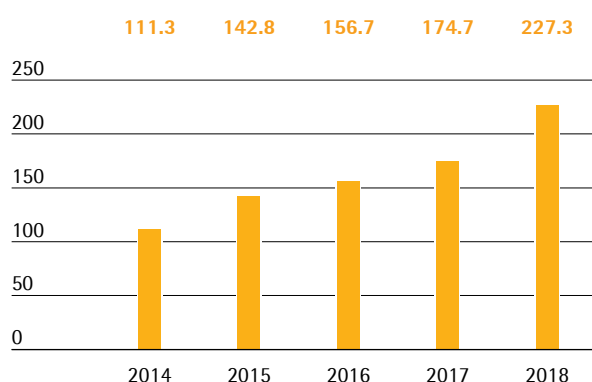
Net Worth and Financial Position

Cash Flow

In the reporting year, Sartorius Stedim Biotech again increased its net cash flow from operating activities. This figure amounted to €227.3 million relative to €174.7 million a year ago, which equates to growth of 30.1% and essentially reflects the improvement in EBITDA.

Net Cash Flow from Operating Activities

€ in millions



Net cash outflows from investing activities rose by 39.2% to €176.5 million. This increase mainly reflects investments for the expansion of our Yauco plant for single-use bags and filters as well as additional membrane casting capacities at the Göttingen site. Thus, the Sartorius Stedim Biotech Group financed its operational investments entirely from operating cash flows. The ratio of capital expenditures relative to sales in 2018 was 14.6% (previous year: 12.6%).

As we did not make any acquisitions in the reporting year, net cash flow from investing activities and acquisitions | divestitures likewise stood at –€176.5 million. The prior-year figure of –€194.9 million included acquisition-related expenses of €68.1 million from the purchase of Umetrics.

Net cash flow from financing activities, which includes payment of dividends of €43.2 million for fiscal 2017, totaled –€59.6 million. This compares to a cash inflow of €16.6 million in the previous year, which essentially reflected financing of above-mentioned acquisition.

Cash Flow Statement Summary

| € in millions | 2018 | 2017 |
|--|--------|--------|
| Net cash flow from operating activities | 227.3 | 174.7 |
| Net cash flow from investing activities and acquisitions | –176.5 | –194.9 |
| Net cash flow from financing activities | –59.6 | 16.6 |
| Cash and cash equivalents | 24.0 | 32.6 |
| Gross debt | 149.6 | 159.7 |
| Net debt | 125.7 | 127.1 |

Consolidated Statement of Financial Position

The balance sheet total of the Sartorius Stedim Biotech Group increased by €167.6 million to €1,571.5 million between year-end 2017 and the reporting date on December 31, 2018.

This increase is predominantly attributable to higher carrying amounts of property, plant and equipment and growth-driven higher inventories and trade receivables.

Non-current assets rose from €913.1 million in 2017 to €1,018.9 million in 2018, primarily due to investments in our production capacities.

Current assets amounted to €552.5 million compared to €490.8 million in the prior year. This increase was mainly driven by the aforementioned buildup in working capital.

Key Working Capital Figures
 in days

| | | 2018 | 2017 |
|-------------------------------------|-------|------|------|
| Days inventories outstanding | | | |
| Inventories | x 360 | 75 | 62 |
| Sales revenue | | | |
| Days sales outstanding | | | |
| Trade receivables | x 360 | 65 | 70 |
| Sales revenue | | | |
| Days payables outstanding | x 360 | 46 | 39 |
| Trade payables | | | |
| Sales revenue | | | |
| Net working capital days | | | |
| Net working capital ¹⁾ | x 360 | 94 | 93 |
| Sales revenue | | | |

¹⁾ Sum of inventories and trade receivables less the trade payables

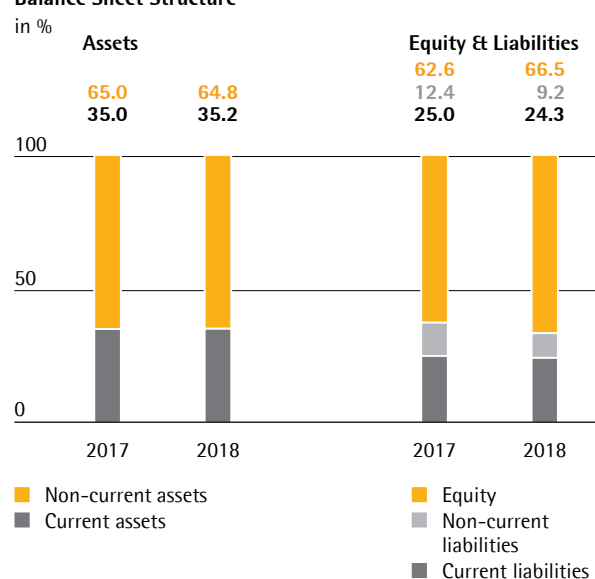
Driven by strong earnings, equity of the Sartorius Stedim Biotech Group grew from €879.5million in 2017 to €1,044.9million in 2018. Its equity ratio improved to 66.5% (December 31, 2017: 62.6%) and thus continued to remain at a comfortable level.

Current and non-current liabilities were largely unchanged at €526.6million, after €524.5million in the previous year.

Overall, gross debt decreased to €149.6million as of December 31, 2018, compared with €159.7million for the year ended December 31, 2017. Net debt as of the reporting date was at €125.7million relative to €127.1million a year ago. This figure excludes the liability for the remaining purchase price for acquisitions amounting to €8.7million in 2018.

Calculation of net debt

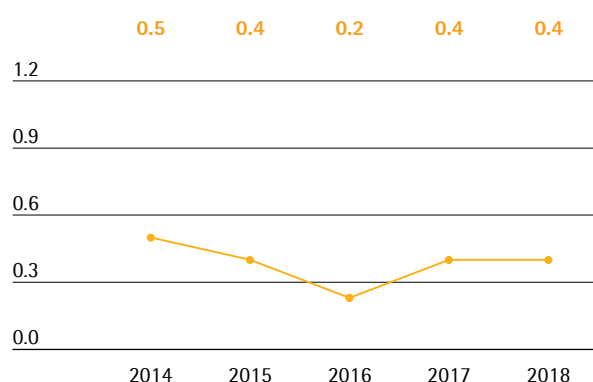
| € in millions | 2018 | 2017 |
|---------------------------|-------|-------|
| Non-current | | |
| Loans and borrowings | 43.1 | 46.3 |
| Finance lease liabilities | 15.0 | 15.8 |
| Current | | |
| Loans and borrowings | 89.8 | 95.9 |
| Finance lease liabilities | 1.7 | 1.7 |
| Gross debt | 149.6 | 159.7 |
| Cash and cash equivalents | 24.0 | 32.6 |
| Net debt | 125.7 | 127.1 |

Balance Sheet Structure


Net Debt to Underlying EBITDA

Regarding the debt financing potential of the Sartorius Stedim Biotech Group, the ratio of net debt to underlying EBITDA represents a key management indicator. This ratio stayed at the previous year's level of 0.4, in line with our expectations.

Ratio of Net Debt¹⁾ to Underlying EBITDA



¹⁾ The net debt excludes the liability for the remaining purchase price for acquisitions;
 2018: €8.7 million, 2017: €46.5 million, 2016: €49.6 million,
 2015: €47.5 million, 2014: €42.8 million

As of December 31, 2018, the total volume of all available capital and guaranteed credit lines amounted to €340 million. Of this amount, Sartorius Stedim Biotech has drawn on €117.5 million, leaving available credit of €221.5 million at the end of 2018. All in all, this ensures that all Group companies have sufficient funds to successfully finance their business operations and new capital expenditures.

We are utilizing hedging transactions to counteract the fluctuations in foreign-exchange rates to which the Group is exposed on account of its worldwide business operations. At the end of 2018, foreign-exchange contracts amounted to €145 million on a reported basis, with a market value of €1.5 million

Financing | Treasury

Our financing strategy aims to ensure our solvency at all times, limit risks associated with financing instruments and optimize our cost of capital. Sartorius Stedim Biotech covers its financing needs through a combination of operating cash flows and the assumption of short-, medium- and long-term financial liabilities.

In December 2014 Sartorius AG has entered into a syndicated revolving credit line agreement of €400 million with a maturity that was extended until December 2021. Since then, Sartorius Stedim Biotech has been utilizing a credit line with a volume of up to €310 million provided by Sartorius AG.

Furthermore, the Group has a long-term loan agreement with the Kreditanstalt für Wiederaufbau (KfW) for a current volume of €6.3 million relating to investments in production capacities and diverse bilateral credit lines of approximately €23 million in total.

The above-mentioned financing comprises instruments with both fixed and variable interest.

Product and Sales

Sartorius Stedim Biotech sells products and services for the entire added value chain in biopharmaceutical production and upstream process development. The portfolio includes cell lines, cellculture media, bioreactors, a wide range of products for separation, purification and concentration, and systems for the storage and transport of intermediate and finished biological products.

New Products with a Focus on Automation

As an industry leader in fermentation, during the reporting year Sartorius Stedim Biotech introduced a fully automated bioreactor system based on single-use technologies. It can be used to portray the development of bioprocesses even on an especially small scale. The ambr 250 system allows customers to execute up to 24 fermentation processes simultaneously during the development and optimization of processes.

In 2018, we launched an automation platform for the process control of single-use bioreactors and various filtration systems, which we developed with Siemens. With this offer, we extend our expertise as a partner for our customers' automation solutions.

In the year under review, Sartorius Stedim Biotech and the U.S. bioprocess company Repligen reached an agreement that helps customers more quickly establish single-use solutions for continuous bioprocessing. At the core of the collaboration is the idea that through continuous processing, the production of biopharmaceuticals can be intensified and system output increased in order to reduce manufacturing costs.

Modified Relationship with Lonza for Cell Culture Media Business

Since late 2012, Sartorius Stedim Biotech has been working on cell culture media with the life science company Lonza. The contract then signed by the two companies granted Sartorius Stedim Biotech the exclusive sales and marketing rights for particular cell culture media and buffers developed and manufactured by Lonza that are used in biopharmaceutical production processes. During the reporting year, Sartorius Stedim Biotech and Lonza modified the contract by mutual accord. Starting in 2019, Sartorius Stedim Biotech will retain current and future Lonza media and buffers in its portfolio of products for cell-based development and manufacture; mutual exclusivity, however, no longer applies. The

new agreement offers both contractual partners greater leeway with regard to advancements and strategic positioning.

Sales Activities Expanded

Sartorius Stedim Biotech markets its product portfolio directly through its own field sales representatives. Sales activities for key accounts are coordinated and supported by global key account management. During the reporting year, we expanded key account management and added more customers to the corresponding programs.

As part of our strategy to position our product portfolio as a platform, we entered into various agreements during the reporting year. The life science group Abzena plc selected Sartorius Stedim Biotech as its preferred partner for equipping its integrated contract development and manufacturing organization (CDMO) systems in Bristol, Pennsylvania and San Diego, California. ABL Europe, a subsidiary of the U.S. contract production and laboratory research company ABL Inc., chose Sartorius Stedim Biotech as its primary supplier of single-use systems.

Production and Supply Chain Management

Sartorius Stedim Biotech operates a very well developed production network around the world. The largest production sites are located in Germany, France and Puerto Rico. Sartorius Stedim Biotech also manufactures in the United Kingdom, Switzerland, Tunisia, India, the United States and, since 2018, in China.

Expansion of Production Capacity

A new manufacturing facility for filters opened in Göttingen during the reporting year. Sartorius Stedim Biotech also expanded production capacity at its Yauco site in Puerto Rico, where membrane filters and sterile bags are manufactured primarily for the U.S. market. For products manufactured in Yauco, the company also began operating a warehouse in Florida, USA, which helps ensure smooth supply to customers on the U.S. mainland.

A new logistics center, which opened in Aubagne in 2018 and covers 12,000 square meters, provides the capacity that the company will need in the years to come. An additional cleanroom in Aubagne also boosts the site's production output.

The topping-off ceremony for the development and production center for cell lines was held in Ulm, Germany, in the year under review. With 6,000 square meters of usable floor space, the new center will be about twice as large as the current facility in Laupheim, near Ulm. The company plans to move operations from Laupheim to Ulm in late 2019.

In addition, Sartorius Stedim Biotech has begun producing sterile single-use bags in Beijing (China), where it plans to manufacture additional products.

Sustainability

Sustainability information for the Sartorius Stedim Biotech Group is not reported. In accordance with the provisions of Article L.225-102-1 IV of the French commercial code, Sartorius Stedim Biotech is exempted from presenting this information, because it is included in the non-financial statement established and published by the controlling company, Sartorius AG, as per applicable German regulations.

Opportunity and Risk Report

Principles

Every business activity entails opportunities and risks, which have to be managed. The skill with which this is done goes a long way in determining the future development of a company's shareholder value.

It is not the task of risk management to eliminate all risks: rather, our approach is to intentionally take a certain measure of risk in our business activities in order to be successful in unlocking opportunities. However, in this endeavor, it is important to keep risks contained within acceptable limits and to control them carefully. Through appropriate guidelines, we ensure that risk assessments are taken into account in the decision-making processes from the very beginning.

Sartorius Stedim Biotech has decided to make the identification and the management of risks and opportunities a cross-functional component of Group management. In this context, Sartorius Stedim Biotech's risk management is integrated into the Sartorius Group organization. Our risk management organization reflects a global functional matrix organization in which individuals heading a functional area are each responsible for their own management of opportunities and risks. The Finance & Controlling department is responsible for the organization of the respective reporting process, including the further development of the Group's risk management system.

Managing Opportunities

Our opportunity management centers on the analysis of target markets and sector environments, as well as the assessment of trends, both of which give strong indications as to future business opportunities. The identification of the potential for development in this context is one of the key roles of the relevant managers and initially takes place at the local rather than the central level. The market-facing functions, such as marketing and product management in the individual divisions, play a leading role in this respect. The central Business Development unit supports these areas with market monitoring, data analysis and the implementation of strategic projects.

As part of strategy reviews, the members of the Group Executive Committee regularly meet with the managers having operational responsibility to discuss short-, medium- and long-term opportunity potential for the various business areas. The subsequent steps of prioritizing the opportunities and evaluating them

from a business management perspective, deriving strategic measures and allocating resources proceed in accordance with a standardized decision-making process that applies throughout the Group. If the opportunities are short-term in nature, they are considered in annual budget planning. Medium- and longer-term opportunities are tracked systematically as part of strategic planning.

As a supplier for the pharmaceutical industry, Sartorius Stedim Biotech operates in a future-oriented and high-growth sector. The significant opportunities generated by the various market and technology trends are described in detail in the sections entitled "Sector Conditions" and "Outlook for the Sector" on pages 22 et seq. and pages 47 et seq., respectively.

Our assessments rank the company as one of the global market leaders in many subsegments and product areas. We believe the high quality of our products, our strong brand recognition and our established customer relationships give Sartorius Stedim Biotech strong opportunities to continue extending our market leadership. The corresponding strategies and the growth opportunities and initiatives based on them are discussed in the section on the strategy of the Group, which begins on page 19.

Strict management of processes and costs provides opportunities to further increase our profitability. Key target areas in this respect include continued enhancements of our procurement chain and ongoing efforts to optimize production, which we present on page 32.

Risk Management

Organization

The overall responsibility for the maintenance of an effective risk management system ensuring comprehensive and consistent management of all material risks rests with the Audit Committee. The Finance & Controlling Department is responsible for coordinating and developing this system and for consolidated risk reporting, while the particular functional areas are responsible for identifying, analyzing and reporting individual risks. This includes the assessment of their potential impact and the decision on taking the appropriate countermeasures.

The Audit Committee monitors the effectiveness of the risk management system. Furthermore, while carrying out their statutory audit mandate for the annual financial statements and consolidated financial statements, the independent auditors examine whether the early warning system in place is capable of prompt identification of risks that could jeopardize the future of the company. Finally, the Internal Audit Department regularly reviews the risk management process and system.

Insurance

We have taken out insurance policies to cover a wide range of risks where possible and economically advisable. These insurance policies include coverage against product liability, property damage, business interruption, transport, material and pecuniary damages and other risks, and provide comprehensive coverage for legal costs. An independent department working in conjunction with an external insurance broker regularly reviews the nature and extent of our insurance protection and makes any adjustments as necessary.

When choosing our insurers, we particularly consider the credit rating of these entities as potential contractual partners, as well as aim to achieve a high degree of diversity in order to mitigate the related risks.

Risk Management System and Risk Reporting

Sartorius has implemented a global guideline (Risk Management Handbook), which includes definitions of the framework, the structural organization, processes, risk reporting and monitoring and controls of the effectiveness of the risk management system. The handbook is based on the internationally recognized COSO (Committee of Sponsoring Organizations of the Treadway Commission) standard. There are also a number of other sources that contain stipulations for handling risks, including the articles of association and rules of procedure of the Group companies and other internal guidelines.

The prescribed reporting process in the risk categories subsequently described establishes the rules for the ongoing review of and information on risk situations. If any specific risks are discernible, these are documented with respect to their assessment, probability of occurrence and measures to be taken to eliminate such risks or to mitigate their impact.

We have an urgent reporting procedure in place to ensure that when a new or emerging significant risk to our net worth, financial position and profitability is identified, the audit committee receives all of the necessary details without undue delay.

To classify risks appropriately, we have defined four main categories: external risks, operating risks, financial risks and compliance risks. Each main category is divided into several subcategories that are described in the following sections.

| Probability of Occurrence | |
|---------------------------|-----------|
| Remote | < 10% |
| Possible | 10% - 50% |
| Probable | 50% - 90% |
| Very likely | > 90% |

| Significance | |
|------------------|--------------------|
| in millions of € | Impact on Earnings |
| Insignificant | < 10 |
| Moderate | 10 - 50 |
| Significant | 50 - 100 |
| Critical | > 100 |

Explanation of Principal Risks and Opportunities

External Risks

General Risks

In principle, our ability to foresee and mitigate the direct and indirect effects of risks entailed by life in general is limited, but we proactively take measures, whenever feasible, to ensure that we can respond appropriately and at short notice or are insured against any damage entailed by such risks that include, for instance, natural catastrophes and their associated damage to commercially significant and critical infrastructure.

Our largest sites in Germany and France do not face major risks from natural catastrophes, while e.g. our production plant in Puerto Rico is exposed to the risk of severe hurricanes. We control this risk by applying high security standards to the buildings and explicitly consider this risk in our warehousing and production network strategy.

Political developments, such as the referendum on the United Kingdom's leaving the European Union ("Brexit")

or the change in politics in the USA, can have an impact on the Group's business. Such developments may involve changes to the tax system or customs duties, as well as impacts on the exchange rate of the euro to the British pound or the U.S. dollar (for more on the subject of exchange rates, see the section below on Exchange Rate Risks).

In the U.K., we run various manufacturing and sales entities with a significant business volume. Any development that has a negative impact on the trading between the U.K. and other countries could therefore lead to a corresponding decrease in Group's earnings. The further developments are being closely observed so that measures can be taken to reduce such risks, as necessary. The tax reform implemented in the United States in 2017 has so far reduced the Group's tax payments; however, other measures might result in a tax burden.

Business Cycle Risks

The nature of our various business areas means that Sartorius Stedim Biotech as a whole is insulated to a certain extent from the full force of wider cyclical effects. If economic developments prove more positive than expected, this, in turn, can additionally stimulate stronger growth.

Operational Risks and Opportunities

Our supply chain extends from procurement to production to sales and distribution. Problems within this workflow can have consequential effects, including delays in deliveries. The global supply chain management system we have instituted throughout our production processes to prevent such problems largely minimizes the associated risks by analyzing and controlling all of the operations involved. The strongly international alignment of our organization opens up a whole series of opportunities too. The various risks and opportunities encountered within our supply chain are explained in detail below.

Procurement Risks and Opportunities

We purchase a wide range of raw materials, components, parts and services from suppliers and are consequently exposed to the risks of unexpected delivery bottlenecks and/or price increases.

Over the past years, we have implemented powerful tools and robust processes in our Materials Management unit to manage risks and critical materials.

These means enable us to meet the needs of our customers with respect to delivery reliability and transparency. This can represent a competitive advantage. Important measures in this respect are to maintain security stock and to define alternative suppliers according to our second-supplier policy. We moreover conduct regular supplier reviews and also use early warning systems.

Risks from raw material prices play a rather subordinate role in our business. On the one hand, the proportion of raw materials in our production costs is comparatively low. On the other hand, we purchase a wide range of materials so that price increases for certain materials do not represent any significant impacts.

Opportunities can arise in the area of procurement when our growth enables us to increase order quantities and thereby strengthen our position with our suppliers, such as by receiving price discounts or preferential treatment as a "preferred customer." In addition, we maintain a list of preferred suppliers in order to enter into long-term business relationships with key suppliers to our mutual benefit.

Increased globalization of our supplier pool holds the prospect of purchasing on more favorable terms, moreover, and there is also a possibility of our expanded purchasing activities in the international markets leading us to identify suppliers with special product and technical expertise that could eventually enhance our own competitive edge.

Production Risks and Opportunities

Based on our core technology expertise, we ourselves manufacture a large proportion of the products that involve a high level of vertical integration. Other products, such as reusable fermenters and bioreactors, are manufactured in collaboration with suppliers so that some of the production risks are transferred to external third parties. When we manufacture products ourselves, we also bear the associated risks of capacity bottlenecks or overcapacity, production downtimes, excessive reject rates and high levels of tied-up working capital, as well as dependency on individual manufacturing sites.

We contain and reduce these risks by planning production capacities carefully, using versatile machines, semi-automated individual workstations and flextime work schedules, and by continuously monitoring production processes. Moreover, our global manufacturing network enables us to compensate for any capacity bottlenecks by shifting production to other

regional plants and to minimize our dependency on individual local production plants. Furthermore, we have taken out policies for business interruption insurance to compensate for any possible losses due to production downtimes.

We consider it an opportunity that our investments in infrastructure and production resources, among other things, have given us high flexibility in our manufacturing operations and that we are capable of meeting our customers' requirements and regulatory standards with respect to business continuity concepts. In addition, this approach ensures that our individual production sites can concentrate on specific manufacturing technologies, gaining added efficiency as a result. Our international manufacturing network also makes it possible to capitalize on the cost advantages offered by individual sites. Furthermore, continuous improvements in production, such as simplifying processes and increasing levels of automation, help drive manufacturing efficiency even higher.

Sales and Distribution Risks and Opportunities

We use a variety of channels to sell and distribute our products around the world. The potential risks entailed are unexpected changes in the demand structure, growing price pressure and non-compliance with supply agreements concluded with customers. We employ targeted market analyses to identify emerging demand trends in individual segments early on so that we have time to respond appropriately. Our technical innovations and the fact that a wide range of our products is used in validated production processes in the biopharmaceutical industry, reduce our exposure to the risk of growing price pressure. We have reduced our risk exposure in the area of logistics in recent years by setting up and using central warehouses to optimize distribution logistics.

Opportunities arise in the area of sales and distribution when the increasing breadth of our product range puts us in a position to sell new products to existing customers. Our business relationships, most of which are established for the long term, and our global presence provide further opportunities.

Sartorius Stedim Biotech sources its key customers from the pharmaceutical, chemical and food industries. These customers are usually relatively large organizations that have been in existence for some time and have strong credit ratings. Accordingly, the Group has had low to zero credit losses over the past years, and its overall credit risk continues to be at a

very low level. Most of our business areas have a highly diversified customer base, so the Group as a whole is not dependent on individual key accounts to any significant degree.

Competitive Risks and Opportunities

Sartorius Stedim Biotech has a leading competitive position in most of its markets. Some of our competitors are larger than us, and most share our status as a globally operating company. As we serve a large number of customers from highly regulated sectors like the pharmaceutical and food industries, and the technology barriers to market entry are substantially high, we regard the probability of new competitors emerging within the short term as low.

The fact that many of our products are used in validated bioprocesses reduces the risk of losing significant market share within a short timeframe. At the same time, it is also more difficult for us to quickly force out the competition that serves customers in this area.

Changes in the competitive environment, for example, a further consolidation in the markets, can pose opportunities. We have been continuously making acquisitions in recent years to reinforce our market position and open up new potential synergies.

Quality Risks and Opportunities

Our customers use Sartorius Stedim Biotech products in a wide range of critical production processes, including the manufacture of pharmaceuticals, foods and chemicals, and in research and development laboratories. The main risk encountered in these areas is non-compliance with agreed quality criteria, which can lead to losses for our customers, or their customers, for which we may be made liable through compensation claims.

We employ rigorous quality checks and advanced production methods and processes, such as cleanroom technology, to ensure that our products satisfy the most stringent quality standards and high regulatory requirements. These manufacturing methods and processes are subject to constant review under our continuous improvement processes, moreover, and are optimized as requirements evolve. Our successful completion of a host of annual audits by customers and implementation of quality systems compliant with ISO 9001 and, where applicable, with ISO 13485 document the high level of quality achieved in Sartorius products and processes. Irrespective of these

measures, we also maintain significant insurance coverage against product liability risks. Sartorius Stedim Biotech has established a traceability system that enables us to recall an entire production batch immediately, if necessary, and minimize any adverse consequences in the event of defects being discovered in a product.

We have installed a complaints management system to deal with customer requests and to ensure full documentation.

In the sectors we address, quality requirements are growing more and more stringent all the time, not least as a result of increasing requirements on protection of medical patients and on product safety by regulatory authorities. Increasing and changing requirements typically entail the risk that a new requirement might be overlooked or be difficult to achieve, but we regard this first and foremost as an opportunity that opens up new market prospects. The reason is that challenging quality demands represent a considerable barrier to entry for potential new competitors and provide stimulus for further technical innovation to which we actively respond. Moreover, we actively seek to draw up new requirements through our work on professional committees, membership in industry associations and standards committees, and are able to identify emerging requirements these at an early stage and prepare ourselves accordingly.

R&D Risks and Opportunities

We devote a considerable share of our resources to research and development. Potential risks in this area may arise from development results that diverge from market needs and application requirements and from exceeding planned development deadlines. Our advanced project management, intensive R&D controlling and early involvement of our customers in the development process substantially limit these R&D risks. We ensure that product developments are always reviewed very promptly with regard how well they meet the customers' needs so products can be adapted accordingly as needed. Patents and continuous tracking of the technologies and competitors relevant to us secure our technology and marketing position.

On the other hand, the R&D sphere also offers a number of potential opportunities. Our intensive collaboration with partners that rank among the global market leaders in their own fields opens up the opportunity for us to jointly develop products with an especially high level of innovation. In areas such as membrane technology and plastics technology, as well as sensorics and bioprocess engineering, in turn, the

expertise of our own specialists puts us at the very forefront of global research and development, presenting us with an opportunity to turn this technical knowledge into potential sales and an even stronger position on the market. The combination of different innovative activities in a separate Corporate Research Department further enables us to identify and benefit from promising developments and emerging trends at universities, startups and at our customers' plants.

Acquisition Risks and Opportunities

By nature, acquisitions provide many opportunities, such as sales growth, extension of our product portfolio and development of new markets. By contrast, the purchase and sale of companies or parts of companies entail a number of typical risks, such as incorrect valuation assumptions or insufficient usage of anticipated synergy effects. To prevent these risks, we take various measures, such as performing a standard due diligence review of important areas and carrying out comprehensive analysis of the market concerned. In addition, we involve external consultants and experts in the purchase or sales process as required. We especially focus on drafting transaction contracts so that they adequately counter such risks, especially by clauses assuring specific characteristics or by contractual warranty or guarantee provisions, as well as agreements on mechanisms for adjustment of the purchase price and on liability clauses. Immediately after an acquisition has taken place, an integration phase is initiated in which any potential risks can likewise be detected as early as possible and prevented or minimized by taking the appropriate counteractions.

Personnel Risks and Opportunities

As an innovative technology group, Sartorius Stedim Biotech employs a large percentage of highly qualified people. We counter the risks of a possible scarcity of required specialists, especially those in key positions and of demographic change by offering performance-related remuneration models, targeted continuing professional development options, further attractive social benefits, continuous education and training for junior staff members within our organization and interesting people development opportunities.

The success of these measures is apparent in the low attrition rates of recent years and the many years of seniority our people accumulate on average. Employment contracts in certain cases contain a clause prohibiting any move to a direct competitor.

Opportunities for Sartorius Stedim Biotech primarily arise in that it can further qualify its staff by offering its own training courses and retain such staff over the long term, thus covering company needs for qualified personnel particularly well.

IT Risks and Opportunities

The business processes of the Sartorius Group are supported by IT applications and systems. Failure or

other impairment of the relevant IT systems or (cyber)attacks can considerably disrupt the smooth functioning of the companies business processes and lead to manipulation or to uncontrolled loss or leakage of knowledge or data.

We minimize this risk by continuously investing in the setup and operation of secure IT systems and applications and by continuously further developing and implementing our concepts and security measures based on the International Standard ISO 27001, Information Security Management System. In addition, we incorporate the results of regular audits and vulnerability assessments carried out by external companies specializing in IT security.

Protection of our data against misuse is ensured by specific authorization and authentication policies based on the assignment of rights limited to a "need-to-know" basis for performing certain tasks, and the application of such policies is reviewed at regular intervals.

We protect our systems against failure and data loss by regular data backups, recovery testing based on rolling disaster scenarios and risk-based use of redundant IT infrastructures. Multi-factor authentication solutions enable us to prevent malware threats.

We are convinced that the threat of cyberattacks is growing worldwide, both in number and intensity. This is why we are continuously extending and strengthening our activities: We are improving our efforts by further automating management of authorizations and reducing the potential for data misuse, among other measures. We inform our staff in a targeted way about possible threats and risks, involving our employees by providing them with simple but effective options for decentralized defense and reporting suspicious emails to IT for checking.

By extending our means for competent and fast response to cyberattacks including other IT security incidents, we supplement our organizational basis for running the Sartorius system and applications at the lowest possible risk across the entire landscape.

Financial Risks and Opportunities

The global nature of the Sartorius Stedim Biotech Group's operations means that its business activities are inevitably exposed to financial risks. The most significant of these, are exchange rate risks, interest rate risks and liquidity risks, all of which are described below and addressed in detail in the Notes to the

Consolidated Financial Statements. Vice versa, financial risks, most notably exchange rate risks and interest rate risks, are balanced by opportunities of approximately equal magnitude.

Exchange Rate Risks

As a consequence of its global business activities, the Sartorius Group is exposed to foreign currency fluctuations. Since we generate around two-thirds of consolidated sales revenue in foreign currencies and, of this figure, approximately two-thirds of this total revenue in U.S. dollars or in currencies pegged to the U.S. dollar, we are positively or negatively impacted by currency effects, especially when converting the currencies of balance sheet items and profit or loss items, respectively. Besides the U.S. dollar, other key currencies are the South Korean won, the Chinese renminbi, the Swiss franc and the British pound.

Our global production network thus enables us to offset the lion's share of sales revenues received in foreign currency within the Group against costs likewise incurred in foreign currency. For example, we manufacture many of our products for the North American market locally, and are not disadvantaged in competition with our U.S. rivals, insofar as this general currency risk is concerned. We continuously monitor the portion of our foreign currency sales revenue that remains after we have settled our costs, so-called net exposure

In order to evaluate and steer the remaining risk based on the expected net exposure for the next 12 months and take into consideration hedging transactions already executed, we are continuously calculating our risk exposure with a cash flow at-risk model. We use this basis to decide on whether to use additional derivative financial instruments, especially spot, forward and swap transactions, to adjust for maximum loss. Hedging transactions are set up by one group of staff and monitored by another, separate group.

Interest Rate Risks and Opportunities

We have concluded fixed interest agreements for approximately two-thirds of our outstanding loans to eliminate the risk posed by variable interest payments. The remaining portion of the financial instruments outstanding on the reporting date is subject to variable interest based on the market rate. We monitor interest rate trends and our interest rate exposure constantly and have the facility to arrange for hedging transactions where we consider it necessary and economically advisable to do so for individual loans. As of December 31, 2018, we did not have any interest

rate derivatives in our portfolio of financial instruments.

Liquidity Risks and Opportunities

The Sartorius Stedim Biotech Group's liquidity is managed centrally in order to minimize liquidity risks and optimize liquidity allocation within the organization. For this purpose, various long- and short-term financial instruments are utilized. Regarding the maturities of our loans, we make it a policy to take a risk-averse approach.

In addition to a 300 MEUR credit line provided by Sartorius AG that can be accessed and repaid at short notice, we have a number of bilateral working capital credit lines for individual Group companies in place. Furthermore, we use cash pooling agreements between selected Group companies as the primary tool to manage liquidity within the Group.

Compliance Risks

Regulatory Risks

Our role as a supplier to the biopharmaceutical industry and health care providers means that Sartorius Stedim Biotech can also be affected by underlying developments in these areas. The possibility of the regulatory authorities (FDA, EMA) adopting a more restrictive approach to the approval of new medications remains the principal source of risk in this context. Such a move would reduce the number of new pharmaceutical products to be marketed and would consequently downgrade future prospects for Sartorius Stedim Biotech over the medium term.

Environmental Risks

Sartorius Stedim Biotech has established an environmental management system that encompasses, and is integrated into, all divisions and covers a whole series of environmentally relevant regulations to minimize environmental risks. This management system has been certified for compliance with ISO 14001 at a number of the company's relatively large manufacturing sites. The respective company organizational units ensure at the particular sites that the laws and regulations relating to environmental protection are observed and that further technical possibilities for limiting environmental risks are identified on an ongoing basis.

The increasing importance of sustainability considerations in many industries represents an opportunity. That is why this aspect is a key element in our supplier selection process for assessing the suitability of a particular company as a business partner.

Litigation Risks

Litigation risks for Sartorius Stedim Biotech can arise from pending or forthcoming legal disputes or from administrative proceedings. All judicial or extrajudicial

disputes are attended to by the company's own attorneys and legal experts, who engage external lawyers as needed.

At present, there are no pending or discernible legal disputes or proceedings that lack any cost coverage allowances in the statement of financial position or that could have a substantial negative impact on Group.

Assessment of the Overall Risk Situation and Risk Outlook

Where feasible, we adopted countermeasures and/or arranged for balance sheet measures during the reporting year to cover all discernible risks within the Sartorius Stedim Biotech Group, and those of a defined probability of occurrence, that had the potential to damage our net worth, financial situation and profitability.

For the purposes of this report, we have assessed the probability of occurrence of the risks shown below and, in the adjacent columns, classify their particular significance for the entire Group.

| Risk Category | Probability of Occurrence | Significance |
|--------------------------------|---------------------------|---------------|
| External risks | | |
| General risks | Possible | Significant |
| Business cycle risks | Possible | Moderate |
| Operating risks | | |
| Procurement risks | Remote | Significant |
| Production risks | Remote | Significant |
| Sales and distribution risks | Possible | Significant |
| Competitive risks | Remote | Moderate |
| Quality risks | Remote | Significant |
| Research and development risks | Remote | Significant |
| Acquisition risks | Possible | Significant |
| Personnel risks | Remote | Moderate |
| IT risks | Possible | Significant |
| Financial risks | | |
| Exchange rate risks | Probable | Moderate |
| Interest rate risks | Probable | Insignificant |
| Liquidity risks | Remote | Moderate |
| Compliance risks | | |
| Regulatory risks | Possible | Significant |
| Environmental risks | Remote | Moderate |
| Litigation risks | Possible | Moderate |

After thorough analysis of the entire risk situation and according to our current review, there are no discernible risks at present that could jeopardize the continued existence of the Group.

Similarly, based on our current review, there are no discernible risks that could jeopardize the future existence of the Group.

Material Events after the Reporting Date

No material events, of any nature, occurred after the reporting date.

Internal Control Procedures

Introduction

The objectives defined by the Chairman for the internal control system of Sartorius Stedim Biotech are as follows:

- Prevent risks that would endanger the quality of the assets of Sartorius Stedim Biotech or even its existence;
- Ensure that the executive management activities, the transactions completed and the conduct of employees comply with the guidelines defined by executive management, applicable laws and regulations, the fundamental values, standards and internal rules of the business and the ethical codes and conventions of the healthcare industry;
- Ensure that accounting and financial information and management data provided to the executive management of the company accurately reflect the operations of Sartorius Stedim Biotech;
- Prevent risks arising from operations, errors or fraud, especially in the accounting and financial area.

Scope of Internal Control

The internal control system described covers the parent company and its affiliates.

Components of Internal Control

Environment for Internal Control

The core of any business is its people (their individual attributes, including integrity, ethical values and expertise) and the environment in which they operate. They are the engine that drives the organization and the foundation that supports the company.

Risk Assessment Process – Risk Mapping

The company must be aware of, and deal with, the risks it faces. It must set itself objectives and integrate them into its sales, production, marketing, financial and other activities so that the organization operates in concert. It must also establish mechanisms to identify, analyze and manage the related risks.

Control Activities

These control activities are undertaken at every level of the Group to ensure that internal control is efficient: checking the accuracy, completeness, authorization, validation and recording of transactions and ensuring that different people discharge different duties so as to reduce the risk of errors or fraud.

Information and Communication

The availability of accurate, reliable and complete information is essential both to achieve business objectives and to enable proper reporting to all parties concerned in compliance with the applicable laws and regulations.

Monitoring, Control and Management

Responsibilities and authorities must be defined and understood at all levels of a company for internal control to function effectively. Duties must be assigned in such a way that a person's work is always checked and approved by a different person. Where the size of the local unit concerned permits, responsibility for initiating, authorizing, recording and processing transactions must always be assigned to different individuals.

Unit management is responsible for maintaining internal checks and internal control at all times.

Internal Controlling Roles

Executive Management

The Chairman and Chief Executive Officer is responsible for the internal control system and management at all levels. He is also responsible for the development, operation, monitoring and management of the internal control and controlling systems and for providing the necessary assurances that these steps have been implemented.

Audit Committee

The Audit Committee is responsible for carrying out any necessary reviews and evaluations of the internal controlling procedures, including those relating to financial information, and also assists with the preparation of the Group's consolidated financial statements. For further information about the Audit Committee, see page 69.

Risk Management

The Sartorius Stedim Biotech Group is inevitably exposed to a wide variety of risks by the nature of its operations around the world. Accordingly, an internal risk management system has been set up to help identify, assess and manage these risks efficiently. Within this risk management system, an ad hoc committee comprised of representatives of different departments regularly studies current issues of risk management. This enables the committee to provide executive management with an overview of the risk to which the company is exposed, enabling it to take appropriate action when required.

Internal Auditing Department

The Internal Auditing Department is in charge of monitoring the effectiveness and suitability of risk management and the internal control system in Sartorius Stedim Biotech Group companies, as well as compliance of all activities and processes with internal and external rules and standards. It provides independent auditing and consulting services that focus primarily on compliance with all relevant legal provisions and the improvement of business processes at the company. To ensure the independence of the internal auditors, the Audit Committee receives, at least once a year, a report from the Internal Auditing Department on the work they have done (according to the audit plan established by this committee) and their findings with regard to Group affiliates.

Finance and Controlling Departments

The Finance and Controlling Departments track and monitor operations and projects to optimize the Group's profitability and cash flow, providing both internal and external stakeholders with reliable information.

These two departments define the Group's accounting rules and methods and its principle financial processes (five-year business plan, budget, etc.) as well as reporting tools, in order to monitor the day-to-day business.

Procedures for Preparing the Group Financial Statements and Other Accounting and Financial Information

The accounts of affiliates are prepared in accordance with the Group's accounting policies. The data is then adjusted, where necessary, to produce company accounts that comply with the applicable local legal and tax provisions. Integrated consolidation software is used both for management reporting purposes and to produce the Group financial statements.

Since 2013, the Group has decided to implement a hard close process as of November 30 in order to anticipate and improve the annual audit.

Accounting Standards

The consolidated financial statements are prepared in accordance with IFRS accounting standards as currently adopted by the European Union. The consolidated financial statements comply with accounting rules and methods as detailed in the Notes to the Consolidated Financial Statements.

Roles of the Group's Finance and Controlling Departments

The Finance and Controlling Departments check the quality of the reporting packages submitted by affiliates, focusing primarily on the following elements: checking corporate data and consolidated adjustments entered locally, inter-company eliminations, the accounting treatment of non-recurring transactions for the reporting period, and verifying principal movements between the opening and closing balance sheets to prepare the cash flow statement.

The Finance Department also verifies the results of procedures, including currency translation, intercompany eliminations, etc.

Key points of review include the preparation and validation of the statement of changes in shareholders' equity and the cash flow statement.

Financial Information and Reporting

The Group's rules and procedures in relation to financial reporting and accounting are set out in the Accounting and Reporting Manual. Application of and compliance with these principles, rules and procedures are the direct responsibility of the finance director of each affiliate. They must ensure that information provided via the Management Information System complies fully with all applicable disclosure requirements.

Executive Management reviews the effectiveness of the internal controlling of financial reporting regularly. In particular, it verifies that transactions have been recorded consistently, in accordance with IFRS international accounting standards as applied by the Group and as set out in the Accounting and Reporting Manual, in order to ensure the pertinence of transactions and assets recognized within the times set.

Internal Control in 2018

From an internal control perspective, the Group focused on the following this year:

Code of Conduct and Anti-Corruption Code

Employees can consult the Sartorius Code of Conduct and the Sartorius Anti-Corruption Code; initial trainings have been completed and implemented as a controlled operation.

These codes are subject to reviews and revisions as required by amended legislation. In addition, all employees of the company and of the Group are made aware of these codes and must ensure compliance with them on a daily basis.

Corporate Transactions

The company complies with the Recommendation of the Autorité des Marchés Financiers of November 3, 2010, and with the AFEP-MEDEF Code, as revised in June 2018. Thus, transactions involving the purchase or sale of the company's securities or financial instruments are prohibited during the periods between the date on which managers, persons considered managers under the law, and any person having regular or occasional access to privileged information are aware of precise information on the course of business or prospects that, if made public, could have a significant influence on the price and the date on which the information is made public.

In addition they are also prohibited for a period of:

- fifteen calendar days prior to the date of publication of the annual and semi-annual financial statements of the company,
- fifteen calendar days, preceding the date of publication of the quarterly information, this day included.

At the beginning of each calendar year, the company prepares and distributes a schedule setting out the periods during which trading in the company's securities is prohibited, specifying that the days indicated do not prejudice the existence of other closed periods resulting from the knowledge of specific information that directly or indirectly concerns the company, which, if it were made public, could have a significant influence on the share price of the company's shares.

In accordance with the recommendations of the AFEP-MEDEF Code and the Autorité des Marchés Financiers Recommendation No. 2010-07 of November 3, 2010, hedging transactions of any kind, on the company's shares with regard to stock options are prohibited.

Mid-Term Prospects

The Group will continue to work on Internal Control issues by strengthening its approach to risk mapping and risk management. This process will be based on elements of the AMF Internal Control Reference Framework. In addition, the process of defining mandatory minimum standards of internal controls applying to all Group companies has been followed by publishing the Group Internal Controls Handbook and will be pursued further in 2019.

Forecast Report

Continued High Growth in the Biopharmaceutical Industry

Strong, long-term trends drive growth in the pharmaceutical industry, which is almost entirely independent of business cycles. Market observers estimate that growth of the world's pharmaceutical market will reach between 3% and 6% per year during the period up to 2022. Experts forecast that the biopharma segment of the pharmaceutical market, which has been enjoying particularly strong growth for years, will continue to outperform the market. Forecasts anticipate average annual growth of around 8% or 9% through 2022, which would increase market volume from a current €217 billion to €300 billion. The share of biological medications and vaccines in the total revenue generated by the global pharmaceutical market is expected to continue rising.

In coming years, the most dynamic market will likely be China. Although the biopharmaceutical market there is still in its infancy, the country offers especially high growth potential owing to favorable policy conditions, an increasing number of domestic biotech companies and rising demand for advanced biopharmaceuticals. Experts also anticipate considerable growth in the United States and Europe, driven in particular by a growing need for medications for aging societies as well as the rising number of chronically ill and multi-morbid patients. In addition, more and more medications are being approved. For example, biopharmaceuticals are increasingly being used in yet-to-be fully explored therapeutic areas and in the treatment of rare diseases that have so far been untreatable. Innovative types of therapy for regenerative medicine and new substance classes, such as antibody-drug conjugates (ADCs), are increasing the

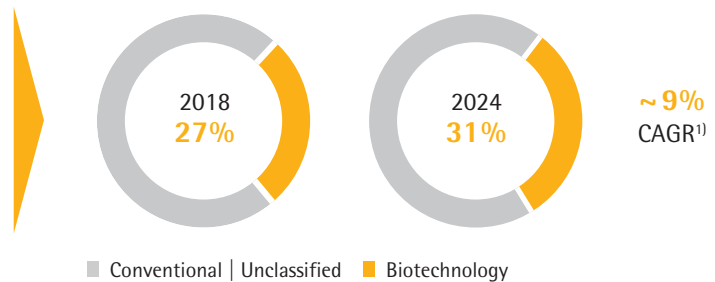
number and range of approved biopharmaceuticals as well as necessitating investments in innovative production technologies.

This relatively young biopharmaceuticals segment is driving sector growth with its high innovative power, as reflected in the strong research and development pipelines. Of the estimated 10,000+ medications in R&D pipelines, over 40% are based on biological manufacturing processes. These include more than 950 biosimilars and 550 biobetters, which are generic versions of biologic agents with better efficacy or fewer side effects than the original compounds.

Biosimilars are contributing increasingly to the growth of the biotechnology market. Governments in emerging and developing countries are fostering the creation of domestic production capacities intended to meet the rising demand for medications, which in turn encourages the founding of new biotech companies. The biosimilars market in industrialized countries is also likely to expand considerably in coming years due to the expiration of patents for high-volume biopharmaceuticals and an increasing number of approved biosimilars. In addition, the regulatory, patent-law and marketing challenges that have slowed progress in the past will likely decrease gradually. Against this backdrop, current estimates indicate that by 2022, the market could grow by an annual average of 30% and amount to a volume of around €18 billion.

Biopharma: A Growing Market

- Growing & aging population
- Increasing access to healthcare
- Strong R&D pipelines
- Emerging biosimilars market



¹⁾ Evaluate Pharma®: World Preview 2018, Outlook to 2024; June 2018; CAGR 2018 to 2024

The biopharmaceutical industry must meet a growing demand for medications while also producing an increasing number of approved medications and new types of therapy. For these reasons, industry observers expect that worldwide bioreactor capacities will continue to expand at a similar pace in the years to come. At the same time, the industry faces rising cost pressure. This increases the significance of innovations for boosting flexibility and cost efficiency in biopharmaceutical research and production. In the future, the biopharmaceutical market will shift away from a low number of especially high-volume medications that account for a majority of total production volume towards an expanding range of products for smaller groups of patients. What's more, technological progress leads to ongoing improvements in the productivity of biopharmaceutical production processes. According to the research and consulting institute BioPlan, manufacturers will therefore likely rely increasingly on single-use technologies for the commercial production of many new medications. Particularly in the case of relatively small batches, single-use technologies already ensure more cost-effective production than conventional stainless-steel units. To master these challenges, more and more pharmaceutical companies are relying on digitalization and automation as well as innovative software solutions for controlling and optimizing their processes.

of the biopharmaceutical industry and the business-friendly climate. In Europe, experts anticipate growth of 3.3% – although especially in this region, macroeconomic and political uncertainties constitute a risk. As before, market observers expect the highest growth rates in Asian countries such as China and India, in which the individual labor markets are likely to grow around 7.2% to 8.7% in 2019.

Sources: IQVIA Institute: 2018 and Beyond: Outlook and Turn-ing Points, March 2018; Evaluate Pharma: World Preview 2018, Outlook to 2024, June 2018; BioPlan: 15th Annual Report and Survey of Biopharmaceutical Manufacturing Capacity and Production, April 2018; Frost & Sullivan: 2018 Mid-year Report: Forecast and Analysis of the Global Market for Laboratory Products, May 2018; BCC Research: Biosimilars: Global Markets, March 2018

Labor Market Remains Stable while Macroeconomic Risks Increase

According to Frost & Sullivan, global demand for laboratory products is likely to continue increasing, with growth of 3.8% in 2019. Budget increases for academic and public research institutes should fuel growth. Risks could arise from a large downturn in global economic growth, because the demand for laboratory products is subject to macroeconomic effects. The U.S. market is expected to generate growth of 3.5%, in part due to the sustained strength

Future Business Development

Our 2019 guidance reflects the sector environment and economic trends, as well as the opportunities and risks outlined in this Annual Report. All forecasted figures are given in constant currencies; potential acquisitions are not considered.

Sartorius Stedim Biotech expects continued profitable growth in 2019. Consolidated sales revenue is projected to increase by about 7% to 11%. This forecast reflects the changes to the sales agreement with the Lonza group for cell culture media. Without these changes, sales growth would probably be approximately 3 percentage points higher.

Regarding profitability, management forecasts that the company's underlying EBITDA margin will increase by slightly more than one percentage point over the prior-year figure of 28.2%. Of this figure, approximately half a percentage point is expected to be an operational increase, whereas the remainder will result from changes to the IFRS accounting rules.¹⁾

The ratio of capital expenditures to sales revenue is projected to be around 11%, down from the year-earlier figure of 14.6%. Key investment projects include completing expansion of the Puerto Rican site for single-use products and increasing production capacities in Germany.

Concerning financial position, management projects that the ratio of net debt to underlying EBITDA will be slightly below the previous year's level of 0.4 reported.

In spite of countermeasures already taken, a disorderly exit of the United Kingdom from the EU may have a certain impact on our supply chain, yet a reliable forecast of possible effects cannot be given at the present time.

¹⁾ IFRS 16 required to be applied as of 2019 regulates accounting of leasing contracts. Ultimately, this will result in the disclosure of longer-term lease payments as depreciation and, accordingly, to a somewhat higher EBITDA, among other things. This will not entail any material changes concerning the Group's relevant net profit or earnings per share.

Financial Statements of the Parent Company Sartorius Stedim Biotech S.A. as of December 31, 2018

Financial Statements of the Parent Company

Sartorius Stedim Biotech S.A. is the parent company of the Group. The company is a mixt holding Company. The company from now on is managing investments of the Group and reals estates for the French Companies.

In 2018, sales revenue generated at Sartorius Stedim Biotech S.A. was €K 1,999 relative to €K 2,198 in 2017. The operating profit is €K -2,371 versus €K -3,197K in 2017. The net financing income totaled €K 48,576 versus €K 47,108 in 2017.

The net profit for 2018 is €K 49,521 compared to €K 49,463 in 2017.

Appropriation of the Net Profit

The ASM will suggest to appropriate the net profit of €49,521,306 for the reporting year of 2018. as follows:

- The following amount is to be added to this balance:
Year-earlier profit carried forward: €34,345,883
- This would yield a distributable profit of €83,867,189
- Total amount of dividends to be disbursed to shareholders: €52,540,761
- Balance resulting from disbursement: €31,326,428

The remaining amount of €31,326,428 is to be carried out to the next year.

Dividends of the last three financial years (information updated as of 1st January 2018)

The table below makes the list of the amount of the dividend per share distributed, since 2015, as well as the tax provisions applicable.

| Exercise | Dividend ¹⁾ | Amount eligible for the 40% abatement | Amount not eligible for the 40% abatement | Dividend per shares ¹⁾ |
|----------|------------------------|---------------------------------------|---|-----------------------------------|
| 2017 | 42,402,887 | 42,402,887 | 0 | 0.46 € |
| 2016 | 38,713,209 | 38,713,209 | 0 | 0.42 € |
| 2015 | 30,734,476 | 30,734,476 | 0 | 2.00 € |

¹⁾ Prior deduction of social contribution on the dividend paid to physical person.

Proposition of dividend for the 2018 financial year

The Board of Directors has decided to propose to the 26 March 2019 Annual Shareholders' Meeting a net dividend of 0,57 euros, per share for the 2018 financial year in comparison with €0.46 for 2017.

The dividends are distributed to the shareholders in ratio with the proportion of the capital held by each shareholder.

The dividend will be paid on 2 April 2019.

Dividend distribution policy

The company follows a policy of dividend distribution linked on one part to the Group's profit over the financial year concerned and on another one to the Group's predictable evolution and profitability.

The 3 April 2018 Shareholders' Meeting voted a net dividend of 0.46 euro per share. The payment of the dividend was paid on 11 April 2018.

Dividends and interim dividends paid and unclaimed are prescribed after five years in favor of the State, from their date of payment (article 2277 of the Civil Code).

Elements likely to have an impact in the event of a public offer

Pursuant to article L. 225-100-3 of the French Commercial Code, an element is likely to have an impact in the event of a public offer: the first shareholder of Sartorius Stedim Biotech S.A. holds a significant percentage of its capital and voting rights.

Sartorius Stedim Biotech S.A. Share Capital

Share Capital as of December 31, 2018

As of 31 December 2018, the share capital amounts to eighteen million four hundred and thirty-six thousand thirty-eight euros (€18,436,038). It is divided into twenty two million one hundred and eighty thousand one hundred and ninety (92,180,190) shares worth twenty cents euros (€0,20) each, all fully subscribed and paid up (Heading I, Article 6 of the bylaws), all of which are entitled to the dividend for the financial year 2018, with the exception of shares held by the Company.

| Date | Nature of the transaction | Share par value | Share capital increase | Share premium | Number of new shares | Number of shares after the transaction | Share capital after the transaction |
|------------------------------|---|-----------------|------------------------|---------------|----------------------|--|-------------------------------------|
| 1 st half of 2012 | Exercise of share subscription options | 0.61 | 5,098.0 | 173,446.0 | 8,358.0 | 17,034,306 | 10,390,926.6 |
| 2 nd half of 2012 | Exercise of share subscription options | 0.61 | 4,270.0 | 202,300.0 | 7,000.0 | 17,041,306 | 10,395,196.6 |
| Year 2013 | Exercise of share subscription options | 0.61 | 610.0 | 8,620.0 | 1,000.0 | 17,042,306 | 10,395,806.6 |
| Year 2014 | Exercise of share subscription options | 0.61 | 9,541.6 | 134,834.0 | 15,642.0 | 17,057,948 | 10,405,348.2 |
| Year 2014 | Reduction of Capital: Cancellation of Treasury Shares | 0.61 | -1,036,213.1 | | -1,698,710.0 | 15,359,238 | 9,369,135.1 |
| Year 2014 | Increase of Capital: nominal value change | 1.00 | 5,990,102.8 | | | 15,359,238 | 15,359,238.0 |
| Year 2015 | Exercise of share subscription options | 1.00 | 8,000.0 | 174,880.0 | 8,000.0 | 15,367,238 | 15,367,238.0 |
| Year 2016 | Reduction of Capital: Cancellation of Treasury Shares | 1.00 | -1,642,095.0 | | -1,642,095.0 | 13,725,143 | 13,725,143.0 |
| Year 2016 | Increase of Capital: new actions created | 1.00 | 1,638,222.0 | | 1,638,222.0 | 15,363,365 | 15,363,365.0 |
| Year 2016 | Increase of Capital: nominal value change | 0.20 | 3,072,673.0 | | 3,072,673.0 | 92,180,190 | 18,436,038.0 |
| Year 2017 | | | | | | 92,180,190 | 18,436,038.0 |
| Year 2018 | | | | | | 92,180,190 | 18,436,038.0 |

Sartorius Stedim Biotech S.A. Shareholdings as of December 31, 2018

Situation of Sartorius Stedim Biotech S.A. Shareholdings

| Shareholders | Shares | Voting rights |
|---------------------------------|--------------|---------------|
| More than 50% | Sartorius AG | Sartorius AG |
| More than 10% but less than 50% | None | None |
| More than 5% but less than 10% | None | None |

Over the past three years, the ownership of Sartorius Stedim Biotech share capital has been distributed as follows:

| Shareholders | December 31, 2016 | | | December 31, 2017 | | | December 31, 2018 | | |
|----------------------------------|-------------------|--------------------|--------------------|-------------------|--------------------|--------------------|-------------------|--------------------|--------------------|
| | Number of shares | % of share capital | % of voting rights | Number of shares | % of share capital | % of voting rights | Number of shares | % of share capital | % of voting rights |
| Sartorius AG | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% |
| Single voting rights | | | | | | | | | |
| Double voting rights | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% |
| Single voting rights | | | | | | | | | |
| Double voting rights | 0 | 0.0% | 0.0% | | | | | | |
| Total Sartorius Group | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% | 68,450,400 | 74.3% | 84.5% |
| Treasury shares | | | | | | | | | |
| Personnel and other shareholders | | | | | | | | | |
| General public | 23,729,790 | 25.7% | 15.5% | 23,729,790 | 25.7% | 15.5% | 23,729,790 | 25.7% | 15.5% |
| Single voting rights | 22,439,112 | 24.3% | 13.9% | 22,439,112 | 24.3% | 13.9% | 22,439,112 | 24.3% | 13.9% |
| Double voting rights | 1,290,678 | 1.4% | 1.6% | 1,290,678 | 1.4% | 1.6% | 1,290,678 | 1.4% | 1.6% |
| Total shares | 92,180,190 | 100.0% | 100.0% | 92,180,190 | 100.0% | 100.0% | 92,180,190 | 100.0% | 100.0% |

Legal Disclosure of Thresholds Crossed

No legal disclosure of thresholds crossed has been registered during the fiscal year under study.

| | Shares | % Issued Capital | Voting rights | % Voting rights |
|---------------------------|-------------------|------------------|-------------------|-----------------|
| Sartorius AG | 68,450,400 | 74.30 | 68,450,400 | 84.56 |
| Total Sartorius AG | 68,450,400 | 74.30 | 68,450,400 | 84.56 |

Control of the Company as of December 31, 2018

Sartorius AG holds, directly or indirectly, 74.3% of the share capital and 84.5% of the outstanding voting rights. Treasury shares are without voting rights.

Staff Shareholdings

None

Treasury Shares Held by Sartorius Stedim Biotech S.A.

None

Unpaid Capital

None

Authorized but Unissued Capital

None

Securities Not Representative of the Share Capital

None

Authority granted by the Annual Shareholders' Meeting to the Board of Directors still valid.

DELEGATION GRANTED FOR INCREASE IN CAPITAL BY THE SHAREHOLDER'S MEETING TO THE BOARD OF DIRECTORS

Delegation of competence

| Object - Duration | Limit | Use in 2018 |
|---|---|-------------|
| Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right to the allotment of the debt instruments, with preferential subscription rights of the shareholders. | The limit is €2,000,000 corresponding to the maximum nominal amount of the increase of the share capital and to the maximal nominal amount of the debt instruments and €500,000,000 on the maximum overall limit of the maximum nominal amount of the debt instruments. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through public offerings. | The limit is deducted on the overall limit of €2,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through private placements as set forth in article L411-2 II of the French Monetary and Financial Code. | The limit is deducted on the overall limit of €2,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to increase the number of shares and/or securities giving access to the share capital of the Company to be issued in the event of a share capital increase with or without preferential subscription rights of the shareholders. | The limit amount 15% of initial issue of shares. It is deducted on the overall limit of €2,000,000 (increase of the share capital) | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the Company, as consideration for securities tendered through public exchange offers initiated by the Company, without preferential subscription right of the shareholders. | The limit is deducted on the overall limit of €500,000,000 (increase of the share capital) and on the overall limit of €2,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to increase the share capital through the capitalization of reserves, earnings or premiums or any other sum upon which capitalization would be permitted. | The limit is €2,000,000 (corresponding to the maximum nominal amount of the increase of the share capital); It is a independent limit. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders and reserved for members of saving plans. | The limit is €2,000,000 corresponding to the maximum nominal amount of the increase of the share capital; it is an independent limit. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |

Other Securities Giving Access to the Share Capital

None

Stock Options

None

Share Capital Dilution

None

Share Subscription Options Granted to Each Senior Executive of the Company and Options Exercised by Them in Fiscal 2018

None

Share Subscription Options Granted to the Ten Top Non-senior Executive Beneficiaries and Options Exercised by Them in the 2018 Fiscal Year

None

Options Exercised During the Fiscal Year

All options have been exercised in 2015. The stock option plans are now expired.

| in € | 2017 | 2016 | 2015 | 2014 | 2013 |
|--|-------------|-------------|-------------|-------------|-------------|
| Dividend per share for the fiscal year | 0.46 | 0.42 | 2.00 | 1.30 | 1.20 |
| Number of shares | 92,180,190 | 92,180,190 | 15,367,238 | 15,359,238 | 15,343,596 |
| Dividend corrected per share¹⁾ | 0.46 | 0.42 | 2.00 | 1.30 | 1.20 |

¹⁾ Compared to the number of shares as of December 31, 2016

Share Subscription Plan

The stock option plans are detailed in the tables below. The authority delegated to the Board of Directors for setting up a new plan has recently expired. The Board of Directors no longer has any such delegated authority to set up any new plan.

Share Subscription Warrants

Sartorius Stedim Biotech S.A. has not issued any share subscription warrants.

Pledging of Shares

No Sartorius Stedim Biotech S.A. shares were pledged.

Pledging of Assets

None

Senior Executives

Information on Sartorius Stedim Biotech S.A. senior executives and a list of the positions they hold or have held over the past five years are included in the Corporate Governance report.

Directors' Meeting Attendance Fees

Directors' meeting attendance fees are calculated on an annual basis. The method of calculating these fees remains the same. It is as follows.

The directors receive directors' meeting attendance fees whose amount and allocation are established by the Board of Directors in consideration of the limits set by the ASM:

- Each Director receives a fixed remuneration of €25,000 per year, to be paid after the annual financial statements have been adopted by the Annual Shareholders' Meeting and which falls due for payment after the Annual Shareholders' Meeting. The chairman of the Board receives twice this amount. Furthermore, members of the Board receive an attendance fee of €1,200 per meeting and reimbursement of its expenses in addition to the annual remuneration.
- For their membership of any committee each Director receives a lump-sum amount of €4,000 per full year of membership in addition to the attendance fee of €1,200. Insofar as they hold the chair, instead of this, they receive a lump-sum amount of €8,000 per full year that they hold the chairperson in addition to the attendance fee. The remuneration for the activities on any committee is

due together with the remuneration under the terms of previous Subsection hereof.

- Any value-added tax is reimbursed by the corporation, insofar as the members of the Board are entitled to invoice the corporation separately for the value-added tax and they exercise this right.
- All these resolutions will not be applied for the Directors that got an executive top management activity at the group level. In this context, the executive corporate officers will not receive any remuneration for their membership.

A total of €268,800 is paid in directors' meeting attendance fees for 2018.

| | | Base fixed salaries € in K | Annual incentive € in K | Long Term Incentive € in K | Other € in K | Stock options € in K | Departure Indemnity € in K | Directors' meeting attendance fees € in K |
|--------------------------------------|----------------|------------------------------------|-------------------------------|----------------------------------|-----------------|----------------------------|----------------------------------|---|
| Total 2017 | 3,056.0 | 832.0 | 363.0 | 1,846.0 | 15.0 | 0.0 | 0.0 | 0.0 |
| Total 2018 | 2,522.0 | 863.0 | 455.0 | 1,189.0 | 15.0 | 0.0 | 0.0 | 0.0 |
| Joachim Kreuzburg ¹⁾ 2017 | 3,056.0 | 832.0 | 363.0 | 1,846.0 | 15.0 | 0.0 | 0.0 | 0.0 |
| Joachim Kreuzburg ¹⁾ 2018 | 2,522.0 | 863.0 | 455.0 | 1,189.0 | 15.0 | 0.0 | 0.0 | 0.0 |

¹⁾ For more details please refer to the Chapter Corporate Governance on pages 57 to 87.

Independent Auditors

The independent auditors for Sartorius Stedim Biotech S.A. are:

- KPMG S.A., represented by John Evans. Alternate auditor: Salustro Reydel.
- Deloitte & Associés, represented by Vincent Gros.

Payment Terms for Trade Payables & Receivables

| | Article D. 441-1 st : Invoices received but not paid at the date of the end of the exercise whose term has expired | | | | | | Article D. 441-2 nd : Invoices sent but not paid at the date of the end of the exercise whose term has expired | | | | | |
|--|---|---------------------------------|---------------|---------------|-------------------|---------|---|---------------------------------|---------------|---------------|-------------------|--------|
| | 0 day (indicative) | 1 à 30 days | 31 at 60 days | 61 at 90 days | 91 days and after | Total | 0 day (indicative) | 1 à 30 days | 31 at 60 days | 61 at 90 days | 91 days and after | Total |
| (A) Repartition of late payment | | | | | | | | | | | | |
| Number of concerned invoices | 31 | | | | | 31 | 4 | | | | | 4 |
| Total Amount of concerned invoices (Including all taxes) | 795,703 | | | | | 795,703 | 78,818 | 0 | | | | 78,818 |
| Percentage of Total amount of purchases including taxes for the exercise | 15% | | | | | 15% | | | | | | |
| Percentage of sales including taxes for the exercise | | | | | | | 3% | | | | | 3% |
| (B) Invoices excluded from (A) relating to disputed and contentious Receivables non recorded | | | | | | | | | | | | |
| Number of invoices excluded | 0 | | | | | 0 | | | | | | |
| Total amount of excluded invoices including taxes | 0 | | | | | 0 | | | | | | |
| (C) Reference payment terms used (Contractual or statutory period - article L. 441-6 or article L. 441-3 of Commerce Code) | | | | | | | | | | | | |
| Payment terms used for the payment term calculation | | Contractual time limit: 30 days | | | | | | Contractual time limit: 30 days | | | | |
| | | Legal time limit: | | | | | | Legal time limit: | | | | |

Five-Year Financial Results of the Parent Company Sartorius Stedim Biotech S.A.








| € in K | 2014 | 2015 | 2016 | 2017 | 2018 |
|--|------------|------------|------------|------------|------------|
| Share capital at end of period | | | | | |
| Share capital (capital stock) | 15,359 | 15,359 | 18,436 | 18,436 | 18,436 |
| Number of shares outstanding | 15,359,238 | 15,359,238 | 92,180,190 | 92,180,190 | 92,180,190 |
| Transactions and financial performance | | | | | |
| Sales revenue (excl. VAT) | 1,465 | 1,593 | 1,843 | 2,198 | 1,999 |
| Profit before tax, employee profit sharing plan, amortization, depreciation and provision expenses (and reversals) | 25,967 | 29,343 | 59,635 | 55,840 | 54,135 |
| Income tax | 468 | -653 | 4,543 | 5,552 | 3,316 |
| Contribution to employee profit-sharing plan | 0 | 0 | 0 | 0 | 0 |
| Net profit | 24,845 | 29,312 | 54,324 | 49,463 | 49,521 |
| Dividends paid or proposal of dividend | 18,412 | 19,967 | 30,734 | 38,713 | 42,403 |
| Earnings per share | | | | | |
| EPS after tax and employee profit-sharing, but before amortization, depreciation and provision expenses | 1.66 | 1.95 | 0.60 | 0.55 | 0.55 |
| EPS after tax and employee profit-sharing, amortization, depreciation and provision expenses | 1.62 | 1.91 | 0.59 | 0.54 | 0.54 |
| Dividend per share | 1.20 | 1.30 | 0.33 | 0.42 | 0.46 |
| Personnel | | | | | |
| Workforce size | 0 | 0 | 0 | 0 | 0 |
| Personnel costs | 0 | 0 | 0 | 0 | 0 |
| Social security costs | 0 | 0 | 0 | 0 | 0 |

Corporate Governance Report

03

The Board of Directors and its Committees

The Board of Directors

| | Name | Mandate | Age | Independent ¹⁾ | First appointment | Expiration of current mandate ²⁾ | Audit Committee member | Remunerations & Nominations Committee member |
|---|--------------------|--------------------------------------|-----|---------------------------|-------------------|---|------------------------|--|
|  | Joachim KREUZBURG | Chairman and Chief Executive Officer | 53 | | 2007 | 2019 | | |
|  | Susan DEXTER | Director | 63 | • | 2015 | 2021 | | |
|  | Anne-Marie GRAFFIN | Director | 57 | • | 2015 | 2021 | | |
|  | Lothar KAPPICH | Director | 61 | | 2017 | 2019 | • | • |
|  | Liliane de LASSUS | Director | 75 | *) | 2006 | 2019 | • | • |
|  | Bernard LEMAITRE | Director | 80 | | 1978 | 2019 | • | • |
|  | Henri RIEY | Director | 57 | • | 2007 | 2019 | • | • |

¹⁾ In accordance with the recommendation N° 8 of the AFEP-MEDEF code

²⁾ Directors are appointed until the date of the Annual General Shareholders' Meeting called to approve the financial statement of the previous fiscal year ending.

^{*)} In accordance with the criteria exposed within the recommendation N°8.5.6 of the AFEP-MEDEF code, Ms. Liliane de Lassus lost her independency qualification on May 19th, 2018, date of the twelfth anniversary of her first appointment as director of the company.

The company is administered by a Board of Directors composed of seven members, three of whom are independent. The directors are appointed for a three-year period.

The organization of the works of the Board and its composition must be suited to the shareholding structure, to the size and the nature of the activity of Sartorius Stedim Biotech S.A. and the particular circumstances it can face.

Composition of the Board of Directors as of 31 December 2018:

For historical reasons due to the shareholding structure of the Company, the composition of the Board of Directors and its Committees reflected the search by our reference shareholder of a long lasting balance between the Directors representing these shareholders, the Independent Directors and the executives.

Our reference shareholder takes its own responsibility towards the other shareholders, direct and distinct from the Board of Directors' one. He takes particular care to avoid possible conflicts of interests in the transparency of the information provided to the market and to fairly take all interests into account.

The Board of Directors should consider what would be the desirable balance in its membership and that of the Committees it has established, in particular in the representation of women and men, nationalities and diversity of skills by taking measures appropriate to guarantee to the shareholders and to the market that its missions are carried out with the necessary independence and objectivity. It makes public in the Reference Document the objectives, methods and results of its politics on these subjects.

Joachim Kreuzburg

Chairman and Chief Executive Officer

Date of birth: 22 April 1965

Nationality: German

First appointment: 29 June 2007

Mandate renewed: 05 April 2016

Appointed until: date of the Annual General Shareholders' Meeting in 2019 to approve the financial statements for the fiscal year ending 31 December 2018

Number of Sartorius Stedim Biotech Shares held: 6

Other current directorships and positions within the Group:

Chairman of the Executive Board (Vorstand) of Sartorius AG;
 Chairman of the Supervisory Board of Sartorius Stedim Biotech GmbH;
 Managing Director of Sartorius Lab Holding GmbH;
 Managing Director of Sartorius Corporate Administration GmbH;
 Managing Director of Sartorius Weighing Technology GmbH;
 Managing Director of SWT Treuhand GmbH;
 Managing Director of SI Weende-Verwaltungs-GmbH;
 Managing Director of SI Grone 1-Verwaltungs-GmbH;
 Managing Director of SIV Grone 2 GmbH;
 Managing Director of Sartorius Ventures GmbH;
 Chairman of the Board of Directors of Sartorius Stedim North America Inc.;
 Member of the Board of Directors of IntelliCyt Corporation;
 Member of the Board of Directors of Essen Instruments, Inc.;
 Chairman of the Board of Directors of Sartorius Stedim Filters Inc.;
 Member of the Board of Directors of Sartorius Stedim Japan K.K.;
 Member of the Board of Directors of Sartorius Stedim Lab Ltd.;
 Member of the Board of Directors of Sartorius Stedim BioOutsource Ltd.;
 Member of the Board of Directors of Denver Instrument (Beijing) Co. Ltd.;
 Chairman of the Board of Directors of Sartorius North America Inc.;
 President and Chairman of the Executive Committee of Sartorius Stedim FMT S.A.S.

Past directorships (held during the past five years) within the Group:

Vice Chairman of the Supervisory Board of Sartorius Stedim Biotech GmbH;
President of VL Finance S.A.S.;
Member of the Board of Directors of kSep Holdings, Inc.;
Member of the Board of Directors of ViroCyt, Inc.;
Member of the Board of Directors of Sartorius Hong Kong Ltd.;
Member of the Board of Directors of Sartorius Scientific Instruments (Beijing) Co. Ltd.;
Member of the Board of Directors of Sartorius Japan K.K.;
Member of the Board of Directors of Sartorius Biohit Liquid Handling Oy.

Other current directorships and positions outside the Group:

Member of the Supervisory Board (Aufsichtsrat) of Carl Zeiss AG, Germany;
Vice Chairman of the Supervisory Board (Aufsichtsrat) of Ottobock SE & Co. KGaA, Germany;
Member of the Administrative Board (Verwaltungsrat) of Ottobock Management SE, Germany;
Member of the regional Advisory Board (Regionalbeirat) of Commerzbank AG, Germany;
Member of the Economic Advisory Board (Wirtschaftsbeirat) of Norddeutsche Landesbank, Germany.

Past directorships (held during the past five years) outside the Group:

Member of the Advisory Board (Beirat) of Hameln Group GmbH, Germany;
Chairman of the Advisory Board (Beirat) of Otto Bock Holding GmbH & Co. KG, Germany.

Educational and professional background:

Diplom-Maschinenbau-Ingenieur, Dr. rer. pol. (University degree in mechanical engineering, doctorate in economics)

1992–1995 Research associate at the Institute for Solar Energy Research in Hamelin, Germany

1995–1999 Research associate at the Faculty of Economics and Management at the University of Hanover, Germany

Since
1 May 1999 Sartorius AG, Goettingen, Germany
Most recent position before promotion to the Executive Board: Vice President, Finances and Investor Relations

Since
11 Nov. 2002 Member of the Executive Board of Sartorius AG, Goettingen, Germany

1 May 2003, to
10 Nov. 2005 Spokesman (Sprecher) of the Executive Board of Sartorius AG, Goettingen, Germany

Since
11 Nov. 2005 CEO and Executive Board Chairman of Sartorius AG, Goettingen, Germany; currently responsible for Operations, Corporate Strategy, Human Resources, Legal Affairs & Compliance, Corporate Communications

Liliane de Lassus

Non-executive member

Date of birth: 29 December 1943

Nationality: French

First appointment: 19 May 2006 ¹⁾

Mandate renewed: 05 April 2016

Appointed until: date of the Annual General Shareholders' Meeting in 2019 to approve the financial statements for the fiscal year ending 31 December 2018

¹⁾ The mandate of Mrs Liliane de Lassus is continuous until today. Mrs Liliane de Lassus has been nominated member of the Board of Directors of Stedim S.A. on 19 May 2006, company which changed its name in Sartorius Stedim Biotech S.A. on 29 June 2007 when Sartorius AG acquired it.

Number of Sartorius Stedim Biotech shares held: 6

Other current directorships and positions outside the Group:

None

Educational and professional background:

Ph.D. in organic chemistry (1972), MBA (1966)

Masters' degree in Sanskrit (1969)

| | |
|-----------|---|
| 1969–1977 | Scientific employee in charge of research at the French CNRS (National Center for Scientific Research), later at the University of California, Berkeley (California, USA) |
| 1977–1981 | PSA – Automobiles Citroën Head of department; in charge of overall manufacturing planning and programming |
| 1981–1985 | Renault Automation (Robotics) Vice President of Strategic Planning |
| 1985–1989 | CEO and Chairman of the Board of a high-tech start-up company specializing in artificial intelligence (Cognitech) |
| 1989–2005 | Consultant in human resources management for company executives, especially in a multi-cultural environment |
| 2005–2007 | CEO of Stedim Biosystems |
| 2007–2008 | Executive Vice President of Sartorius Stedim Biotech |
| 2008–2018 | Managing Director of L2 L Conseil SARL (management consulting services; management of human resources) |

Bernard Lemaître

Non-executive member

Date of birth: 16 December 1938

Nationality: French

First appointment: 27 September 1978 ²⁾

Mandate renewed: 05 April 2016

Appointed until: date of the Annual General Shareholders' Meeting in 2019 to approve the financial statements for fiscal year ending 31 December 2018

²⁾ The mandate of Mr Bernard Lemaître is continuous until today. Mr Bernard Lemaître has been nominated member of the Board of Directors of Stedim S.A. on 27 September 1978, company which changed its name in Sartorius Stedim Biotech S.A. on 29 June 2007 when Sartorius AG acquired it.

Number of Sartorius Stedim Biotech shares held: 6

Other current directorships and positions outside the Group:

President of Financière de La Seigneurie S.A.S., La Ciotat;
Member of the Board of Directors of Senova Systems Inc., USA;
Member of the Board of Directors of Sycovest Asset Management, Paris;
Member of the Supervisory Board of Azulis Capital S.A., Paris;
Member of the Supervisory Board of Solon Ventures Ltd., London;
Member of the Supervisory Board of Qualium Investments S.A.S., Paris.

Educational and professional background:

| | |
|-----------|--|
| 1979–2007 | Founder, CEO and Chairman of Stedim S.A. |
|-----------|--|

Lothar Kappich

Non-executive member

Date of birth: 15 February 1957

Nationality: German

First appointment: 14 September 2017

Appointed until: the 2019 Annual General Shareholders' Meeting approving the financial statements for the fiscal year ending 31 December 2018

Number of Sartorius Stedim Biotech shares held: 1

Other current directorships and positions within the Group:

Chairman of the Supervisory Board of Sartorius AG.

Past directorships (held during the past five years) within the Group:

Member of the Supervisory Board of Sartorius AG.

Other current directorships and positions outside the Group:

None

Past directorships outside the Group:

Managing Director of ECE Projektmanagement GmbH & Co. KG, Germany.

Educational and professional background:

Doctorate (Dr. rer. pol.) in economics (subject of the doctoral dissertation: Theory of International Business Activity)

1988 - 1990 Controller in the Central Controlling Department from Schering AG in Berlin

1990 - 2017 ECE Projektmanagement G.m.b.H. & Co. KG in Hamburg, latest position: Managing Director of ECE's HR & Corporate Services as well as Managing Director of numerous subsidiaries at the ECE group

2007 - 2017 Member of the Supervisory Board of Sartorius AG, Goettingen

Since 2017 Chairman of the Supervisory Board of Sartorius AG, Goettingen

Henri Riey

Non-executive member

Independent Director

Date of birth: 5 November 1961

Nationality: Monegasque

First appointment: 29 June 2007

Mandate renewed: 05 April 2016

Appointed until: date of the Annual General Shareholders' Meeting in 2019 to approve the financial statements for the fiscal year ending 31 December 2018

Number of Sartorius Stedim Biotech shares held: 31

Other current directorships and positions outside the Group:

President of Aidea;

President of Groupe HR S.A.S.;

Director and secretary-treasurer of The Princess Grace Foundation (Monaco)

Educational and professional background:

Diplôme Institut Supérieur de Gestion (France) (degree earned at the French Higher Institute of Business Management "Institut supérieur de gestion")

1985-1988 Fund Manager at Paribas bank

1988-1996 Fund Manager, responsible for the European Equity Fund Management Team at Barclays Bank, France

1996-1999 Head of Research of Barclays Asset Management Europe

1999-2004 Executive Vice President of Barclays Asset Management; in charge of all fund management businesses

2004 - 2013 CFO of Hendyplan S.A.

Anne-Marie Graffin

Non-executive member
Independent Director
Date of birth: 3 May 1961
Nationality: French

First appointment: 7 April 2015
Mandate renewed: 03 April 2018
Appointed until: date of the Annual General Shareholders' Meeting in 2021 to approve the financial statements for the fiscal year ending 31 December 2020

Number of Sartorius Stedim Biotech shares held: 6

Other current directorships and positions outside the Group:

Member of the Supervisory Board of Valneva SE;
Member of the Supervisory Board of Nanobiotix S.A.;
Managing Director of SMAG Consulting SARL.

Past directorships (held during the past five years) outside the Group:

Member of the Board of Directors of Themis Bioscience GmbH.

Educational and professional background:

Graduated from ESSEC (Ecole Supérieure des Sciences Economiques et Commerciales)

| | |
|-------------|---|
| 1984 - 1987 | International Distillers and Vinters, France Products Manager |
| 1988 - 1990 | URGO Laboratories Marketing Manager |
| 1991 - 1995 | RoC S.A (Johnson & Johnson) - Head of International Marketing Group |
| 1998 - 2000 | Sanofi Pasteur MSD - France Products Manager Adults Vaccines |
| 2001 - 2005 | Sanofi Pasteur - Head of range then Europe Adults Vaccines Marketing Director |
| 2006 - 2008 | Sanofi Pasteur MSD - Executive Director Business Management |
| 2009 - 2010 | Sanofi Pasteur MSD - Vice President Business Management |
| Since 2011 | Managing Director SMAG Consulting SARL - Advice Biotech and Medtech Strategy Management |

Susan Dexter

Non-executive member
Independent Director
Date of birth: 11 October 1955
Nationality: American

First appointment: 7 April 2015
Mandate renewed: 03 April 2018
Appointed until: date of the Annual General Shareholders' Meeting in 2021 to approve the financial statements for the fiscal year ending 31 December 2020

Number of Sartorius Stedim Biotech shares held: 6

Other current directorships and positions outside the Group:
None

Past directorships (held during the past five years) outside the Group:
Kalon Biotherapeutics, College Station, Texas, USA - CMO;
BioSense Technologies, Woburn, Massachusetts, USA - Clinical diagnostic technology based on cellular impedance.

Educational and professional background:

Degrees and Certifications: BS in Immunology and Marketing (double major, honors), American University, Washington, D.C., USA
Harvard University Negotiation Course for Lawyers, Harvard University, Cambridge, Massachusetts, USA
Finance for non-financial Managers, Harvard University through Dow Chemical Company internal training program

| | |
|-------------|--|
| 1975 - 1980 | University of Massachusetts Medical School, Research, mammalian cell culture, animal toxicology studies, basic research |
| 1980 - 1986 | Collaborative Research, Biotechnology Sales in emerging markets for bioprocessing supplements and raw materials for biomanufacturing |
| 1986 - 1998 | Celltech Biologics, Lonza Biologics, Business Development-bioprocessing and manufacturing of biotechnology based biotherapeutics |
| 1998 - 2004 | Collaborative BioAlliance, Dow Chemical Company (Dow Biotechnology Contract Manufacturing Services) - Vice President, Business Development for microbial fermentation services, technologies |

| | |
|-------------|---|
| | and implementation of single use bioprocessing technologies |
| 2004 – 2008 | Xcellerex, Inc. (now GE Healthcare), Chief Business Officer; CMO services using fully integrated single-use bioprocessing technology, sales of single use bioprocessing technologies |
| Since 2008 | Latham Biopharm Group, Managing Director; Due Diligence, Acting VP Business Development for multiple CMO's offering contract manufacturing services to the biotechnology life sciences industry, strategic consulting, single-use disposable technology implementation, project management and high-level business development and marketing, Advisor and speak for BioProcess International, Outsourced Pharma |

Registered Addresses

With regards to their social mandates, the members of the Board of Directors and of the General Management are domiciled at the Company's headquarters.

Independent Directors

The Company being controlled by a majority shareholder, the portion of independent administrator board members should be at least a third of the Board. As of 31 December 2018, the Board of Directors of Sartorius Stedim Biotech S.A. is composed of 43% of independent members under the independence criteria defined by the APEF-MEDEF code.

Pursuant to the principles of good corporate governance, the independent members may not be principal shareholders, employees, former Group employees, suppliers or bankers of the Group or major customers, nor may they have any other link likely to impair their judgment.

In accordance with the internal rules of the Board of Directors and in application of the APEF-MEDEF code, the independence of directors is assessed each year with respect to the following criteria.

An independent director:

- May not be an employee or senior executive employee or director of his or her parent company or of one of its consolidated companies and may not have been so during the five previous years (criterion 1);
- May not be a senior executive of a company in which the company directly or indirectly holds a director's position or in which an employee as such or a senior executive of the company (either currently or having been so for less than five years) holds a director's position (criterion 2);
- May not be a significant client, supplier, business banker or investment banker of the company or of its group, for which the company or its group represents a significant part of its business (criterion 3);
- May not have any close family ties with one of the senior executives (criterion 4);
- May not have been a statutory auditor of the company for the five past years (criterion 5);
- May not have been a director of the company for more than twelve years (criterion 6).

In addition to the abovementioned criteria, the Board of Directors analyses other factors, such as the ability to understand the issues and risks, prior to making a decision on whether a director qualifies as independent.

– As part of the Assessment of the Board of Directors, the Board of Directors goes through all the criteria listed above and currently it states that it has three independent directors: Mr. Susan Dexter, Mrs Anne-Marie Graffin, and Mr Henri Riey.

| | Not an employee or executive officer | No cross-directorships | No significant business relationships | No family links | Not a statutory auditor | First appointment | Not a director for over 12 years | Classification adopted |
|--------------------|--------------------------------------|------------------------|---------------------------------------|-----------------|-------------------------|-------------------|----------------------------------|------------------------|
| Joachim Kreuzburg | No | No | Yes | Yes | Yes | 2007 | Yes | Not independent |
| Liliane de Lassus | Yes | Yes | Yes | Yes | Yes | 2006 | No ²⁾ | Independent |
| Bernard Lemaître | Yes | Yes | Yes | Yes | Yes | 1978 | No | Not independent |
| Henri Riey | Yes | Yes | Yes | Yes | Yes | 2007 | Yes | Independent |
| Susan Dexter | Yes | Yes | Yes ¹⁾ | Yes | Yes | 2015 | Yes | Independent |
| Anne-Marie Graffin | Yes | Yes | Yes | Yes | Yes | 2015 | Yes | Independent |
| Lothar Kappich | Yes | No | Yes | Yes | Yes | 2017 | Yes | Not independent |

¹⁾ Particular attention was paid to the situation of Mrs. Susan Dexter, who holds a position at Latham Biopharm Group Inc., a company with whom Sartorius Stedim Biotech GmbH signed a consultancy agreement, which expired on December 31, 2017 and has not been renewed. The Board noted that the relation between these two companies were not significant in terms of their volume. In addition, the possibility for Sartorius to use a panel of providers having similar expertise than Latham Biopharm Group, in a competitive context, rules, out all relationship of dependence.

²⁾ Mrs. Liliane de Lassus first mandate as director of the company occurred on May 19, 2006. As a consequence, the criteria N°6 of the recommendation N°8 of the AFEP MEDEF code (not being a director of the company for more than twelve years) is not fulfilled anymore since May 19, 2018. Therefore, Mrs. Liliane de Lassus correlatively lost her quality of independent director as of that date.

Balanced representation of women and men

Each year, the Board of Directors examines the desired balance in its composition and that of its committees, seeking in particular a balanced representation of men and women, and a wide diversity of skills and nationalities, reflecting as best it can both the highly technical and global nature of the company's business.

Specifically, as regards the threshold of 40% women to be reached under the provisions of Article L. 225 - 18 - 1 of the French Commercial Code, the Board of Directors has put significant effort into searching for skilled, independent and dedicated female directors with a proven level of expertise in biotechnologies or related industries. As of 31 December 2018, the Board of Directors of Sartorius Stedim Biotech S.A. is composed of 43% of women.

Assessment of the Board of Directors

The internal rules of the Board of Directors require that once a year the Board devotes an item on its agenda to discuss its functioning and ensures that a formal assessment is carried out. For this purpose, in December 2018, members of the Board completed a questionnaire on the following topics:

- the Board's composition;
- the mode and structure of governance;
- the effectiveness of the Board of Directors;
- the Board's working methods;
- the areas of competence of the Board's members;
- areas for improvement.

Consistent with last year's efficiency review, the results are satisfactory in terms of flow of information, active participation of each Board members, quality of the Committee's work. Those answers are reflecting the high quality teamwork of Board members and their convergence of views.

Board of Directors' internal rules

The Board of Directors has adopted a set of internal rules that defines and includes rules of operation for this body relating to its powers, members' attendance, operations requiring approval and prior validation with a certain number of triggering thresholds. The directors' charter is included in the Annexe and defines the rights and obligations of directors, in particular regarding the code of ethics and prevention of conflicts of interest.

Staggering of the mandate terms

According to the APEF MEDEF governance code for listed companies, the staggering of terms should be organized in order to avoid renewing a group of mandates and to promote harmonious renewing of the directors' mandates. The afferent rule relating to staggering terms is respected, as such, on December 31, 2018 two memberships have been renewed. Five membership renewals will be submitted to the Annual Combined Shareholders' Meeting on March 26, 2019, as ordinary resolutions.

Plurality of mandates

In accordance with the APEF MEDEF governance code for listed companies, an executive Director can't exercise more than two other mandates of Director in listed companies outside its group, including foreign companies. It should in addition collect the notice from the Board before accepting a new Director mandate in a listed company.

Moreover, an administrator can't exercise more than four other mandates in listed companies outside its group, including foreign companies. This recommendation is applied during the nomination or the renewal of the administrator's mandate.

Procedures established and followed by the Committees are set up within their respective internal regulations.

Committees' members are appointed by the Board of Directors. A special attention is paid by the Board of Directors to the nomination of renewal of the mandate of the Chairman of the Audit Committee, upon recommendation of the Remunerations and Nominations Committee.

Other Information

The Board of Directors met six times during fiscal 2018.

In accordance with the bylaws of Sartorius Stedim Biotech S.A. Company, each Director owns personally at least one share of the company.

All Directors fulfill the below mentioned thresholds with regards to numbers of mandates in listed companies:

- For the executive Directors: maximum of two mandates in companies not belonging to the group,
- For non-executive Directors: maximum of four mandated in companies not belonging to the group.

To the company's knowledge, within the last five years, no member of the Board of Directors:

- has been convicted of fraud during the last five years or has been subject to any official public investigation or sanction by statutory regulatory authorities;
- has been associated in his | her capacity of manager in any bankruptcy, receivership or liquidation for the past five years ;
- has been disqualified by a court from acting in the capacity of a member of an administrative, management or supervisory body of an issuer or from acting in the capacity of a management executive or conducting the business of any issuer for the past five years.

To the company's knowledge, no family relationships exist among the members of the company's Board of Directors.

Furthermore, to the company's knowledge, there is no conflict of interest between any duty of the members of the Board of Directors and their private interests and | or other duties. A Director must inform the Board as soon as he | she is aware of any conflict of interests, or even the possibility of a potential conflict, and must refrain from any participation in discussions on the relevant subject matter and from voting on any associated resolutions.

To the company's knowledge, no settlement or agreement has been reached with shareholders, clients, suppliers or others to appoint a member of the Board of Directors.

To the company's knowledge, there is no service contract linking a Board member to the Sartorius Stedim Biotech Group and granting him or her benefits.

Measures taken to ensure that control is not done in an abusive way are the following:

- Three members of the Board of Directors out of seven are Independent Directors.
- One independent member of the Board out of three is a member and Chairman of the Audit Committee.
- One independent member of the Board out of three is a member of the Remunerations and Nominations Committee.

Conditions for Preparation and Organization of the Work of the Board of Directors

Internal Rules and Regulations

The procedures governing the organization and functioning of the Board of Directors are defined by the Internal Rules and Regulations of the Board which is published on the website of Sartorius Stedim Biotech S.A. as of the publication of this particular report.

The Internal Rules and Regulations currently applicable has been adopted on 5 April 2016 and will be updated by the Board of Directors in 2019 in order to bring it into line with the new provisions of the AFEP MEDEF governance code for listed companies.

The Board of Directors deals with all matters concerning the proper operation of the company and takes decisions on subjects that concern it.

Its Missions

The main missions of the Board of Directors are as follows:

- The Board of Directors shall define the company's strategic goals and assess them from an overall perspective at least once a year, as proposed by the CEO, and ensure that these goals are implemented. It shall also appoint the corporate officers responsible for managing the company in pursuit of this strategy and review all delegations of authority;

- The Board of Directors shall review the management of the Group and monitor the quality of information provided to shareholders and to the market through the financial statements or when material events occur, especially about the company's shareholdings;
- The Board of Directors is responsible for approving all strategic investment projects and any transaction, in particular acquisitions or disposals, likely to materially affect the company's results, the structure of its balance sheet or risk profile;
- The Board of Directors will beforehand decide for each significant transaction outside the scope of the announced strategy;
- The Board of Directors shall deliberate prior to making any changes to the management structure of the company, and shall be informed of the principal organizational changes;
- The Board of Directors shall examine the corporate and consolidated accounts and approve the management report and the sections of the annual report dealing with corporate governance and those setting out the company's policies with respect to remuneration and stock options;
- Although it is not a modification with a social purpose, the Board of Directors must seize the Shareholders' Meeting if the transaction concerns a preponderant share of the assets or the activities of the group;
- The Board of Directors shall convene annual shareholders' meetings and propose changes to the articles of association.

The missions mentioned above summarize the internal bylaws of the Board of Directors.

Activity Report of the Board of Directors for Fiscal 2018

The Board of Directors met six times during the fiscal year. The average attendance was 97.62%.

The Board reviewed and approved the corporate and consolidated accounts for 2017.

The Board of Directors considered and debated on the following at its meetings:

- Strategic direction and major Group projects.
- The annual, half-year and quarterly financial statements.
- Budgets presented by executive management.
- Information on the financial structure and cash flow items.
- Significant off-balance sheet commitments.
- Risk indicators for the Group.
- New internal organization.
- Stock market performance.
- Self-assessment of the Board members.
- Elements of remuneration due or attributed.
- Renewal of the mandates of members of the Board of Directors
- Follow up on recent acquisitions.

Information to be Provided to Directors

Before each Board Meeting, Directors receive a report on the agenda items that require prior consideration, in due time and following notification.

Preliminary figures of the annual and interim statements are generally sent to all Directors at least one week before the meeting of the Audit Committee, which is always held on the day of or on the day before the Board meeting.

In addition to Board meetings, the Chairman regularly informs the Directors of any event or development that may have a material impact on Group operations or on any information previously communicated to the Board.

The members of the Board of Directors receive a copy of each press release published by the Company. The Directors may, at any time, request further information from the Chairman of the Board, who shall assess the relevance of the request.

The Audit Committee and the Remuneration and Nominations Committee are responsible for studying and making preparations for the Board's main deliberations in order to improve the Board's efficiency.

Under no circumstances do these Committees relieve the Board which has the only legal power of decision nor are allowed to cause division within its college which is and stays responsible of the accomplishment of its missions. The Committees don't replace but are an emanation of the Board of Directors facilitating its work.

The Committees of the Board may consult, in the performance of their functions, any of the main company's executive members after having informed the Chairman of the Board of Directors and subject to reporting back to the Board.

The Committees of the Board may request external technical studies relating to matters within their competence, at the expense of the Company, after having informed the Chairman of the Board of Directors or the Board of Directors itself and subject to reporting back to the Board.

In the event the Committees solicit the services of external counsels (e.g. the Remuneration Committee for the purpose of gathering information related to systems and levels of remunerations applicable within the main markets), the Committees shall ensure the objectiveness of the sought advice.

Each Board of Directors meeting is preceeded with at least one meeting of one of the two committees, depending on the addressed topics. The Committees remain accountable to the Board of Directors and address to them their position, advice and recommendations.

Procedures established and followed by the Committees are set up within their respective internal regulations.

Committees' members are appointed by the Board of Directors. A special attention is paid by the Board of Directors to the nomination of renewal of the mandate of the Chairman of the Audit Committee, upon recommendation of the Remunerations and Nominations Committee.

The Audit Committee

The Audit Committee assists the Board of Directors in areas relating to accounting policy, reporting, internal and external control, financial communication and management of the risks to which the company is exposed.

Audit Committee duties

Regarding accounting policy and internal control, the Audit Committee has the following duties:

- To proceed as soon as possible, and in any event prior to examination of the annual parent company financial statements and, where appropriate, the consolidated financial statements by the Board of Directors, with the review of all the financial, interim and annual parent company and, where appropriate, consolidated financial statements, including their notes and, where appropriate, the management report presented by the Board of Directors to the General Meeting of Shareholders called to approve the financial statements for the year ended and to present its observations to the Board of Directors. During the examination of the financial statements, the Committee pays particular attention to significant transactions that could have given rise to a conflict of interests;
- To ensure the pertinence of the selected methods and accounting procedures chosen by the company and to check their proper application;
- To check the accounting treatment of any significant transaction made by the company;
- To ensure that the internal procedures for data collection and control are sufficient to ensure the quality and reliability of the annual parent company financial statements and, where appropriate, the company's consolidated financial statements;
- To examine the scope of the consolidated companies and, where appropriate, the reasons for which any companies are not included.

Regarding external control, the Audit Committee has the following duties:

- To submit to the Board of Directors recommendations concerning the Statutory Auditors in view of their appointment or renewal by the General Meeting of Shareholders, to analyse and issue an opinion on the definition, extent and timetable of their assignment and their fees. For this purpose, the Committee steers the selection procedure for the Statutory Auditors and submits to the Board of Directors a recommendation on the Statutory Auditors proposed for appointment by the General Meeting of Shareholders. The Committee proposes to the Board the selection procedure and, in particular, whether a call for tender should be issued. It supervises the call for tender and approves the specifications and the selection of the companies consulted, taking care to select the "best bid" and not the "lowest bid";
- To ensure the independence of the Statutory Auditors.

Regarding risk analysis and prevention, the Audit Committee has the following duties:

- To analyse all disputes, including fiscal, that may have a significant impact on the parent company financial statements and, where appropriate, the company's consolidated financial statements, or its financial position;
- To examine the company's exposure to significant financial risks. The Committee examines the risks and significant off-balance sheet commitments and assesses the importance of malfunctions or weaknesses that it is made aware of and informs the Board, as appropriate;
- To review the conclusions of internal audit reports;
- To verify the satisfactory application of internal controls and information reporting procedures.

Regarding financial communication, the Audit Committee's duties include reviewing the company's financial communication projects relating to the annual and interim parent company financial statements, as well as quarterly sales.

Given the extent of its remit, the Audit Committee consults with the Statutory Auditors, but also with the Finance, Accounts and Treasury Directors. These meetings may be held, at the Committee's request, without the Company's executive bodies being present.

Composition of the Audit Committee

The Audit Committee comprises at least three members chosen by the Board of Directors for their accounting and finance expertise, of whom one must be an independent member.

The independence criteria retained by the Audit Committee's internal rules are based on those proposed by the recommendations of the AFEP MEDEF code and the Ethics code and adapted to suit the company's size, organization and means.

Audit Committee's internal rules

The Audit Committee has adopted a set of internal rules designed to provide a framework for its duties and operation and, in particular, to ensure the implementation and application of independence criteria for its members. It also includes the conditions for remuneration of the latter.

As of 31 December 2018, the Audit Committee has four members:

- Mr. Henri Riey, Chairman of the Committee since December 5, 2007
- Mr. Lothar Kappich
- Mrs. Liliane de Lassus
- Mr. Bernard Lemaître

The Chairman of the Audit Committee is independent. The Board of Directors will meet after the Annual Shareholders Meeting called to approve the financial statement of fiscal year ending on December 31, 2018 in order to modify the membership of the Audit Committee and ensure that the composition of the Committee complies with the recommendation N°15.1 of the AFEP MEDEF code for the fiscal year 2019.

The Chairman of the Board of Directors, who is also the CEO of the Group, is a permanent guest of the Audit Committee, but has no voting rights.

The Audit Committee met five times during fiscal 2018.

Remunerations and Nominations Committee

Remunerations and Nominations Committee duties

The Remunerations and Nominations Committee's purpose is to assist the company's Board of Directors in setting the remuneration policy for corporate officers and, in particular, relating to incentive mechanisms (allocation of stock options and bonus shares) that the company may implement.

During the year, the Remunerations and Nominations Committee may consult all the company's executive members, after it has informed the Chairman of the Board of Directors, and must report on this to the Board.

The Remunerations and Nominations Committee's duties also include assisting the Board of Directors with the appointment of new Board members.

Composition of the Remunerations and Nominations Committee and functioning

As of 31 December 2018, the Remunerations and Nominations Committee has four members:

Mr. Lothar Kappich, appointed member of the Remunerations of Nominations Committee during the Board of Directors meeting that took place on October 10, 2017, was appointed Chairman of the Committee by its members during the meeting held on 15 February 2018.

- Mrs. Liliane de Lassus
- Mr. Henri Riey
- Mr. Bernard Lemaître

One of the four members of the Remunerations and Nominations Committee is independent. The Board of Directors will meet after the Annual Shareholders Meeting called to approve the financial statement of fiscal year ending on December 31, 2018 in order to modify the membership of the Remunerations and Nominations Committee and ensure that the composition of the Committee complies with the recommendations N°16.1 and N°17.1 of the AFEP MEDEF code for the fiscal year 2019.

The Remunerations and Nominations Committee met once in fiscal 2018.

Report on Corporate Governance

1. Regulated Agreement

Continued agreement

The Company has decided to continue the services agreement between the Company and Sartorius AG, made effective retroactively since January 1st, 2015 and adopted by the Annual Shareholders meeting of April 4th, 2017, and covering the recharge of services of Mr. Joachim Kreuzburg.

Consequently, and in accordance with the provisions set out in Article L. 225-40-1 of the French Code of Commerce, the Company will proposed to its Board of Directors of February 7th, 2019 and further to its Annual Shareholders meeting of March 26th, 2019 to approve the continuation of the said agreement (as below detailed).

The said agreement contains the following modalities:

Nature: General assistance and administrative services

Purpose: formalization of the recharges between the Company and its parent company.

Amounts:

For Mr. Joachim Kreuzburg

Year 2017: 759 996 €

Year 2018: 674.216€

This regulated agreement has been rejected by the Annual Shareholders Meeting of April 3rd, 2018.

New agreement as of 31 December 2018

None.

2. Regulated commitments concerning Mr. Joachim Kreuzburg

There are certain commitments described in this section that are regarded as regulated under French Regulation.

Such commitments were subscribed by Sartorius AG in accordance with the global remuneration policy of the Group; 20% of their total amounts are re-charged to the Company.

These commitments were rejected by the Annual Shareholders meetings of 4 April 2017 and of April 3rd, 2018, but were approved by the Board of Directors on 16 February 2017. They will be submitted to the approval of the Annual Shareholders meeting called on March 26, 2019 to approve the financial statements for the fiscal year ending 31 December 2018.

These commitments subscribed by the German parent company comply with the German law.

Earlier departure severance

The service contract of Joachim Kreuzburg includes a severance pay cap of a maximum of two annual salaries to cover cases in which Sartorius AG Executive Board membership is terminated prematurely.

Non-competition clause

Joachim Kreuzburg has a post-contractual non-competition obligation, which is in accordance with German law. This obligation will last for two years after an Executive Board member has left the Group. During this time, if the non-competition clause is not waived or terminated, this Executive Board member may claim half of his most recent annual remuneration received from the company.

Pension commitments

Mr. Joachim Kreuzburg benefit from a supplementary pension scheme that is applicable under German Law. These commitments and their modalities are exhaustively described in the section Remuneration Report of this annual report

Corporate Governance Code / AFEP MEDEF

Since fiscal 2008, the Sartorius Stedim Biotech S.A. Board of Directors decided to follow the AFEP-MEDEF recommendations, as revised in June 2018, as the reference code for corporate governance (see www.medef.fr).

The AFEP-MEDEF Corporate Governance Code (the "Code") defines a set of regulations for good and responsible corporate governance. It follows the "comply or explain" principle that is implemented in most countries of the European Union. If a listed company does not comply with a recommendation of this Code, it must explain this in its corporate governance report.

In accordance with article 27.1 of the Corporate Governance Code for listed companies in effect from the presented date (the "Code"); listed companies referring to the code are required to precisely identify, in their Registration Document, the application of these recommendations. In case on non-application of one of the-se provisions, companies are required to provide a comprehensible, relevant and circumstantial explanation according to the rule "apply or explain". It is recommended by the AMF (recommendation n°2014-08 of 22 September 2014) that companies indicate in a specific table each recommendation that are not applied and the related explanations.

SPECIFIC TABLE ON RECOMMENDATIONS OF THE AFEP MEDEF CODE FOR THE GOVERNANCE OF LISTED COMPANIES

| ARTICLE | DEVIATIONS OF THE PROVISIONS OF THE CODE | EXPLANATIONS |
|---------|--|---|
| 3.2 | <p>Disclosure of the option selected</p> <p>It is essential for the shareholders and third parties to be fully informed of the choice made between separation of the offices of Chairman and Chief Executive Officer and maintenance of these positions as a single office.</p> | <p>The Board of Directors has opted for the Chairman's functions meeting of the Board Committee and as Chief Executive Officer in order to simplify the company operational management and increase its effectiveness.</p> <p>This organization turned out to be a factor of efficient governance considering the organization of the Sartorius Stedim Biotech Group. Mr Joachim Kreuzburg is Chairman of the Board and CEO of Sartorius AG mother company of the group. He is on one hand bound to the controlling shareholder and on the other hand very involved in the business affairs of the Group which he particularly knows and experienced.</p> <p>Also, the Board of Directors is proceeding to an annual evaluation of its functioning to identify the improvements that could be made. The result of the evaluation shows that this organization is well suited for the interests of the company.</p> |
| 10.3 | <p>Non-executive directors meeting</p> <p>It is recommended that the non-executive directors meet periodically without the executive or "in-house" directors. The internal rules of operation of the Board of Directors must provide for such a meeting once a year, at which time the evaluation of the Chairman's, Chief Executive Officer's and Deputy Chief Executive's respective performance shall be carried out, and the participants shall reflect on the future of the company's executive management.</p> | <p>Board meetings are organized in the presence of the executive members to maintain the same degree of information between the members of the Board and strengthen the open and transparent collective character.</p> <p>According to the Code AFEP-MEDEF planning that the non executive members have to meet annually without the presence of the executive or internal members, the internal rules of the Board mentions the possibility for the non executive members to organize this kind of meeting. The concerned Directors have duly been made aware of this possibility but did not express their wish to hold such a meeting during the past year. They will be reminded of such a possibility during the first Board of Directors Meeting that will occur for the next financial year.</p> |
| 15.1 | <p>Independent directors within the Audit Committee</p> <p>The proportion of independent directors on the audit committee (excluding the directors representing employee shareholders and directors representing employees, who are not taken into account) should be at least equal to two-thirds, and the committee should not include any executive director.</p> | <p>On December 31, 2018, 25% of the Audit Committee members are independent (i.e. one member out of four). This is the direct consequence of the loss of the status of independent of one of its member during the fiscal year ended on December 31, 2018.. The Audit Committee is chaired by the independent administrator: Mr. Henri Riey. The Board of Directors will met after the Annual Shareholders Meeting called in March 26, 2019 to approve the financial statement of fiscal year ending on December 31, 2018 in order to modify the membership of the Audit Committee and ensure that the composition of the Audit Committee complies with the recommendation of the AFEP MEDEF code for the fiscal year 2019., in addition to ensuring the efficiency of their respective work, requiring a high expertise in finance or accounting matters.</p> |
| 15.3 | <p>Examination deadline of the accounts between the Audit Committee and the Board</p> <p>The appointment or extension of the term of office of the audit committee's Chairman is proposed by the appointments/nominations committee, and should be specially reviewed by the Board.</p> | <p>For practical reasons, connected in particular to the presence within the Committee of a majority of non resident members, the meetings of the Audit committee usually take place the same days as those of the Board of Directors. Taking into consideration this obligation, and in order to give to the Audit committee the possibility of achieving completely its missions, the internal rules of the Board mentions that any documents and useful information must be communicated to the Board by the Chairman and Chief Executive Officer upfront and in a sufficient delay. The files are like this transmitted to the members of the Audit Committee with a sufficient upstream delay and at the latest three days before every meeting of the Committee or of the Board allowing them to have a sufficient delay for the examination of the statements before these meetings.</p> <p>Therefore, each member of the said committee is spending the necessary time to examine each topic and is duly enabled to require such information if needed.</p> |

| ARTICLE | DEVIATIONS OF THE PROVISIONS OF THE CODE | EXPLANATIONS |
|--------------|--|---|
| 16/17 | THE COMMITTEE IN CHARGE OF REMUNERATIONS AND NOMINATIONS | |
| 16.1/17.1 | Independent directors within the Remunerations and Nominations Committee It must mostly consist of independent directors It is recommended that the Chairman of the committee should be independent. | <p>The Board of Directors decided to create a Remunerations and Nominations Committee. To this end the committee in charge of remunerations and nominations has updated its internal regulations to state clearly its new functions.</p> <p>On December 31, 2018, 25% of the Remunerations and Nominations Committee members are independent (i.e. 1 member out of four). This is the direct consequence of the loss of the status of independent of one of its member during the fiscal year ended on December 31, 2018. The Board of Directors will meet after the Annual Shareholders Meeting called in 2019 to approve the financial statement of fiscal year ending on December 31, 2018 in order to modify the membership of the Remunerations and Nominations Committee and ensure that the composition of such Committee complies with the recommendations of the AFEP MEDEF code for the fiscal year 2019. It is further mentioned that the Chairman and Chief Executive Officer of the Board is not a member of the committee.</p> <p>For historical reasons related to the company share options, the composition of the specialized committee was reflecting the research by our shareholder in order to reflect a balance between the directors representing the shareholders and the independent directors.</p> <p>Lothar Kappich has been appointed Chairman of the Remunerations and Nominations Committee of the Sartorius Stedim Biotech Group for management and coherency reasons: Although Mr. Lothar Kappich is non-independent, he is also the Chairman of the remunerations committee of the Sartorius Group AG</p> |
| 19. | ETHICAL RULES FOR DIRECTORS | |
| | The director should be a shareholder personally and hold a fairly significant number of shares to the received attendance fees: by default if he does hold the shares upon assuming his functions, he must use the acquired attendance fees when acquired. | <p>The Board of Directors has implemented these ethic principles within its internal regulations, in particular within the Director Charter, which is attached to the internal regulations.</p> <p>Beyond the application of Article L 225 - 25 of the French Code of Commerce, the Board of Directors has left until now the freedom to each director to invest significantly or not within the company.</p> |
| 21. | TERMINATION OF EMPLOYMENT CONTRACT IN THE EVENT OF BECOMING A COMPANY OFFICER | |
| 21.1 | When an employee is appointed as a company officer, it is recommended to terminate his or her employment contract with the company or with a company affiliated to the group, whether through contractual termination or resignation | <p>This recommendation is not applied since there are no Directors under an employment contract with the company.</p> <p>According to German law, it is not necessary to change such an employment contract when a person becomes a Managing Director of the company he/she works for. It should also be considered that the Sartorius Stedim Biotech Group is controlled by a German majority shareholder, and the biggest group company is a German company; therefore, in this respect German rules and regulations are very common in the whole group and have to be observed at the respective group level.</p> |

| | | |
|--------|--|---|
| 24. | COMPENSATION OF COMPANY OFFICERS | |
| 24.3.2 | <p>Annual variable compensation of executive officers</p> <p>The rules for fixing this compensation must be consistent with the annual review of the performances of the executive officers and the corporate strategy. They depend on the director's performance and the progress made by the company</p> | <p>Mr Joachim Kreuzburg is representing the Group Sartorius AG, his compensation policy is deliberated and decided at the level of the headquarter Sartorius AG.</p> <p>The performance action elements are detailed in the document reference within the parts of the corporate governance Report and the internal control within this Registration Document. It is also reminded that this variable compensation is exclusively allocated by Sartorius AG, and, as such, performance rules are decided under the German legislation, through applicable governance codes.</p> |
| 24.5.1 | <p>Departure of company officers</p> <p>It is not acceptable that directors whose company has failed or who have personally failed may receive benefits upon departure.</p> | <p>Severance payments for Joachim Kreuzburg are defined in the course of the remuneration program at the headquarter level Sartorius AG, and are capped. Further and detailed information related to these severance payments are available within the report on the remuneration of the Executive Director inserted in this Registration Document.</p> |

Shareholders' Meeting

Convening

Annual (or Ordinary) General Shareholders' Meetings are those convened to take all decisions that do not result in a revision of the bylaws. Extraordinary General Shareholders' Meetings are those called to decide or authorize direct or indirect revisions to the bylaws. Special Meetings bring together the holders of a specific class of share to consider revisions to the rights of this class of share. Decisions made at the General Meetings are binding for all shareholders, even those who are absent, dissenting or legally incapable or incapacitated (Company bylaws, Heading 3, Article 13). General Meetings are convened by the Board of Directors or, by default, the independent auditors or a person thus empowered. General Meetings are held at the registered office or any other place stated in the notice of convocation (Company bylaws, Heading 3, excerpt of Article 14). The forms and timescale of the notice of convocation are governed by French laws.

Agenda

The notices and letters of call shall indicate the indications required by the law, particularly the agenda, the company electronic address where written questions of Shareholders may be sent and, eventually the mention of the obligation to collect the opinion or the prior approval of the mass of securities Shareholders giving access to the share capital.

The meeting may only deliberate on the matters placed on the agenda. It may, however, remove one or more directors at any time.

One or more shareholders representing the percentage of share capital required by law may, under the conditions and time limits set forth by law, require the inclusion on the agenda of draft resolutions.

In accordance to the Articles R 225-71 to R 225-74 of the Commercial Code, requests made by the Shareholders to register draft resolutions on the agenda and written questions are sent to the Headquarters by registered letter with recorded delivery beginning on the publication of the Meeting announcement and until 25 days before the General Meeting, or in a delay of 20 days beginning on the publication of the Meeting announcement, when this one is published more than 45 days before the General Meeting (date of reception of the request by the company will be taken into account).

The request of a new item on the agenda must be motivated. The request to register draft resolutions is provided with the text of draft resolutions, which may have a short explanation of reasons. These requests are subject to justification of possession or representation of required Share capital, in accordance to regulatory rules (Company bylaws, Heading 3, Article 14, excerpt of point 2).

Moreover, in accordance to the Articles L. 2323-67 paragraph 2 of the Labor Code, requests of draft resolutions made by the Work Council, to be added on the agenda, are sent in the next 10 days following the publication of the Meeting announcement. (Company bylaws, Heading 3, Article 14, excerpt of point 2).

If the meeting has been unable to make a valid decision due to a lack of the required quorum, the second meeting and, where appropriate, the second meeting adjourned are called at least ten days in advance in the same form as the first meeting (Company bylaws, Heading 3, Article 14, excerpt of point 1).

Admission to Meetings – Powers

Every shareholder has the right to attend General Meetings and to participate in the discussions, in person or by proxy, regardless of the number of shares held, on simple proof of identity and the ownership of shares. The right to participate in a General Meeting is subject to the condition that the shares must be recorded, in the name of the shareholder or the shareholder's appointed broker, either in the nominative share accounts held by the company or in the bearer share accounts held by the authorized broker, by zero hours, Paris time, on the second working day prior to the meeting. The recording or registration of the shares in the bearer share accounts held by the authorized broker must be confirmed by a share certificate provided by the broker. This share certificate must be attached to the postal voting form, the proxy form or the application for an admission pass, issued in the name of the shareholder or on behalf of the shareholder represented by the appointed broker. A certificate must also be supplied to shareholders who wish to attend the General Meeting in person but who have not received an admission pass by zero hours, Paris time, on the second working day prior to the meeting.

A Shareholder may be represented by another Shareholder, his or her spouse or by the partner with who he or she signed a Civil Partnership. Furthermore, he or she may be represented by any other moral or physical person of his choice in accordance to the Articles L. 225-106 to -106-3 of the Commercial Code; in that aim, the representative must present valid proof of proxy.

The legal representatives of shareholders who are legally incapable or incapacitated and individuals representing corporate shareholders take part in meetings, whether or not they are shareholders (Company bylaws, Heading 3, Article 14, point 3).

All Shareholders may also have a postal voting, using a registration form and sent to the company according to the law and regulations; to be acceptable this registration must be received by the company three days before the date of the Meeting.

In case of remote voting using an electronic vote, or a proxy vote given by electronic signature, this vote is made according to the conditions of the current regulations (Company bylaws, Heading 3, Article 14, point 4).

All legal documents relative to legal information for shareholders are made available to them at the registered office of the company.

Delegation granted for increase in capital by the Shareholders' meeting to the Board of Directors

Delegation of competence

| Object - Duration | Limit | Use in 2018 |
|---|---|-------------|
| Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right to the allotment of the debt instruments, with preferential subscription rights of the shareholders. | The limit is €2,000,000 corresponding to the maximum nominal amount of the increase of the share capital and to the maximal nominal amount of the debt instruments and €500,000,000 on the maximum overall limit of the maximum nominal amount of the debt instruments. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the company and/or securities giving the right of the allotment of debt instruments, without preferential subscription rights of the shareholders - through public offerings. | The limit is deducted on the overall limit of €2,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the Company and/or securities giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders - through private placements as set forth in article L411-2 II of the French Monetary and Financial Code. | The limit is deducted on the overall limit of €2,000,000 (increase of the share capital) and on the overall limit of €500,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to increase the number of shares and/or securities giving access to the share capital of the Company to be issued in the event of a share capital increase with or without preferential subscription rights of the shareholders. | The limit amount 15% of initial issue of shares. It is deducted on the overall limit of €2,000,000 (increase of the share capital) | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital of the Company, as consideration for securities tendered through public exchange offers initiated by the Company, without preferential subscription right of the shareholders. | The limit is deducted on the overall limit of €500,000,000 (increase of the share capital) and on the overall limit of €2,000,000 (debt instruments). | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to increase the share capital through the capitalization of reserves, earnings or premiums or any other sum upon which capitalization would be permitted. | The limit is €2,000,000 (corresponding to the maximum nominal amount of the increase of the share capital); It is a independent limit. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |
| Ability to issue shares and/or securities giving access to the share capital giving the right to the allotment of debt instruments, without preferential subscription rights of the shareholders and reserved for members of saving plans. | The limit is €2,000,000 corresponding to the maximum nominal amount of the increase of the share capital; it is an independent limit. | None |
| Granted for a period of 26 months as from 03 April 2018 | | |

Remuneration of the Members of the Board

Information about the Remuneration of the Executive Director

The Executive Director, Joachim Kreuzburg, is at the same time Chairman of the Executive Board of the major shareholder of Sartorius Stedim Biotech S.A. He receives his fixed and variable remuneration from the major shareholder Sartorius AG. A portion of his remuneration is charged to Sartorius Stedim Biotech S.A., reflecting his role as Executive Director of the Company. A portion of his total remuneration is charged to the SSB Group for his management services based on his proportional work for Sartorius Stedim Biotech (please refer also to section "Related Parties" of the "Financial Statements and Notes"). This allocation key is applied to all components of his remuneration.

Remuneration of the Executive Director who is chairman of the Executive Board of the major shareholder Sartorius AG (Joachim Kreuzburg)

General and Fixed Remuneration

The total amount of the remuneration of the chairman of the Executive Board of Sartorius AG, as well as for other members of this Executive Board, reflects the scope of the responsibilities of the Executive member concerned, the Executive member's personal performance, the company's economic situation and sustainable progress. In addition, this amount is benchmarked with those at peer companies and with the vertical remuneration structure within the company as well as at peer companies. Remuneration is comprised of both fixed non-performance-based components and of variable performance-based components, and is reviewed annually to ensure that it remains appropriate. The variable performance-based remuneration components consist of those to be paid annually and of multi-year components intended to have a long-term incentive. Fixed non-performance-based remuneration is paid in the year in which it is granted. For 100% target achievement, the variable annual and long-term performance-based components generally represent half of total remuneration, which excludes pension commitments under a defined benefit plan as well as fringe benefits. The targets set for the performance-based remuneration refer to financial key figures of the Sartorius Group in which the Sartorius Stedim Biotech Group is fully consolidated. Specifically, Sartorius Stedim Biotech represents approx. 80% of the business and assets of the Sartorius Group. Therefore, the development of

Sartorius Stedim Biotech has a significant influence on the financial results of the Sartorius Group and thus on the variable remuneration of Sartorius AG's Executive Board members. However, all components of the remuneration described below refer to parameters and financial key figures of the Sartorius Group in total.

Variable Remuneration

The variable portion of this remuneration contains components that are paid annually (subordinate targets measured against sales revenue | order intake, underlying EBITDA and ratio of net debt to EBITDA) and components determined by multi-year assessment (measured against (i) consolidated net profit and (ii) the phantom stock plan).

The components to be annually paid and the elements determined by multi-year assessment each make up one half of the target achievement that is possible. A cap is provided for all variable components to be paid.

Of the total that can be awarded for 100% target achievement, the subordinate targets of the components to be annually paid are weighted as follows:

- sales revenue | order intake 15%;
- EBITDA 20%;
- ratio of net debt to underlying EBITDA 15%.

The subordinate targets constituted by (i) consolidated net profit and (ii) the phantom stock plan as components determined by multi-year assessment are each weighted at 25%.

a) Annually paid variable remuneration

The portion of the variable remuneration that is to be paid annually depends on the degree to which the target is achieved, which the Supervisory Board of Sartorius AG defines by setting each individual subordinate target. Thus, target achievement is subdivided into the previously mentioned three subordinate targets, which are each separately paid.

Sales Revenue | Order Intake

If the degree of target achievement is below 90%, no remuneration is paid. If 90% is achieved, 50% of the sum awarded is paid out. Thereafter, payment increases linearly up to a target achievement of 104%, at which a maximum of 120% of the sum awarded is paid out. The degree of payment of 120% constitutes the cap for this subordinate target at the same time.

Underlying EBITDA

If the degree of target achievement is below 70%, no remuneration is paid. If 70% is achieved, 70% of the sum awarded is paid out. Thereafter, payment increases linearly up to a target achievement of 120%, at which a maximum of 120% of the sum awarded is paid out. The degree of payment of 120% constitutes the cap for this subordinate target at the same time.

Ratio of Net Debt to underlying EBITDA

No remuneration is paid if the ratio of net debt to underlying EBITDA achieved is below the lower limit defined. If this defined value is achieved, 50% of the sum awarded is paid out. Thereafter, payment increases linearly up to a target achievement of 120%, at which a maximum of 120% of the sum awarded is paid out. The degree of payment of 120% constitutes the cap for this subordinate target at the same time.

b) Variable remuneration with multi-year components

On the one hand, components determined by multi-year assessment depend on the degree to which the target is achieved, which the Supervisory Board of Sartorius AG defines by setting the subordinate target constituted by consolidated net profit. On the other hand, these multi-year components depend on the value of the monetary sum ascribed to the Executive Board member at the beginning of each year.

Consolidated Net Profit

For this subordinate target, the basis for assessment is the consolidated net profit after non-controlling interest excluding amortization (amortization of the value of intangible assets, such as customer databases or patents, which results from purchase price allocation within the scope of business combinations pursuant to IFRS 3). Target achievement for assessing annual variable remuneration is based on the average taken over a period of three fiscal years, beginning with the present fiscal year.

To smooth the amounts to be paid out, a partial payment amounting to 50% of the target achievement for a fiscal year will be effected. Any overpayments as a result of these partial payments will be offset in the following year against other remuneration components (fixed or variable). No partial payment will be made in the year prior to an Executive Board member's resignation. Full account is thus taken of any negative results, and the effects thereof continue to have an impact on the remuneration of the Executive Board member concerned even after he or she has left the company. If a defined minimum value is attained, payment of the awarded sum will increase linearly from 0% to a maximum of 120% of the subordinate target achievement value defined by the Supervisory Board. The degree of payment of 120% constitutes the cap for this subordinate target at the same time.

Phantom Stock Plan

Through the issue of shadow shares, called phantom stocks, the Executive Board members are treated as if they were owners of a certain number of shares in Sartorius AG, without, however, being entitled to receive dividends. The development of the value of these phantom stocks are linked with the development of the Sartorius share; both increases and decreases in the share price are taken into account. Later, this phantom stock is valued based on the share price at the time and its equivalent is paid out, provided that the associated conditions are met. Phantom stocks cannot be traded and does not entail any share subscription rights.

According to the Sartorius phantom stock plan, each Executive Board member is credited at the beginning of every year with phantom stock units valued at an agreed monetary sum. The value of these phantom stocks can be paid out only as an entire annual tranche. Payment can be requested, at the earliest, after a period of four years and no later than after eight years.

An Executive Board member is entitled to receive payment for phantom stock units only if the share price at the time of the payment request has appreciated at least 7.5% per year relative to the time the phantom stock was assigned or if the share price outperformed the TecDAX® as a comparative index. The phantom stock plan rules out subsequent changes to the parameters used for comparative stock valuation.

The amount to be paid is capped at a maximum of 2.5 times the share price at the time the phantom stocks were assigned, based in each case on the actual annual tranche concerned.

Assignment of this phantom stock and payment of its monetary equivalent depend on the mean value calculated from the average prices of the preference share in the closing auction of Xetra trading on the Frankfurt Stock Exchange over the last 20 days of trading of the previous year or over the last 20 days of trading prior to submission of the payment request. This serves to compensate for any short-term fluctuations in the share price.

Payment for phantom stocks is blocked for the four weeks preceding the scheduled publication date of quarterly and preliminary year-end results and for 20 days of trading on the stock exchange following the actual publication of quarterly and preliminary year-end results. These blackout periods are intended to prevent Executive Board members from profiting from potential insider knowledge.

Pension Commitments

According to the company's remuneration policy, Executive Board members of Sartorius AG receive performance-related benefit commitments under a defined benefit plan when reappointed for the first time. In addition to including a basic pension, these commitments provide for the Executive Board member to make his own contribution from his variable earnings and for the company to match this contribution by a bonus amount. An Executive Board member may choose to receive such defined benefits in the form of a monthly retirement pension for old

age or as a one-time payment to cover the member's retirement pension for old age and invalidity as well as in the form of survivor's benefits for the surviving spouse and children of the decedent.

Beyond such commitments, Joachim Kreuzburg is additionally entitled under a former company pension scheme to receive performance-based retirement benefits based on the salary of a German federal civil servant classified as grade 10 of salary class B for ministry officials according to the Federal Civil Service Remuneration Act [Bundesbesoldungsgesetz]. Such benefits are paid in the form of a retirement pension for old age and invalidity as well as in the form of survivors' benefits for the surviving spouse and children of the decedent.

After a member has turned 65, this shall be considered the regular age limit at which this member shall automatically be entitled to receive all such benefits.

Other Remuneration Components

The remuneration system provides that the Supervisory Board of Sartorius AG at its discretion may grant an Executive Board member special compensation based on that member's exceptional performance.

Severance Caps

The service contracts include a severance pay cap of a maximum of two annual salaries to cover cases in which Sartorius AG Executive Board membership is terminated prematurely. Potential amounts have to be paid by Sartorius AG.

Non-competition Clause

All Executive Board members of Sartorius AG have a post-contractual non-competition obligation, which is in accordance with German law. This obligation will last for two years after an Executive Board member has left the Group. During this time, if the non-competition clause is not waived or terminated, this Executive Board member may claim half of his most recent annual remuneration received from the company. It should be noted that this indemnity is paid by Sartorius AG. However, an allocation of this indemnity is recharged by Sartorius AG to Sartorius Stedim Biotech S.A at the date of its payment

Fringe Benefits

The members of the Executive Board of Sartorius AG are each entitled to use a company car, reclaim expenses incurred on business travel and to be covered by accident insurance and D&O insurance as fringe benefits in addition to receiving the remuneration components mentioned. The D&O insurance provides for the application of a deductible or excess in the amount required by law.

Share-based Payment

The remuneration policy for Executive Board members of Sartorius AG does not provide for the transfer of Sartorius AG shares as compensation for members. An exception to this was made in December 2014 for Joachim Kreuzburg in connection with his third appointment as a member of the Executive Board and its Chairmann and CEO.

Consequently he holds now 100,000 shares of each class of shares. The shares transferred are subject to a holding period that will end on November 10, 2019. Should Joachim Kreuzburg leave the company after November 11, 2017, and before November 11, 2019, at his own request, Joachim Kreuzburg shall be required to transfer half of the shares granted to him back to the company. The amount resulting since December 16, 2014, for the shares granted are to be spread as an employee benefits expense over the full vesting period and recognized as such in profit or loss. In fiscal 2018, an amount of €202 K was accordingly recognized in the accounts of Sartorius Stedim Biotech.

Information about the Remuneration of the Non-Executive Directors

The remuneration for non-executive board members is defined in the Board of Directors internal rules of Sartorius Stedim Biotech S.A. and comprises fixed remuneration, meeting attendance fees and reimbursement of out-of-pocket expenses. Members also serving as a member of a committee of the Board receive higher fixed remuneration.

Tables Summarizing the Remuneration and Options and Shares Granted to Each Executive Board Member

Joachim Kreuzburg
(Chairman of the Board and Chief Executive Officer)

| € in K | Year 2018 | Year 2017 |
|--|--------------|--------------|
| Remuneration due | 2,522 | 3,056 |
| Valuation of options granted during the reporting period | 0 | 0 |
| Valuation of the performance of shares granted in previous years | 0 | 0 |
| Total | 2,522 | 3,056 |

The amount cross-charged by the company Sartorius AG to the Sartorius Stedim Biotech Group concerning Joachim Kreuzburg is €1.348 K, the amount charged to Sartorius Stedim Biotech S.A. is submitted to the vote of the Annual Shareholders' Meeting in accordance with the AFEP-MEDEF code and amounted to €674 K.

Pension Commitments

| in T€ | Expected pension | Present value of obligation | | Service cost (IFRS) | |
|-----------------------|------------------|-----------------------------|--------------|---------------------|------------|
| | p. a. | 31.12.2018 | 31.12.2017 | 2018 | 2017 |
| Dr. Joachim Kreuzburg | 243 | 3,385 | 2,989 | 257 | 258 |
| | 243 | 3,385 | 2,989 | 257 | 258 |

Summary of the Remuneration for Each Executive Board Member

Joachim Kreuzburg¹⁾
(Chairman of the Board and Chief Executive Officer)

| € in K | Year 2018 | | Year 2017 | |
|--------------------------------|--------------|--------------|--------------|--------------|
| | Amounts due | Amounts paid | Amounts due | Amounts paid |
| Fixed remuneration | 863 | 863 | 832 | 832 |
| Variable remuneration | | | | |
| Annually paid | 455 | 363 | 363 | 418 |
| Long-term incentive | 1,189 | 588 | 1,846 | 602 |
| Exceptional remuneration | | | | |
| Director's attendance fees | | | | |
| Benefits in kind ²⁾ | 15 | 15 | 15 | 15 |
| Total | 2,522 | 1,829 | 3,056 | 1,867 |

¹⁾ Joachim Kreuzburg receives his salary from Sartorius AG for his duties performed for the entire Sartorius Group. His remuneration is determined annually by the Supervisory Board of Sartorius AG.

²⁾ Company Car

Table on Directors' Meeting Attendance Fees and Other Remuneration Received by Non-executive Board Members

| € in K | Year 2018 | Year 2017 |
|----------------------------|--------------|--------------|
| Liliane de Lassus | | |
| Director's attendance fees | 47.4 | 52.2 |
| Other remuneration | | |
| Bernard Lemaître | | |
| Director's attendance fees | 47.4 | 52.2 |
| Other remuneration | | |
| Arnold Picot | | |
| Director's attendance fees | 0.0 | 41.8 |
| Other remuneration | | |
| Henri Riey | | |
| Director's attendance fees | 50.2 | 55.0 |
| Other remuneration | | |
| Susan Dexter | | |
| Director's attendance fees | 36.2 | 37.4 |
| Other remuneration | | |
| Anne-Marie Graffin | | |
| Director's attendance fees | 36.2 | 36.2 |
| Other remuneration | | |
| Lothar Kappich | | |
| Director's attendance fees | 51.4 | 38.2 |
| Other remuneration | | |
| Total | 268.8 | 313.0 |

Performance Shares Available for Each Board Member

Performance shares available for each corporate officer

| Performance shares available for each corporate officer ¹⁾ | Date of the plan | Number of shares available during the reporting period | Acquisition conditions |
|---|------------------|--|------------------------|
| Joachim Kreuzburg | | Not applicable | |
| Liliane de Lassus | | Not applicable | |
| Bernard Lemaître | | Not applicable | |
| Lothar Kappich | | Not applicable | |
| Henri Riey | | Not applicable | |
| Susan Dexter | | Not applicable | |
| Anne-Marie Graffin | | Not applicable | |
| Total | | | |

¹⁾ The performance shares are bonus shares allocated to the Board members within the framework of the L225-197-1 articles and following of the commercial law, and which are subjected to additional requirements laid down by the recommendations AFEP/MEDEF of October 2008.

The performance shares are bonuses allocated pursuant to the provisions of Articles L225-197-1 and the following of French commercial law and according to the recommendations of the AFEP/MEDEF Code.

Performance Shares Granted to Board Members

There is no performance share program in place for the board members of Sartorius Stedim Biotech S.A.

The information provided in the table below refers to the phantom stock plan of Sartorius AG. This plan only relates to Joachim Kreuzburg who is Chairman of the Executive Board of Sartorius AG.

| Performance shares granted by the AGM during the reporting period to any corporate officer by the issuer or any other company of the Group | Date of the plan | Number of shares granted during the year | Valuation of the shares according to the consolidated accounts methodology | Date of acquisition | Date of availability | Performance conditions |
|--|------------------|--|--|---------------------|----------------------|------------------------|
| Joachim Kreuzburg | | 2,685 | 239 | Jan. 1, 2018 | Jan. 1, 2022 | |
| Liliane de Lassus | | | | | | |
| Bernard Lemaître | | | | | | |
| Lothar Kappich | | | | | | |
| Henri Riey | | | | | | |
| Susan Dexter | | | | | | |
| Anne-Marie Graffin | | | | | | |
| Total | | 2,685 | 239 | | | |

| | 2018 € in K | 2017 € in K |
|------------------------------|----------------|----------------|
| Total | 743 | 1,404 |
| Phantom Stocks | 239 | 163 |
| Sartorius AG shares granted | 504 | 1,241 |
| Dr. Joachim Kreuzburg | 743 | 1,404 |
| Phantom Stocks | 239 | 163 |
| Sartorius AG shares granted | 504 | 1,241 |

| | Number of phantom stock units | Subscription price in € | Fair value when granted on Jan. 1 of the particular year € in K | Fair value at year-end on Dec. 31, 2017 € in K | Fair value at year-end on Dec. 31, 2018 € in K | Paid out € in K | Change in fair value in 2017 | Exercisable |
|---|-------------------------------|-------------------------|--|---|---|--------------------|------------------------------|-------------|
| Dr. Joachim Kreuzburg | | | | | | | | |
| Tranche of phantom stock units for 2014 | 8,032 | 21.01 | 169 | 422 | 0 | 422 | 0 | |
| Tranche of phantom stock units for 2015 | 7,360 | 24.70 | 182 | 454 | 454 | 0 | 0 | no |
| Tranche of phantom stock units for 2016 | 3,484 | 57.41 | 200 | 225 | 359 | 0 | 134 | no |
| Tranche of phantom stock units for 2017 | 2,950 | 70.51 | 208 | 163 | 281 | 0 | 118 | no |
| Total tranches previous years | 21,826 | | 759 | 1,264 | 1,094 | 422 | 252 | |
| Tranche of phantom stock units for 2018 | 2,685 | 80.32 | 216 | 0 | 239 | 0 | 23 | no |
| Total | 24,511 | | 975 | 1,264 | 1,333 | 422 | 275 | |

Stock Options Granted During the Reporting Period to the Board Members by the Issuer or Any Other Company of the Group

Not applicable

Stock Options Exercised During the Reporting Period by Each Board Member

Not applicable

Stock Options Granted | Historical Information

Not applicable

**Stock Options Granted to the Top Ten
Non-corporate Officers and Exercised by Them**

Not applicable

Additional Information about the Executive Board Members

| Corporate officer | Employment contract | | Additional pension plan | | Indemnities or compensation due with regard to termination of contracts or positions | | Non-competition clause indemnities | |
|---------------------------------------|---------------------|----|-------------------------|----|--|----|------------------------------------|----|
| | Yes | No | Yes | No | Yes | No | Yes | No |
| Joachim Kreuzburg CEO and Chairman | [1] | | [3] | | 2,400 | | 600 | |

[1] Joachim Kreuzburg has a service contract (without social security components) with Sartorius AG for his duties performed as members of the Executive Board of the major shareholder Sartorius AG. This is standard practice in Germany. The contract includes a cap regarding potential severance payments at the maximum of a two years annual remuneration. Furthermore there is a post-contractual non-competition clause obligation, that will last for two years after an Executive Board member has left the Group. During this time, if the non-competition clause is not waived or terminated, this Executive Board member may claim half of his most recent annual remuneration received from the company.

[3] Additionally there is a general pension plan in place at the Sartorius AG level for Joachim Kreuzburg. The level of their entitlement to benefits paid under a company pension plan depends on their respective tenure

Independent Auditors' Fees

Principal Independent Auditors

KPMG S.A.

480, avenue du Prado – CS 90021 – 13272 Marseille Cedex 08 – France

Represented by John Evans.

First commissioned by the Annual General Shareholders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General Shareholders' Meeting to approve the 2020 financial statements.

Member of the Compagnie régionale de Versailles.

Deloitte et Associés

10, Place de la Joliette – Les Docks – Atrium 10.4 – BP 64529 - 13567 Marseille Cedex 02 – France

Represented by Vincent Gros.

First commissioned by the Annual General Shareholders' Meeting on 19 May 2006.

Date commission expires: 2024 Annual General Shareholders' Meeting to approve the 2023 financial statements.

Member of the Compagnie régionale de Versailles.

Independent Auditors' Fees

| € in K | KPMG | | | | Deloitte | | | |
|--|------------|---------------|------------|---------------|------------|---------------|------------|--------------|
| | 2018 | | 2017 | | 2018 | | 2017 | |
| Audit | | | | | | | | |
| Independent audit, certification, parent company & consolidated financial statements | | | | | | | | |
| Parent company | 61 | 7.4% | 76 | 9.3% | 54 | 37.2% | 58 | 35.2% |
| Subsidiaries | 754 | 92.0% | 753 | 91.9% | 91 | 62.8% | 97 | 58.8% |
| Services directly related to audit services | | | | | | | | |
| Parent company | | | | | | | | |
| Subsidiaries | | | | | | | | |
| Subtotal | 815 | 99.5% | 829 | 101.1% | 145 | 100.0% | 155 | 93.9% |
| Other services | | | | | | | | |
| Legal, tax, corporate | 5 | 0.5% | 47 | 5.8% | 0 | 0.0% | 0 | 0.0% |
| Information technology, other | 0 | 0.0% | 0 | 0.0% | 0 | 0.0% | 10 | 6.1% |
| Subtotal | 5 | 0.5% | 47 | 5.8% | 0 | 0.0% | 10 | 6.1% |
| Total | 820 | 100.0% | 876 | 107% | 145 | 100.0% | 165 | 100% |

Substitute Independent Auditors

SALUSTRO REYDEL

3, cours du Triangle – Immeuble Le Palatin – 92939
Paris-La Défense Cedex – France

Represented by Jean-Claude Reydel.

First commissioned by the Annual General Shareholders' Meeting on 7 April 2015.

Date commission expires: 2021 Annual General Shareholders' Meeting to approve the 2020 financial statements.

Member of the Compagnie régionale de Versailles.

| Ernst & Young | | | | Other | | | | Total | | | |
|---------------|--|------|--|-------|--------|------|-------|-------|--------|-------|-------|
| 2018 | | 2017 | | 2018 | | 2017 | | 2018 | | 2017 | |
| | | | | | | | | | | | |
| | | | | | | | | | | | |
| | | | | | | | | 115 | 10.2% | 134 | 10.4% |
| | | | | 60 | 37.8% | 88 | 35.7% | 905 | 80.6% | 938 | 72.7% |
| | | | | | | | | | | | |
| | | | | | | | | | | | |
| | | | | | | | | | | | |
| | | 0 | | 60 | 37.8% | 88 | 35.5% | 1,020 | 90.8% | 1072 | 83.1% |
| | | | | | | | | | | | |
| | | 1 | | 51 | 32.4% | 140 | 56.5% | 56 | 5.0% | 188 | 14.6% |
| | | 0 | | 47 | 29.7% | 20 | 8.0% | 47 | 4.2% | 30 | 2.3% |
| | | 1 | | 98 | 62.2% | 160 | 64.5% | 103 | 9.2% | 218 | 16.9% |
| | | 1 | | 158 | 100.0% | 248 | 100% | 1,123 | 100.0% | 1,291 | 100% |

Consolidated Financial Statements and Notes

04

Statement of Profit or Loss and Other Comprehensive Income

| | Notes | 2018 12 months € in K | 2017 12 months € in K |
|--|-------|-----------------------------|-----------------------------|
| Sales revenue | [9] | 1,212,152 | 1,081,033 |
| Cost of sales | | -582,589 | -526,208 |
| Gross profit on sales | | 629,563 | 554,825 |
| Selling and distribution costs | | -215,208 | -195,241 |
| Research and development costs | | -60,616 | -53,210 |
| General administrative expenses | | -67,004 | -61,685 |
| Other operating income and expenses | [11] | 13,500 | -22,987 |
| Earnings before interest and taxes (EBIT) | | 300,234 | 221,702 |
| Financial income | [12] | 5,349 | 9,500 |
| Financial expenses | [12] | -21,035 | -10,589 |
| Financial result | | -15,685 | -1,089 |
| Profit before tax | | 284,549 | 220,613 |
| Income taxes | [13] | -74,590 | -56,849 |
| Net profit for the period | | 209,959 | 163,763 |
| Attributable to: | | | |
| Equity holders of Sartorius Stedim Biotech | | 208,052 | 161,088 |
| Non-controlling interest | [22] | 1,907 | 2,676 |
| Earnings per share (€) | [15] | 2.26 | 1.75 |
| Diluted earnings per share (€) | [15] | 2.26 | 1.75 |

Other Comprehensive Income

| | Notes | 2018 12 months € in K | 2017 12 months € in K |
|---|-------|-----------------------------|-----------------------------|
| Net profit for the period | | 209,959 | 163,763 |
| Cash flow hedges | [36] | -11,547 | 26,299 |
| of which effective portion of changes in fair value | | -20,016 | 26,901 |
| of which reclassified to profit or loss | | 8,469 | -602 |
| Income tax on cash flow hedges | [19] | 3,463 | -7,889 |
| Foreign currency translation differences | | 6,433 | -27,034 |
| Items that are or may be reclassified subsequently to profit or loss | | -1,651 | -8,624 |
| Remeasurements of the net defined benefit liabilities | [23] | 469 | -667 |
| Income tax on remeasurements of the net defined benefits liabilities | [19] | -160 | 481 |
| Items that will not be reclassified to profit or loss | | 309 | -186 |
| Other comprehensive income after tax | | -1,342 | -8,810 |
| Total comprehensive income | | 208,617 | 154,953 |
| Attributable to: | | | |
| Equity holders of Sartorius Stedim Biotech | | 206,673 | 152,335 |
| Non-controlling interest | | 1,943 | 2,619 |

Statement of Financial Position

| | Notes | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|---|-------|-------------------------|-------------------------|
| Non-current assets | | | |
| Goodwill | [16] | 384,695 | 386,045 |
| Other Intangible Assets | [16] | 177,492 | 172,906 |
| Property, plant and equipment | [17] | 435,980 | 331,295 |
| Financial Assets | | 5,644 | 11,808 |
| Other Assets | | 632 | 18 |
| Deferred tax assets | [19] | 14,490 | 10,989 |
| | | 1,018,932 | 913,060 |
| Current assets | | | |
| Inventories | [20] | 252,002 | 185,604 |
| Trade receivables | [28] | 220,231 | 211,616 |
| Other financial assets | [29] | 22,036 | 20,049 |
| Current tax assets | | 12,950 | 24,465 |
| Other assets | | 21,334 | 16,561 |
| Cash and cash equivalents | [27] | 23,975 | 32,552 |
| | | 552,529 | 490,845 |
| Total assets | | 1,571,461 | 1,403,905 |
| Equity | | | |
| Equity attributable to SSB S.A. shareholders | | 1,036,398 | 872,028 |
| Issued capital | [21] | 18,436 | 18,436 |
| Capital reserves | | 231,526 | 231,526 |
| Retained earnings (including net profit) | | 786,436 | 622,066 |
| Non-controlling interest | [22] | 8,476 | 7,426 |
| | | 1,044,874 | 879,454 |
| Non-current liabilities | | | |
| Pension provisions | [23] | 35,595 | 35,446 |
| Other provisions | [24] | 2,877 | 3,103 |
| Loans and borrowings | [30] | 43,125 | 46,250 |
| Finance lease liabilities | | 14,976 | 15,829 |
| Other financial liabilities | [31] | 8,889 | 30,392 |
| Deferred tax liabilities | [19] | 39,150 | 42,987 |
| | | 144,612 | 174,007 |
| Current liabilities | | | |
| Provisions | [24] | 12,283 | 9,558 |
| Trade payables | [32] | 154,568 | 117,959 |
| Loans and borrowings | [30] | 89,817 | 95,933 |
| Finance lease liabilities | | 1,717 | 1,658 |
| Other financial liabilities | [33] | 14,312 | 34,964 |
| Employee benefits | | 39,335 | 31,895 |
| Current tax liabilities | | 33,070 | 31,791 |
| Other liabilities | | 36,873 | 26,686 |
| | | 381,975 | 350,444 |
| Total equity and liabilities | | 1,571,461 | 1,403,905 |

Statement of Cash Flows

| | Notes | 2018 12 months € in K | 2017 12 months € in K |
|--|--------------|-----------------------------|-----------------------------|
| Profit before tax | | 284,549 | 220,613 |
| Financial result | [12] | 15,685 | 1,089 |
| Earnings before interest and taxes (EBIT) | | 300,234 | 221,702 |
| Depreciation amortization of fixed assets | [16][17] | 60,914 | 50,678 |
| Increase decrease in provisions | [23][24] | 2,361 | 655 |
| Income taxes paid | [13] | -65,537 | -49,248 |
| Other non-cash items | | -32,909 | 3,489 |
| Gross cash flows from operating activities | | 265,062 | 227,276 |
| Increase decrease in receivables and other assets | [28][29] | -21,604 | -46,316 |
| Increase decrease in inventories | [20] | -64,962 | -21,091 |
| Increase decrease in liabilities | [31][32][33] | 48,793 | 14,819 |
| Net cash flow from operating activities | | 227,289 | 174,689 |
| Acquisitions of intangible and tangible assets | [16][17] | -176,540 | -126,826 |
| Net cash flow from investing activities | | -176,540 | -126,826 |
| Payments for acquisitions of consolidated subsidiaries and other business operations; net of cash acquired | [8] | 0 | -68,100 |
| Net cash flow from investing activities and acquisitions | | -176,540 | -194,926 |
| Interest received | [12] | 2,569 | 290 |
| Interest paid and other financial charges | [12] | -8,995 | -1,914 |
| Dividends paid to: | | | |
| - Shareholders of Sartorius Stedim Biotech SA | | -42,403 | -38,713 |
| - Non-controlling interest | | -783 | -704 |
| Gross cash flows from financing activities | | -49,612 | -41,041 |
| Proceeds from loans and borrowings raised | [6][30] | 12,829 | 60,994 |
| Repayments of loans and borrowings | [6][30] | -22,811 | -3,333 |
| Net cash flow from financing activities | | -59,594 | 16,620 |
| Net increase decrease in cash and cash equivalents | | -8,845 | -3,617 |
| Cash and cash equivalents at the beginning of the period | | 32,552 | 34,756 |
| Net effect of currency translation on cash and cash equivalents | | 268 | 1,414 |
| Cash and cash equivalents at the end of the period | | 23,975 | 32,552 |

The Notes to the Consolidated Financial Statements are an integral part of these statements.

Statement of Changes in Equity

| € in K | Issued capital | Capital reserves | Hedging reserves | Pension reserves | Retained earnings | Foreign currency translation reserves | Group equity | Non-controlling interest | Total equity |
|---|----------------|------------------|------------------|------------------|-------------------|---------------------------------------|------------------|--------------------------|------------------|
| Balance at Jan. 1, 2017 | 18,083 | 231,526 | -6,961 | -10,983 | 490,275 | 36,064 | 758,004 | 5,551 | 763,555 |
| Net profit for the period | 0 | 0 | 0 | 0 | 161,088 | 0 | 161,088 | 2,676 | 163,763 |
| Cash flow hedges | 0 | 0 | 26,299 | 0 | 0 | 0 | 26,299 | 0 | 26,299 |
| Remeasurements of the net defined benefit liabilities | 0 | 0 | 0 | -667 | 0 | 0 | -667 | 0 | -667 |
| Foreign currency translation differences | 0 | 0 | 0 | 0 | 0 | -26,977 | -26,977 | -57 | -27,034 |
| Deferred taxes | 0 | 0 | -7,889 | 481 | 0 | 0 | -7,408 | 0 | -7,408 |
| Other comprehensive income for the period | 0 | 0 | 18,410 | -186 | 0 | -26,977 | -8,753 | -57 | -8,810 |
| Total comprehensive income | 0 | 0 | 18,410 | -186 | 161,088 | -26,977 | 152,335 | 2,619 | 154,953 |
| Dividends | 0 | 0 | 0 | 0 | -38,713 | 0 | -38,713 | -704 | -39,417 |
| Changes in non-controlling interest | 0 | 0 | 0 | 0 | 40 | 0 | 40 | -40 | 0 |
| Other changes | 353 | 0 | 0 | 0 | 9 | 0 | 362 | 0 | 362 |
| Balance at Dec. 31, 2017 | 18,436 | 231,526 | 11,449 | -11,169 | 612,699 | 9,087 | 872,027 | 7,426 | 879,453 |
| Adjustment on adoption of IFRS 9 | | | | | 321 | | 321 | | 321 |
| Balance at Jan. 1, 2018 | 18,436 | 231,526 | 11,449 | -11,169 | 613,020 | 9,087 | 872,349 | 7,426 | 879,775 |
| Net profit for the period | 0 | 0 | 0 | 0 | 208,052 | 0 | 208,052 | 1,907 | 209,959 |
| Cash flow hedges | 0 | 0 | -11,547 | 0 | 0 | 0 | -11,547 | 0 | -11,547 |
| Remeasurements of the net defined benefit liabilities | 0 | 0 | 0 | 469 | 0 | 0 | 469 | 0 | 469 |
| Foreign currency translation differences | 0 | 0 | 0 | 0 | 0 | 6,396 | 6,396 | 37 | 6,433 |
| Deferred taxes | 0 | 0 | 3,463 | -160 | 0 | 0 | 3,303 | 0 | 3,303 |
| Other comprehensive income for the period | 0 | 0 | -8,084 | 309 | 0 | 6,396 | -1,379 | 37 | -1,342 |
| Total comprehensive income | 0 | 0 | -8,084 | 309 | 208,052 | 6,396 | 206,673 | 1,943 | 208,617 |
| Dividends | 0 | 0 | 0 | 0 | -42,403 | 0 | -42,403 | -783 | -43,186 |
| Other changes | | 0 | 0 | 0 | -221 | 0 | -221 | -110 | -331 |
| Balance at December 31, 2018 | 18,436 | 231,526 | 3,365 | -10,860 | 778,448 | 15,483 | 1,036,398 | 8,476 | 1,044,874 |

Notes to the Financial Statements

1. General Information

Sartorius Stedim Biotech is a leading international supplier of products and services that enable the biopharmaceutical industry to develop and manufacture drugs safely and efficiently. As a total solutions provider, Sartorius Stedim Biotech offers a portfolio covering nearly all steps of biopharmaceutical manufacture. The company focuses on single-use technologies and value-added services to meet the rapidly changing technology requirements of the industry it serves. With its own manufacturing and R&D sites in Europe, North America and Asia and an international network of sales companies, Sartorius Stedim Biotech has a global reach.

Headquartered in Aubagne, France, Sartorius Stedim Biotech S.A. is listed on the Euronext Paris (ISIN code: FR 0000053266).

Sartorius Stedim Biotech S.A.'s ultimate parent company is Sartorius AG, headquartered in Goettingen, Germany, and listed at several German stock exchanges (ISIN codes: 0007165607 ordinary shares, 0007165631 preference shares).

In compliance with the European Regulation 1606/2002 of July 19, 2002, requiring listed companies to use International Accounting Standards, the consolidated financial statements of the Sartorius Stedim Biotech Group for the year ended December 31, 2018, are compliant with the Standards and Interpretations IFRS and IFRIC of the IASB as adopted by the European Union, which are available at the following site:

https://ec.europa.eu/commission/index_fr.

The consolidated financial statements are prepared in euros. Unless otherwise specified, all amounts are disclosed in thousands of euros (abbreviated as € in K). In some cases, the sum of the figures given in this report may not precisely equal the stated totals and percentages may not be exact due to rounding.

These consolidated financial statements were approved by the Board of Directors on February 7, 2019 and will be submitted for approval by the Shareholders' Meeting on March 26, 2019.

2. Effects of New Financial Reporting Standards

The following major new accounting rules were applicable for the first time to the present consolidated financial statements of the Group:

– IFRS 9, Financial Instruments

This standard changes the accounting for financial instruments. For detailed information on these changes and the initial application of IFRS 9, see note 26.

– IFRS 15, Revenue from Contracts with Customers

– Clarifications to IFRS 15, Revenue from Contracts with Customers

This standard defines when and in which amount revenue is recognized. For detailed information on the related changes and the initial application of IFRS 15, see note 9.

The following new accounting rules were applicable for the first time to the present financial statements and had no material impact on the presentation of the company's financial position and financial performance:

– Annual Improvements to IFRSs 2014 - 2016 Cycle, Amendments to IFRS 1 and IAS 28

The amendments are related to the first-time application of IFRS (IFRS 1) as well as the option for venture capital organizations and similar organizations to measure investments in associates or joint ventures at fair value through profit or loss (IAS 28).

– Amendments to IFRS 2, Classification and Measurement of Share-based Payment Transactions

The amendments include clarifications with regard to the accounting for share-based payment transactions.

– Amendments to IAS 40, Transfers of Investment Property

The amendments intend to clarify the guidance regarding transfers to, or from, investment properties.

– IFRIC 22, Foreign Currency Transactions and Advance Consideration

The interpretation addresses a question in relation to IAS 21. It clarifies which exchange rate is used for foreign currency transactions that involve advance payments received or made.

The following standards, interpretations and amendments were not yet applied to the consolidated financial statements of the reporting year as they had not yet been adopted by the EU or their application was not mandatory for 2018:

| Standard Interpretation | Title | Applicable for financial years from ¹⁾ | Endorsement by the EU Commission |
|--|---|---|----------------------------------|
| IFRS 14 | Regulatory Deferral Accounts | January 1, 2016 | No |
| IFRS 16 | Leases | January 1, 2019 | Yes |
| IFRIC 23 | Uncertainty over Income Tax Treatments | January 1, 2019 | Yes |
| Amendments to IFRS 9 | Prepayment Features with Negative Compensation | January 1, 2019 | Yes |
| Amendments to IFRS 28 | Long-term Interests in Associates and Joint Ventures | January 1, 2019 | No |
| Amendments to IFRS 3, IFRS 11, IAS 12 and IAS 23 | Annual Improvements to IFRSs 2015 - 2017 Cycle (issued in Dec. 2017) | January 1, 2019 | No |
| Amendments to IAS 19 | Plan Amendment, Curtailment or Settlement | January 1, 2019 | No |
| Amendments to IAS 1 and IAS 8 | Definition of Material | January 1, 2020 | No |
| Various Standards / Framework | Amendments to References to the Conceptual Framework in IFRS Standards | January 1, 2020 | No |
| Amendments to IFRS 3 | Definition of a Business | January 1, 2020 | No |
| IFRS 17 | Insurance Contracts | January 1, 2021 | No |
| Amendments to IFRS 10 and IAS 28 | Sale or Contribution of Assets between an Investor and its Associate or Joint Venture | n/a | No |

¹⁾ These are required to be applied once they are endorsed by the EU Commission.

The dates mentioned above are those required by the standard themselves (IASB effective dates).

IFRS 16, Leases

The Group did not yet apply IFRS 16, Leases. The application of the standard is mandatory for periods from 2019 onwards. IFRS 16 introduces a standardized accounting model according to which leases are generally to be recognized on the lessee's balance sheet.

A lessee recognizes a right-of-use asset representing its right to use a lease asset, as well as a liability resulting from the lease, which represents its obligation to make lease payments. There are exemptions for short-term leases and leases of low-value assets. Accounting for the lessor is comparable to that of the current standard; i.e., lessors continue to classify leases as financial or operating leases.

The Group has analyzed the effects of this new standard on the consolidated financial statements during the year 2018. The Group did not apply the standard early. The Group does not plan to apply the standard retrospectively in accordance with IAS 8, but rather plans to recognize the effect from initially applying IFRS 16 in retained earnings at the date of

initial application. The Group will use the exemptions for short-term leases and leases of low-value assets and recognize the corresponding lease payments as an expense generally on a straight-line basis over the particular lease term.

IFRS 16 will lead to an increase in fixed assets and financial liabilities. Based on its present level of knowledge, the Group does not expect any significant impact overall on its most important key figures, such as equity ratio or underlying EBITDA.

For example, on the basis of a differentiated analysis of the operating lease commitments (see also note 18) conducted in the course of the preparation of the present consolidated financial statements, total assets would increase by approx. €35 million as of December 31, 2018. This would result in a reduction of the equity ratio by approx. 1.5%. On this basis, the EBITDA margin would increase by about 0.5 - 1% for the year 2018.

3. Significant Accounting Policies

Basis of Preparation

The consolidated financial statements of the Group are based on the principle of the historical cost of acquisition, construction or production, with the exception of the items reflected at fair value, such as derivative financial instruments.

Consolidation

The consolidated financial statements of the Sartorius Stedim Biotech Group include the annual financial statements of all companies, which are controlled directly or indirectly by Sartorius Stedim Biotech S.A. In terms of IFRS 10, Consolidated Financial Statements, the Group Sartorius Stedim Biotech controls an entity when it is exposed to, or has rights to, variable returns from its involvement with the entity and has the ability to affect those returns through its power over the entity.

Such entities are included in the consolidated financial statements from the time when Sartorius Stedim Biotech S.A. or its subsidiaries obtain such control until the date on which control ceases.

Subsidiaries have been included on the basis of their annual financial statements for the same reporting period as the parent company, using uniform Group recognition and measurement methods.

All intragroup assets and liabilities, equity, income, expenses and cash flows relating to transactions between members of the Group are eliminated on consolidation.

Business Combinations

Business combinations are accounted according to the acquisition method. The identifiable acquired assets and assumed liabilities are generally recorded at fair value on the date of acquisition.

For significant acquisitions, the purchase price allocation is carried out with assistance from independent third-party valuation specialists. The valuations are based on the information available at the acquisition date.

The Group determines goodwill at the acquisition date as:

- the fair value of the consideration transferred; and
- the amount recognized for any non-controlling interest in the acquiree; and
- if the business combination is carried out in stages, the fair value of any previously held equity interest in the acquiree; less
- the net recognized amount for the identifiable assets acquired and liabilities assumed.

When the difference is negative, the purchase gain is recognized immediately in income. Expenses directly related to business combinations are recorded in the profit or loss as they are incurred.

Foreign Currency Transactions

The presentation currency of the consolidated financial statements of the Sartorius Stedim Biotech Group is the euro (financial statements presented in thousands of euros). In the financial statements of each company, transactions denominated in foreign currencies have been translated into the functional currency of the subsidiary at the exchange rate applicable on the date of the transaction. Monetary assets and debts denominated in a foreign currency have been translated at the exchange rate on the balance sheet date. Rate gains and losses have been recognized in profit or loss for the period.

Translation of financial statements prepared in foreign currencies

Subsidiaries' financial statements prepared in foreign currencies have been translated pursuant to IAS 21, The Effects of Changes in Foreign Exchange Rates, in accordance with the concept of a functional currency. Foreign subsidiaries have been regarded as independent subdivisions of the Sartorius Stedim Biotech Group. The assets (including goodwill) and liabilities of the entities that have a functional currency different from the presentation currency are translated at the exchange rate prevailing at the balance sheet date. The incomes, expenses, and cash flows of these entities have been translated using the average rate for the year, to the extent that this rate represents an approximate value of exchange rates used as of the date of the transaction in the absence of significant fluctuations. Resulting translation differences are recognized in other comprehensive income.

For long-term loans for which settlement is neither planned nor likely in the foreseeable future, the Group applies the principle of "net investment in a foreign operation." Exchange differences resulting from these loans are recognized in other comprehensive income in accordance with IAS 21.32.

The exchange rates for major currencies against the euro were considered as follows:

| For 1 € | Year-end exchange rates | | Average exchange rates | |
|---------|-------------------------|------------|------------------------|------------|
| | 2018 | 2017 | 2018 | 2017 |
| USD | 1.14500 | 1.19930 | 1.18129 | 1.12955 |
| GBP | 0.89453 | 0.88723 | 0.88465 | 0.87670 |
| JPY | 125.85000 | 135.01000 | 130.40242 | 126.70218 |
| CHF | 1.12690 | 1.17020 | 1.15517 | 1.11173 |
| INR | 79.72980 | 76.60550 | 80.69389 | 73.52751 |
| KRW | 1277.93000 | 1279.61000 | 1299.12464 | 1276.62397 |
| CNY | 7.87510 | 7.80440 | 7.80666 | 7.62790 |

4. Use of Judgments and Estimates

During the preparation of consolidated financial statements, management uses estimates and assumptions based on their best knowledge of the current and future situation. However, actual results may differ from these estimates. These estimates and assumptions are revised on a regular basis, and the impact of changes in estimates is recognized prospectively.

In addition, Group management exercises its judgment in defining the accounting treatment of specific transactions when the existing Standards and Interpretations do not specifically treat the accounting problems concerned.

Assumptions and estimates primarily concern the following topics:

Business Combinations

The accounting for business combinations requires that the assets acquired and liabilities assumed be recorded at their respective fair values on the date the Group obtains control. The application of the acquisition method requires estimates and assumptions to be made, especially concerning the fair values of the acquired intangible assets, property, plant and equipment and the liabilities assumed at the acquisition date, and the useful lives of the acquired intangible assets, property, plant and equipment.

These measurements are based to a large extent on anticipated cash flows. If actual cash flows vary from those used in calculating fair values, this may materially affect the Group's future results of operations.

Impairment of Assets

The book values (carrying amounts) of property, plant and equipment and intangible assets are subject to impairment testing if there is an indication of impairment and at least once a year for assets with an indefinite useful life or not yet available for use in accordance with IAS 36, Impairment of Assets. When an asset is tested, the recoverable amount of the asset is estimated. The recoverable amount of an asset or a cash-generating unit (CGU) is the higher of its fair value – less costs to sell the asset or its CGU – and its value in use. In the event the individual asset's recoverable amount cannot be estimated, the recoverable amount of the asset's CGU is estimated.

If the estimated recoverable amount of an asset (or a CGU) goes below its book value (carrying amount), this carrying amount is reduced to the recoverable amount (impairment allocated in priority to goodwill).

If the causes of the asset impairment are removed, the book value of the asset (or the CGU) is credited to the newly estimated recoverable amount. However, the book value increase is limited to the value that the asset (or CGU) would have had if no asset impairment loss had been recognized in previous financial years.

The calculation of the value in use is generally based on discounted cash flow methods using cash flow projections up to five years. These projections take into account past experience and represent management's best estimate about future sales revenue and cost developments. Cash flows after the planning period are extrapolated using individual growth rates. Key assumptions on which management has based its determination of the value in use include estimated growth rates, weighted average cost of capital and tax rates. These estimates can have a material impact on the respective values and ultimately the amount of any impairment.

Intangible Assets

The capitalization of self-constructed intangible assets also includes a significant level of judgment, e.g. the evaluation of feasibility of a development project, the expected market prospects and the determination of useful lives.

Employee Benefits – Pension Provisions

Obligations for pension and other post-employment benefits are determined in accordance with actuarial valuations. These valuations rely on key assumptions including discount rates, expected salary increases and mortality rates. The discount rate assumptions are determined by reference to yields on high-quality corporate bonds of appropriate duration and currency at the end of the reporting period.

Due to changing market and economic conditions the underlying key assumptions may differ from actual developments and may lead to significant changes in pension and other post-employment benefit obligations.

Such differences are recognized in other comprehensive income in the period in which they occur. For a sensitivity analysis, see note 23, Pension and Employee Benefits Provisions.

Provisions, Contingent Liabilities and Contingent Assets

Provisions are recognized for legal or constructive obligations that exist as of the balance sheet date. To determine the amount of the obligations, certain estimates and assumptions have to be applied, including the determination of the probability and the amount of future outflows of resources. Typically, significant estimates are involved in the determination of provisions related to onerous contracts, warranty costs, asset retirement obligations and legal proceedings.

Income Taxes

The Group operates in various tax jurisdictions and therefore has to determine tax positions under respective local tax laws and tax authorities' views which can be complex and subject to different interpretations of taxpayers and local tax authorities. Deferred tax assets have to be recognized for all deductible temporary differences and unused tax losses to the extent that it is probable that future taxable profit will be available against which the deductible temporary differences and unused tax losses can be utilized. As future developments are uncertain and partly beyond management's control,

assumptions are necessary to estimate future taxable profits as well as the period in which deferred tax assets will be recovered.

Estimates are revised in the period in which there is sufficient evidence to revise the assumption. If management considers it probable that all or a portion of a deferred tax asset cannot be realized, the corresponding amount is not recorded as an asset.

Fair Value Measurement

A number of the Group's accounting policies and disclosures may require the measurement of fair values, for both financial and non-financial assets and liabilities, including Level 3 fair values (unobservable inputs).

If third party information, such as broker quotes or pricing services, is used to measure fair values, then management assesses the evidence obtained from the third parties to support the conclusion that these valuations meet the requirements of IFRS, including the level in the fair value hierarchy in which the valuations should be classified.

When measuring the fair value of an asset or a liability, the Group uses observable market data as far as possible. If the inputs used to measure the fair value of an asset or a liability fall into different levels of the fair value hierarchy, then the fair value measurement is categorized in its entirety in the same level of the fair value hierarchy as the lowest level input that is significant to the entire measurement.

5. Operating Segments

According to IFRS 8, Operating Segments the identification of reportable operating segments is based on the "management approach"; i.e. the segments are defined analogously to the internal financial reporting of an entity. Therefore, an area of activity is to be considered an operating segment if its business activities may result in revenues and expenses, its operating results are regularly reviewed by the entity's chief operating decision maker (= the Executive Members of the Board of Directors) and discrete financial information is available in its internal reporting. Internal control and reporting within Sartorius Stedim Biotech is based on the approach of operating as a "total solution provider" for its customers. Accordingly, there is only one single segment to be identified for Sartorius Stedim Biotech, driven by the product and customer perspective: Biopharm.

The key performance indicator of the operating segment of the Sartorius Stedim Biotech Group is the so-called "underlying EBITDA", as the board monitors this performance measure at a consolidated level and they believe this measure is relevant to an understanding of the Group's financial performance.

EBITDA corresponds to earnings before interest, taxes, depreciation and amortization; "underlying EBITDA" means EBITDA adjusted for extraordinary items. In this connection, extraordinary items are expenses and income that are of an exceptional or a one-time nature and accordingly distort the sustainable profitability of a segment and have a material impact on the net worth, financial position and earnings of the Group. Examples of such items are restructuring

expenses, large Group projects as well as proceeds or losses from the disposal, sale or other transfer of financial assets or of property, plant and equipment, provided that these are not of a recurrent nature.

Underlying EBITDA is not a defined performance measure in IFRS. The Group's definition of underlying EBITDA may not be comparable with similarly titled performance measures and disclosures by other entities.

Segment assets and segment liabilities are not analyzed on a regular basis to the chief operating decision maker and are therefore not part of the segment report.

| € in K | 2018 | 2017 | Biopharm Change | 2018 | 2017 | Group Change |
|--|-----------|-----------|--------------------|-----------|-----------|-----------------|
| Sales revenue | 1,212,152 | 1,081,033 | 12% | 1,212,152 | 1,081,033 | 12% |
| Underlying EBITDA | 342,430 | 294,912 | 16% | 342,430 | 294,912 | 16% |
| as a % of sales revenue | 28.2% | 27.3% | | 28.2% | 27.3% | |
| EBIT | 300,234 | 221,702 | 35% | 300,234 | 221,702 | 35% |
| as a % of sales revenue | 24.8% | 20.5% | | 24.8% | 20.5% | |
| Acquisitions of intangible and tangible assets | 176,945 | 136,734 | 29% | 176,945 | 136,734 | 29% |
| as a % of sales revenue | 14.6% | 12.6% | | 14.6% | 12.6% | |

Reconciliation of Segment Profit or Loss:

| € in K | 2018 | 2017 |
|----------------------------------|----------------|----------------|
| Underlying EBITDA of the segment | 342,430 | 294,912 |
| Depreciation and amortization | -54,918 | -50,645 |
| Extraordinary effects | 12,722 | -22,564 |
| EBIT | 300,234 | 221,702 |
| Financial result | -15,685 | -1,089 |
| Profit before tax | 284,549 | 220,613 |

Supplementary Information by Region

To provide additional information required by IFRS 8, the table below presents the supplementary information by geographical region. The key figures for non-current assets of the geographical areas refer to the company location, whereas sales revenue is reported according to the customer's location.

The non-current assets correspond to property, plant and equipment as well as to intangible assets (including goodwill) of the Group affiliates that are to be allocated to these various regions.

The amount of sales revenue with a single customer does not exceed 5% of the consolidated sales revenue (2018 and 2017).

| € in K | Sales revenue | | Non-current assets | |
|---------------------|------------------|------------------|--------------------|----------------|
| | 2018 | 2017 | 2018 | 2017 |
| EMEA | 508,241 | 460,664 | 816,639 | 756,242 |
| thereof Germany | 144,602 | 130,454 | 340,599 | 296,567 |
| thereof France | 54,977 | 56,955 | 331,371 | 305,508 |
| Americas | 422,625 | 368,817 | 159,484 | 113,722 |
| thereof USA | 396,509 | 339,319 | 159,484 | 113,722 |
| Asia Pacific | 281,286 | 251,552 | 22,044 | 20,281 |
| thereof China | 83,904 | 66,978 | 1,474 | 1,526 |
| thereof South Korea | 72,824 | 76,585 | 7,502 | 7,109 |
| Group | 1,212,152 | 1,081,033 | 998,167 | 890,245 |

6. Statement of Cash Flows

The statement of cash flows shows the impact of cash inflows and outflows on the cash and cash equivalents of the Group. The cash flows are classified by operating, investing and financing activities according to IAS 7 (Statement of Cash Flows).

In this context cash equivalents are assets that can be converted into cash within a short maturity (generally less than three months). The amount considered in the

statement of cash flows is equal to the amount of cash and cash equivalents in the statement of financial position.

The impact resulting from the modified agreement with Lonza in the field of cell culture media posted in 2018 had no impact on the statement of cash flows (see note 11).

The following table summarizes the development of the liabilities arising from financing activities during the reporting period:

| | Balance at Dec. 31, 2016 € in K | cash flows | Currency effects | other non-cash changes | Balance at Dec. 31, 2017 € in K |
|--|---------------------------------------|---------------|---------------------|---------------------------|---------------------------------------|
| Loans and borrowings | 84,052 | 58,347 | -215 | 0 | 142,184 |
| Finance lease liabilities | 18,270 | -686 | -173 | 76 | 17,487 |
| Liability for phantom units in connection with the AllPure acquisition | 5,833 | 0 | -706 | 155 | 5,282 |
| Total financial liabilities from financing activities | 108,155 | 57,661 | -1,094 | 231 | 164,952 |

| | Balance at Dec. 31, 2017 € in K | cash flows | Currency effects | other non-cash changes | Balance at Dec. 31, 2018 € in K |
|--|---------------------------------------|---------------|---------------------|---------------------------|---------------------------------------|
| Loans and borrowings | 142,184 | -9,229 | -12 | 0 | 132,943 |
| Finance lease liabilities | 17,487 | -754 | -40 | 0 | 16,693 |
| Liability for phantom units in connection with the AllPure acquisition | 5,282 | 0 | 250 | 3,206 | 8,739 |
| Total financial liabilities from financing activities | 164,953 | -9,982 | 198 | 3,206 | 158,375 |

7. Scope of Consolidation

The 2018 financial statements of the following entities:

- TAP Biosystems (PHC) Ltd., UK
- TAP Biosystems Ltd., UK
- Distribo GmbH, Germany

were not included in the scope of consolidation, because the figures were of minor importance for assessing the financial position of the Group.

The sales revenue and total assets of the non-consolidated companies are below 1% of the Group figures.

The financial statements of the following companies have been included in the Group financial statements:

| | Ownership in % |
|--|-------------------|
| EMEA | |
| Sartorius Stedim Biotech S.A., Aubagne, France | Parent company |
| Sartorius Stedim Belgium N.V., Brussels, Belgium | 100 |
| Sartorius Stedim Nordics Oy, Helsinki, Finland | 100 |
| Sartorius Stedim Biotech GmbH, Goettingen, Germany | 100 |
| Sartorius Stedim Plastics GmbH, Goettingen, Germany | 100 |
| Sartorius Stedim North America Holding GmbH, Goettingen, Germany | 100 |
| Sartorius Stedim Systems GmbH, Guxhagen, Germany | 100 |
| Sartorius Stedim Cellca GmbH, Laupheim, Germany | 100 |
| Sartorius Stedim UK Ltd., Epsom, UK | 100 |
| Sartorius Stedim BioOutsource Ltd., Glasgow, UK | 100 |
| Sartorius Stedim Lab Ltd., Stonehouse, UK | 100 |
| TAP Biosystems Group Ltd., Royston, UK | 100 |
| TAP ESOP Management Ltd., Royston, UK | 100 |
| The Automation Partnership Cambridge Ltd., Royston, UK | 100 |
| Sartorius Stedim FMT S.A.S., Aubagne, France | 100 |
| Sartorius Stedim France S.A.S., Aubagne, France | 100 |
| Sartorius Stedim Aseptics S.A., Lourdes, France | 100 |
| Sartorius Stedim Ireland Ltd., Dublin, Ireland | 100 |
| Sartorius Stedim Italy S.p.A., Florence, Italy | 100 |
| Sartorius Stedim Netherlands B.V., Amersfoort, Netherlands | 100 |
| Sartorius Stedim Austria GmbH, Vienna, Austria | 100 |
| Sartorius Stedim Poland sp. z.o.o., Kostrzyn, Poland | 100 |
| Sartorius Stedim RUS, St. Petersburg, Russia | 100 |
| Sartorius Stedim Data Analytics AB, Umeå, Sweden | 100 |
| Sartorius Stedim Switzerland AG, Tagelswangen, Switzerland | 100 |
| Sartorius Stedim Spain S.A., Madrid, Spain | 100 |
| Sartorius Stedim Hungaria Kft., Budapest, Hungary | 100 |
| Sartorius Stedim Bioprocess S.A.R.L., M'Hamdia, Tunisia | 100 |
| Americas | |
| Sartorius Stedim Filters Inc., Yauco, Puerto Rico | 100 |
| Sartorius Stedim North America Inc., Wilmington, USA | 100 |
| Asia Pacific | |
| Sartorius Stedim Australia Pty. Ltd., Dandenong South, Victoria, Australia | 100 |
| Sartorius Stedim Biotech (Beijing) Co. Ltd., Beijing, China | 100 |
| Sartorius Stedim (Shanghai) Trading Co. Ltd., Shanghai, China | 100 |
| Sartorius Stedim India Pvt. Ltd., Bangalore, India | 100 |
| Sartorius Stedim Japan K.K., Tokyo, Japan | 100 |
| Sartorius Korea Biotech Co. Ltd., Seoul, South Korea | 69 |
| Sartorius Stedim Malaysia Sdn. Bhd., Kuala Lumpur, Malaysia | 100 |
| Sartorius Stedim Singapore Pte. Ltd., Singapore, Singapore | 100 |

There are no associates or joint ventures included in the scope of consolidation, all companies are consolidated in full.
The ownership rate equals the share in voting rights.

8. Business Combinations

None.

Notes to the Statement of Profit or Loss

9. Sales Revenue

Initial Application of IFRS 15, Revenue from Contracts with Customers

IFRS 15, Revenue from Contracts with Customers, defines a comprehensive model to determine when to recognize revenue and in which amount. It replaces existing guidelines for recognition of revenue, including IAS 18, Revenue; IAS 11, Construction Contracts; and IFRIC 13, Customer Loyalty Programmes.

The Group applied IFRS 15 for the first time as of January 1, 2018 on the basis of the modified retrospective method. Accordingly, the comparative period is presented in line with previous standards. Any effects of the initial application of IFRS 15 are recorded as an adjustment to the opening balance of retained earnings at the date of initial application. Furthermore, upon initial adoption, the Group has been applying IFRS 15 only to contracts that are not considered completed contracts at the date of initial application. Besides, the Group used the practical expedients offered regarding contract modifications that occurred prior to the date of initial application of IFRS 15 and did not retrospectively restate such contracts. As the impact of the new standard on the Group's consolidated financial statements is low, the use of this practical expedient is expected not to have a material impact.

There are no material effects as a result of the application of the new standard. However, as of the date of initial application, there are certain effects with regard to construction contracts in the company's project business. Sales revenue of approximately €5 million, which was recognized in the past, is not recognized over time according to IFRS 15, but rather at a point in time. This results from new criteria introduced by IFRS 15 for revenue recognition over time for contracts on the construction of customer-specific goods. The effects resulted in a balance sheet extension by about €2 million at the date of initial application. There were no effects to be recognized in equity. The effects of applying the modified retrospective method on the opening statement of financial position as of January 1, 2018 are shown in the following table.

| Effects from the application of IFRS 15 on the opening balances of the statement of financial position as of Jan. 1, 2018 | Carrying amount Dec. 31, 2017 € in mn | Adjustments on adoption of IFRS 15 € in mn | Carrying amount Jan. 1, 2018 € in mn ³⁾ |
|---|---|---|--|
| Inventories | 185.6 | 4.1 | 189.7 |
| Amounts due from customers for contract work ¹⁾ | 7.0 | -1.9 | 5.0 |
| Trade receivables | 204.6 | | 204.6 |
| Trade receivables | 211.6 | -1.9 | 209.7 |
| Trade payables payments received for orders ²⁾ | 40.7 | 2.2 | 42.9 |
| Trade payables | 77.2 | | 77.2 |
| Trade payables | 118.0 | 2.2 | 120.2 |

¹⁾ Contract assets according to IFRS 15.

²⁾ Contract liabilities according to IFRS 15.

³⁾ Carrying amount without adjustments for the new credit impairment model according to IFRS 9.

If the Group had applied the previous standards, IAS 18 and IAS 11, respectively, during the reporting period, sales revenue as well as cost of sales would have been higher by about €0.6 million while there would be no impact on the net result. Total assets and total liabilities would have been lower by about €2 million. The effects result from the revised criteria

for revenue recognition over time according to which revenue is now recognized at a point in time rather than over time for one project. The following table presents the impact on the consolidated statement of financial position as of December 31, 2018 that the continued application of the previous standards would have had in comparison with IFRS 15.

| Effects from the application of IFRS 15 on the consolidated statement of financial position as of December 31, 2018 | Carrying amount December 31, 2018 (IFRS 15) € in mn | Adjustments on adoption of IFRS 15 € in mn | Carrying amount December 31, 2018 (IAS 18, IAS 11) € in mn |
|---|---|---|--|
| Inventories | 252.0 | -5.1 | 246.9 |
| Amounts due from customers for contract work ¹⁾ | 3.9 | 2.9 | 6.8 |
| Trade receivables | 216.3 | | 216.3 |
| Trade receivables | 220.2 | 2.9 | 223.1 |
| Trade payables payments received for orders ²⁾ | 51.0 | -2.2 | 48.8 |
| Trade payables | 103.6 | | 103.6 |
| Trade payables | 154.6 | -2.2 | 152.4 |

¹⁾ Contract assets according to IFRS 15.

²⁾ Contract liabilities according to IFRS 15.

Other Disclosures

The revenues from contracts with customers according to IFRS 15 are disaggregated into geographical regions (see Segment Report, note 5). For details on the impairments on trade receivables and contract assets recognized in the reporting period see note 38.

The following table presents the opening balances as of January 1, 2018, and the balances as of December 31, 2018 for the Group's contract liabilities. The balances of trade receivables and contract assets are presented in note 28.

| | Line item in statement of financial position | Carrying amount as of January 1, 2018 € in k | Carrying amount as of December 31, 2018 € in k |
|--|---|--|--|
| Deferred revenue | Other liabilities | 10,153 | 16,235 |
| Payments received on account of orders | Trade payables | 42,949 | 51,000 |
| Contract liabilities (total) | | 53,101 | 67,234 |

There were no extraordinary changes in the carrying amounts of the contract liabilities and contract assets in the reporting period. Revenue in the amount of 41,182 k euros was recognized in the reporting period that was included in the contract liability balance at the beginning of the reporting period.

The Group produces and sells instruments and consumables for customers in the Biopharm segment.

The Group satisfies its performance obligations depending on the goods to be transferred and the promised services. The far majority of the revenues from sales of products is recognized at a point in time when the customer obtains control over the goods. This is typically the case when the significant risks and rewards of ownership of the goods are transferred to the customer. Therefore, the point in time may vary

depending on the agreement with the individual customer.

For complex products that require installation at the customer's site, revenue is recognized upon formal customer acceptance. To a low extent, revenue is recognized over time in the customer-specific project business. In these cases, revenue is recognized according to the project progress which is measured based on the percentage of costs to date compared to the total estimated contract costs. The amount of actual costs incurred to date reflects the progress and the transfer of control to the customer appropriately as the Group has a right to a reimbursement of cost to date plus an appropriate margin, if the project is cancelled by the customer without cause.

Revenue from services is generally recognized when the services are performed or have been performed. When the services are performed continuously over a period of time, the Group recognizes the related revenue over time. In this case, revenue is generally recognized pro rata temporis in relation to the total contract period. Product sales are typically accompanied by the legally required warranty. Any material extended warranties are accounted for as separate performance obligations.

According to the general payment terms, customer payments are due in the short-term, typically within 30 days. To some extent, the Group obtains advance payments, e.g. to avoid credit risks. Therefore, the Group regularly has contract liabilities (payments received on account of orders). In addition, the Group recognizes contract liabilities in connection with service contracts (deferred revenues) when customers pay in advance.

There are no material effects from contracts with significant financing components. The Group uses the practical expedient regarding the existence of a significant financing component. This means that a financing component is only taken into consideration when the length of time between the transfer of goods or services and the receipt of consideration is expected to exceed one year and the effect is material. As of December 31, 2018, the Group has refund liabilities in the amount of 5,755k euros as a result from incentive agreements with customers.

The aggregate amount of the transaction price allocated to the performance obligations that are unsatisfied (or partially unsatisfied) at the end of the reporting period (orders on hand) amounts to 489.0million euros. The Group expects that these unsatisfied performance obligations will be satisfied to a large extent in 2019.

10. Functional Costs

The statement of profit or loss has been presented according to the "cost of sales format", i.e. expenses have been allocated to the relevant functions production, sales & marketing, research & development and general administration.

Expenses relating to cross-functional initiatives or projects are assigned to the respective functional costs based on an appropriate allocation principle.

The caption "cost of sales" includes the costs of products sold and the acquisition costs of merchandise sold. In addition to directly attributable expenses, such as raw materials and supplies, employee benefits expense and energy expenses, cost of sales also includes overhead, which can be allocated to the manufacturing area, and the corresponding depreciation and amortization.

The selling and distribution costs pertain, in particular, to the costs of the sales and marketing function, distribution and market research.

Research and development costs comprise the costs of research and product and process development, unless they are recognized as assets.

The item "general administrative expenses" mainly includes employee benefits expense and the cost of materials of the general administrative area.

All profit and loss items that cannot be allocated to one of the mentioned functional areas are recognized as other income and expenses. This includes essentially effects from translation of transactions in foreign currencies, sale of fixed assets, allowances on trade receivables and reorganization and other non-recurring expenses. Income from grants related to income is recognized as other income, when there is reasonable assurance that the conditions attached to the grants are complied with and the grants will be received. They are recognized systematically as income over the period in which the related costs are recorded.

Operating expenses by nature are presented in the Profit or Loss Statement by nature in note 14.

The material expense and personnel cost are as follows:

Raw Materials and Supplies

This caption consists of the following:

| | 2018 12 months € in K | 2017 12 months € in K |
|----------------------------|-----------------------------|-----------------------------|
| Purchases consumed | 237,889 | 235,762 |
| Cost of purchased services | 54,079 | 48,353 |
| Total | 291,968 | 284,114 |

Personnel Cost

This caption can be broken down as follows:

| | 2018 12 months € in K | 2017 12 months € in K |
|---|-----------------------------|-----------------------------|
| Wages and salaries | 274,115 | 244,438 |
| Social security | 59,877 | 54,481 |
| Expenses for retirement benefits and pensions | 6,861 | 8,088 |
| Total | 340,853 | 307,007 |

11. Other Operating Income and Expenses

| | 2018 12 months € in K | 2017 12 months € in K |
|--|-----------------------------|-----------------------------|
| Currency translation gains | 10,392 | 7,652 |
| Extraordinary income | 35,248 | 0 |
| Income from the decrease in allowances for bad debts | 475 | 780 |
| Income from release of provisions and liabilities | 353 | 973 |
| Income from grants | 4,235 | 8,013 |
| Other income | 1,921 | 1,139 |
| Other operating income | 52,624 | 18,557 |
| Currency translation losses | -10,403 | -14,762 |
| Extraordinary expenses | -22,526 | -22,564 |
| Allowances for bad debts | -2,041 | -801 |
| Other expenses | -4,153 | -3,416 |
| Other operating expenses | -39,124 | -41,543 |
| Total other operating income and expenses | 13,500 | -22,987 |

The item reported as income from grants discloses the grants for expenses (essentially related to research and development projects).

Extraordinary items amounted to €12.7million (net) (previous year: €-22.6million). The extraordinary income in 2018 refers mainly to the derecognition of the liability for the acquisition of the cell culture media business as a consequence of the modified cooperation agreement with Lonza; extraordinary expenses essentially cover one-time expenses for strategic Group projects and integration and acquisition-related items.

12. Financial Result

| | 2018 12 months € in K | 2017 12 months € in K |
|---|-----------------------------|-----------------------------|
| Interest and similar income | 385 | 274 |
| - of which from affiliated companies | 126 | 18 |
| Income from derivative financial instruments | 2,798 | 801 |
| Other financial income | 2,167 | 8,425 |
| Financial income | 5,349 | 9,500 |
| Interest and similar expenses | -3,344 | -3,834 |
| - of which from affiliated companies | -1,378 | -1,814 |
| Expenses for derivative financial instruments | -5,779 | -1,706 |
| Interest expense for pensions | -649 | -585 |
| Other financial expenses | -11,263 | -4,465 |
| Financial expenses | -21,035 | -10,589 |
| Total | -15,685 | -1,089 |

The other financial income in 2017 (other financial expenses in 2018) includes mainly foreign exchange gains (losses) in connection with bank deposits, loans denominated in foreign currencies as well as the Lonza liability. Furthermore, in 2018 an increase in the liability for the phantom units in AllPure has been posted with an amount of approx. €3 mn (see note 31).

The interest expenses to affiliated companies are in connection with the loan granted by the Group's ultimate parent Sartorius AG (see also chapter 41).

13. Income Taxes

| | 2018 12 months € in K | 2017 12 months € in K |
|----------------------|-----------------------------|-----------------------------|
| Current income taxes | -78,330 | -56,173 |
| Deferred taxes | 3,740 | -676 |
| Total | -74,590 | -56,849 |

Current income taxes are determined based on the respective local taxable income of the period and local tax rules. In addition, current income taxes include adjustments for uncertain tax payments or tax refunds for periods not yet assessed. Changes in deferred tax assets and liabilities are included in income taxes except for changes recognized in other comprehensive income or equity.

Income taxes in France are calculated at 33.33% of the estimated taxable profit for the year. For Germany, a rate of approx. 30% was applied to the taxable income. Income generated outside France and Germany is taxed at rates applicable in the corresponding country.

Considering the French and German average tax rates and the impact of other tax legislations, the expected tax rate for the Sartorius Stedim Biotech Group is roughly 26% (29% in 2017). The reduction of 3% in the expected tax rate is mainly in connection with the reduced corporate tax rate in the US. The following table describes the difference between the expected tax expense and the income tax expenses reported for the particular financial year.

| | 2018 12 months € in K | 2017 12 months € in K |
|--|-----------------------------|-----------------------------|
| Expected tax rate | 26% | 29% |
| Expected tax expense | -73,983 | -63,978 |
| Differences from the Group average income tax rate | 9,771 | 8,638 |
| Permanent differences | -11,172 | -1,551 |
| Tax-free income and other tax exemptions | 1,222 | 1,262 |
| Taxes for previous years | 604 | -853 |
| Withholding and similar taxes | -404 | -596 |
| Other | -628 | 228 |
| Total | -74,590 | -56,849 |
| Effective tax rate | -26.2% | -25.8% |

14. Profit or Loss Statement by Nature

| | 2018 12 months € in K | 2017 12 months € in K |
|--|-----------------------------|-----------------------------|
| Sales revenue | 1,212,152 | 1,081,033 |
| Purchases consumed | -237,889 | -235,762 |
| Cost of purchased services | -54,079 | -48,353 |
| Personnel costs | -340,853 | -307,007 |
| Amortization and depreciation | -60,914 | -50,678 |
| Other operating costs | -218,183 | -217,531 |
| Subtotal | -911,918 | -859,331 |
| Operating profit (EBIT) | 300,234 | 221,702 |
| Financial income expenses | -15,685 | -1,089 |
| Income tax | -74,590 | -56,849 |
| Non-controlling interest | -1,907 | -2,676 |
| Net profit after non-controlling interest | 208,052 | 161,088 |

15. Earnings per Share

According to IAS 33, the earnings per share must be determined as follows: The basic earnings per share (basic EPS) are calculated on the basis of the weighted average number of ordinary shares during the period.

| | 2018 | 2017 |
|---|-------------------|-------------------|
| Net profit after tax (€ in K) | 209,959 | 163,763 |
| Group net profit after tax (€ in K) | 208,052 | 161,088 |
| Earnings per share (€) | 2.26 | 1.75 |
| Diluted earnings per share (€) | 2.26 | 1.75 |
| Number of shares (statutory level) | 92,180,190 | 92,180,190 |
| Treasury shares | -3,416 | 0 |
| Weighted average number of shares used in earnings per share calculation | 92,176,774 | 92,180,190 |
| Weighted average number of shares used in diluted earnings per share calculation | 92,176,774 | 92,180,190 |

Notes to the Individual Balance Sheet Items

16. Goodwill and Other Intangible Assets

Goodwill

| | Goodwill € in K |
|---|--------------------|
| Gross book values at Jan. 1, 2017 | 344,777 |
| Currency translation | -4,791 |
| Business combinations | 46,059 |
| Gross book values at Dec. 31, 2017 | 386,045 |
| Impairment losses at Jan. 1, 2017 | 0 |
| Currency translation | 0 |
| Impairment losses | 0 |
| Impairment losses at Dec. 31, 2017 | 0 |
| Net book values at Dec. 31, 2017 | 386,045 |
| | |
| | Goodwill € in K |
| Gross book values at Jan. 1, 2018 | 386,045 |
| Currency translation | -1,350 |
| Business combinations | 0 |
| Gross book values at Dec. 31, 2018 | 384,695 |
| Impairment losses at Jan. 1, 2018 | 0 |
| Currency translation | 0 |
| Impairment losses | 0 |
| Impairment losses at Dec. 31, 2018 | 0 |
| Net book values at Dec. 31, 2018 | 384,695 |

The caption reported as goodwill in the amount of €384,695K is the capitalized difference in assets resulting from business combinations. According to IAS 36, goodwill acquired in a business combination may not be amortized, but rather, must be tested annually for impairment and as soon as there is any indication of asset impairment. The increase recorded in 2017 concerns the acquisition of Umetrics.

For the purpose of impairment testing, goodwill must be allocated to each of the acquirer's cash-generating units (CGUs) that are expected to benefit from the synergies of the combination. The CGU represents the lowest level within the entity at which goodwill is monitored for internal management purposes and may not be larger than a segment. Sartorius Stedim Biotech

Group follows the strategy to be a total solution provider for its customers. Because of the various interdependencies within the business, the lowest level at which goodwill is monitored is that of the Biopharm segment. Therefore, the acquired goodwill is allocated to this group of CGU's.

As in 2017, the impairment test conducted for 2018 measures the recoverable amount on the basis of the value in use of the particular cash-generating unit (Biopharm segment). The cash flow forecasts consider previous experiences and are generally based on Group management's forecasts for a period of four years. The calculations were based on a terminal year growth rate of 2.5% for the years after 2022. This rate is derived from market expectations, which forecast significant growth rates for the targeted biopharmaceutical market. The major growth driver for the Sartorius Stedim Biotech Group will be the aging and increase in population and the improved access to drugs in the emerging markets as well as the ongoing paradigm shift from reusable products to single-use products utilized in bio manufacturing by the biopharmaceutical industry.

The discount rates applied correspond to the weighted capital cost rates and were recognized as follows:

| | 2018 | | 2017 | |
|------------------|---------------|-----------|------------|-----------|
| | Before tax | After tax | Before tax | After tax |
| Biopharm segment | 8.9% | 7.2% | 8.5% | 6.8% |

In 2018, our impairment test did not result in recognition of impairment losses. In this context, various sensitivity analyses based on realistic variations of the assumptions disclosed above did not result in an impairment either. The following variations would theoretically represent the "break-even point":

| | 2018 | 2017 |
|----------------------|--------|--------|
| Discount rates | 22.9% | 20.4% |
| Terminal growth rate | -29.6% | -28.4% |
| Cash flows | -78.5% | -80.2% |

Intangible Assets

| | Patents, licenses and similar rights € in K | Brand name € in K | Customer relationships € in K | Capitalized development costs € in K | Payments on account € in K | Total € in K |
|--|--|----------------------|-------------------------------------|---|----------------------------------|-----------------|
| Gross book values at Jan. 1, 2017 | 74,664 | 11,061 | 121,351 | 68,054 | 36 | 275,165 |
| Currency translation | -2,817 | -47 | -1,006 | -387 | 0 | -4,257 |
| Business combinations | 11,614 | 892 | 14,486 | 0 | 0 | 26,992 |
| Acquisitions | 5,767 | 0 | 469 | 20,788 | 0 | 27,024 |
| Disposals | -34 | 0 | 0 | -2,685 | 0 | -2,720 |
| Transfers | -143 | 0 | 0 | 155 | 0 | 12 |
| Gross book values at Dec. 31, 2017 | 89,051 | 11,906 | 135,300 | 85,924 | 36 | 322,216 |
| Amortization and impairment losses at Jan. 1, 2017 | -27,813 | -68 | -67,710 | -35,556 | 0 | -131,147 |
| Currency translation | 1,298 | 5 | 410 | 139 | 0 | 1,853 |
| Amortization and impairment losses | -7,306 | -98 | -10,072 | -5,176 | 0 | -22,652 |
| Disposals | 34 | 0 | 0 | 2,607 | 0 | 2,641 |
| Transfers | -5 | 0 | 0 | 0 | 0 | -5 |
| Amortization and impairment losses at Dec. 31, 2017 | -33,792 | -160 | -77,372 | -37,986 | 0 | -149,310 |
| Net book values at Dec. 31, 2017 | 55,259 | 11,746 | 57,928 | 47,937 | 36 | 172,906 |

| | Patents, licenses and similar rights € in K | Brand name € in K | Customer relationships € in K | Capitalized development costs € in K | Payments on account € in K | Total € in K |
|--|--|----------------------|-------------------------------------|---|----------------------------------|-----------------|
| Gross book values at Jan. 1, 2018 | 89,051 | 11,906 | 135,300 | 85,924 | 36 | 322,216 |
| Currency translation | 246 | -32 | 172 | -52 | 1 | 335 |
| Business combinations | 0 | 0 | 0 | 0 | 0 | 0 |
| Acquisitions | 20,443 | 0 | 0 | 22,839 | 19 | 43,301 |
| Disposals | -11,010 | 0 | -11,804 | -1,089 | 0 | -23,902 |
| Transfers | 363 | 0 | 0 | 0 | 0 | 363 |
| Gross book values at Dec. 31, 2018 | 99,094 | 11,874 | 123,669 | 107,622 | 55 | 342,313 |
| Amortization and impairment losses at Jan. 1, 2018 | -33,792 | -160 | -77,372 | -37,986 | 0 | -149,310 |
| Currency translation | -444 | 1 | -41 | 28 | 0 | -455 |
| Amortization and impairment losses | -11,519 | -152 | -13,991 | -4,341 | 0 | -30,004 |
| Disposals | 8,183 | 0 | 5,680 | 1,089 | 0 | 14,952 |
| Transfers | -3 | 0 | 0 | 0 | 0 | -3 |
| Amortization and impairment losses at Dec. 31, 2018 | -37,576 | -311 | -85,724 | -41,210 | 0 | -164,821 |
| Net book values at Dec. 31, 2018 | 61,518 | 11,563 | 37,944 | 66,412 | 55 | 177,492 |

Intangible assets acquired are recorded at cost less the accumulated, regular amortization that is calculated according to the straight-line method and any impairment loss. The useful life of an intangible asset is the period during which the Group expects to use the asset.

Amortization of intangible assets is based on the following estimated useful lives:

| | |
|-------------------------------------|-----------------------|
| Software | 2 to 5 years |
| Capitalized R&D expenses | 4 to 6 years |
| Customer relations and technologies | 5 to 15 years |
| Brand name | 5 years to indefinite |

Costs incurred within the scope of the development of new products and methods were capitalized as internally generated intangible assets if the following criteria were met:

- The technical feasibility of completing the intangible assets so that it will be available for use or sale;
- The intention to complete the intangible asset and use or sell it;
- The ability to use or sell the intangible asset;
- The demonstration of how the intangible asset will generate probable future economic benefits;
- The availability of adequate technical, financial and other resources to complete the development and to use or sell the intangible asset;
- The ability to measure reliably the expenditure attributable to the intangible asset during its development.

The capitalized development costs essentially cover the costs that were allocated to the staff involved in R&D, raw materials and supplies, outside services and directly attributable overhead. Intangible assets generated internally are amortized on a straight line basis over their useful lives, which generally do not exceed six years.

If an internally generated intangible asset may not be recognized, the development costs are expensed in the period in which they are incurred. Costs for research activities are reported as expenses in the period in which they are incurred.

The Stedim brand name acquired in 2007 is considered to have an indefinite useful life and is therefore not amortized. There is no foreseeable limit to the period over which the brand name is expected to generate net cash inflows for the Group. The brand name is tested annually for impairment at the level of the "Biopharm segment" cash-generating unit (CGU).

In 2018, the development costs of €22,839 K were recognized as assets (€20,788 K in 2017).

Amortization of intangible assets is allocated to the corresponding functions in the statement of profit or loss. For capitalized development costs, amortization is disclosed in the line "cost of sales".

As a consequence of the modified agreement with Lonza in the cell culture media business, the related intangible assets (technology and customer relationships) have been recognized as disposals (see note 11).

In 2018, impairments were recognized in the amount of €5.8 million. The impairments relate mainly to technologies and customer relationships.

17. Property, Plant and Equipment

| | Land, buildings and improvements € in K | Technical machinery and equipment € in K | Factory and office equipment and other equipment € in K | Payments on account and construction in progress € in K | Total € in K |
|---|--|---|--|--|-----------------|
| Gross book values at Jan. 1, 2017 | 171,308 | 152,063 | 81,784 | 36,704 | 441,860 |
| Currency translation | -4,127 | -2,165 | -2,873 | -4,742 | -13,907 |
| Business combinations | 0 | 0 | 141 | 0 | 141 |
| Acquisitions | 2,317 | 7,438 | 10,995 | 88,958 | 109,708 |
| Disposals | -454 | -2,683 | -3,153 | -1,797 | -8,087 |
| Transfers | 7,494 | 7,209 | 1,710 | -15,944 | 469 |
| Gross book values at Dec. 31, 2017 | 176,538 | 161,862 | 88,603 | 103,179 | 530,183 |
| Depreciation at Jan. 1, 2017 | -52,038 | -78,460 | -49,899 | 1 | -180,396 |
| Currency translation | 1,033 | 764 | 2,246 | 0 | 4,042 |
| Depreciation | -6,947 | -12,687 | -8,393 | 0 | -28,026 |
| Disposals | 436 | 2,553 | 2,978 | 0 | 5,967 |
| Transfers | -405 | 190 | -260 | 0 | -475 |
| Depreciation at Dec. 31, 2017 | -57,921 | -87,641 | -53,327 | 1 | -198,888 |
| Net book values at Dec. 31, 2017 | 118,617 | 74,222 | 35,275 | 103,181 | 331,295 |

| | Land, buildings and improvements € in K | Technical machinery and equipment € in K | Factory and office equipment and other equipment € in K | Payments on account and construction in progress € in K | Total € in K |
|---|--|---|--|--|-----------------|
| Gross book values at Jan. 1, 2018 | 176,538 | 161,862 | 88,603 | 103,179 | 530,183 |
| Currency translation | 386 | -302 | 445 | 3,709 | 4,238 |
| Business combinations | 0 | 0 | 0 | 0 | 0 |
| Acquisitions | 20,018 | 11,247 | 9,137 | 93,242 | 133,644 |
| Disposals | -3,643 | -7,316 | -9,959 | -299 | -21,218 |
| Transfers | 6,035 | 11,899 | 852 | -19,113 | -326 |
| Gross book values at Dec. 31, 2018 | 199,335 | 177,390 | 89,078 | 180,718 | 646,521 |
| Depreciation at Jan. 1, 2018 | -57,921 | -87,641 | -53,327 | 1 | -198,888 |
| Currency translation | 35 | -242 | -133 | 0 | -341 |
| Depreciation | -8,490 | -13,443 | -8,980 | 0 | -30,913 |
| Disposals | 3,320 | 6,712 | 9,257 | 0 | 19,290 |
| Transfers | -11 | 72 | 250 | 0 | 311 |
| Depreciation at Dec. 31, 2018 | -63,067 | -94,542 | -52,933 | 1 | -210,541 |
| Net book values at Dec. 31, 2018 | 136,268 | 82,848 | 36,146 | 180,719 | 435,980 |

The "Property, plant and equipment" caption is recorded at cost, and related assets are depreciated over their estimated useful life using the straight line method. Tangible assets are subject to impairment tests whenever there are indicators of impairment.

Depreciation of fixed assets is based on the following periods of useful life:

| | |
|------------------------------|----------------|
| Buildings | 15 to 50 years |
| Machinery | 5 to 15 years |
| Factory and office equipment | 3 to 13 years |

Depreciation is included in the statement of profit or loss according to use of the assets in the cost of sales, selling and distribution costs, research and development costs, general administrative expenses and other operating expenses.

Borrowing costs are expensed as incurred unless they are directly attributable to the acquisition, construction or production of a qualifying asset and are therefore part of the cost of that asset. A qualifying asset is defined as an asset that takes a substantial period of time (six to twelve months) to get ready for its intended use.

Grants related to assets are deducted from the cost of the related asset.

In 2018, as for fiscal 2017, there were no significant impairment losses to be recognized with regard to property, plant and equipment.

18. Leases

A lease is an arrangement whereby the lessor conveys to the lessee in return for a payment or series of payments the right to use an asset for an agreed period of time. According to IAS 17 a lease is classified as either an operating or a finance lease. A finance lease is a lease that transfers substantially all the risks and rewards incidental to ownership of an asset. All other leases are designated as operating leases.

When the Group is a lessee in a finance lease, the amount equal to the fair value of the leased property, or if lower, the present value of the minimum lease payments is recognized as an asset on the balance sheet and simultaneously recognized as a financial liability. The minimum lease payments essentially consist of the finance charge and the reduction of the outstanding liability, which are measured according to the effective interest method. A leased asset is depreciated on a straight-line basis over the period of its expected useful life or over the shorter lease term.

For an operating lease, the lease instalments to be paid by the lessee are recognized as expenses over the lease term and the lease payments received by the lessor are recognized as income, respectively. The leased asset continues to be recognized on the lessor's balance sheet as property, plant and equipment.

Capitalized property, plant and equipment for assets held under finance leases amount to €14,599 K (2017: €15,672). The cost of acquisition of these assets totals €19,631 K (2017: €19,682 K).

In 2018 rental payments amounting to €14.0 million (2017: €11.3 million) were made for assets leased under operating leases.

For the new lease accounting rules according to IFRS 16 (applicable from 2019 on) please refer to note 2.

The Group's financial obligations from operating lease commitments (future minimum lease payments under non-cancellable leases), are as follows:

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|---------------------------|-------------------------|-------------------------|
| Operate leases | | |
| - due within one year | 11,434 | 8,682 |
| - due within 2 to 5 years | 22,145 | 13,900 |
| - due thereafter | 8,813 | 4,817 |

19. Deferred Tax

| | Deferred Tax Assets | | Deferred Tax Liabilities | |
|---|-------------------------|-------------------------|--------------------------|-------------------------|
| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
| Intangible assets | 1,066 | 783 | 37,746 | 36,630 |
| Tangible assets | 0 | 0 | 5,865 | 6,553 |
| Inventory | 10,236 | 9,143 | 0 | 0 |
| Receivables and other current assets | 230 | 189 | 640 | 5,105 |
| Provisions | 6,384 | 6,649 | 0 | 0 |
| Liabilities | 5,391 | 2,181 | 1,866 | 3,140 |
| Gross amount | 23,307 | 18,945 | 46,117 | 51,428 |
| Carry forward of taxable losses | 0 | 2,334 | 0 | 0 |
| Tax on undistributed earnings of subsidiaries | 0 | 0 | 1,850 | 1,848 |
| Offset | -8,817 | -10,289 | -8,817 | -10,289 |
| Net amount | 14,490 | 10,989 | 39,150 | 42,987 |
| Change | 3,501 | 235 | 3,837 | -14,207 |
| thereof recognized in profit or loss | 3,450 | 292 | 291 | -968 |

Deferred tax assets or liabilities are determined based on temporary differences between the carrying amounts and the tax basis of assets and liabilities (except in special cases provided by IAS 12) including loss carry forwards and tax credits. Measurement is based on the tax rates expected to be effective in the period in which an asset is realized or a liability is settled.

For this purpose, the tax rates and tax rules are used which have been enacted or substantively enacted at the reporting date. Deferred tax assets are recognized for deductible temporary differences and tax losses and unused tax credits only to the extent that it is probable that the Group will have future taxable income against which they can be charged.

Deferred Tax Assets

On the balance sheet date, the Group had unused tax loss amounts carried forward of €11.6 million to be deducted from future taxable profits (€18.2 million in 2017). A deferred tax amount was reported on approx. €0.0 million of these losses (€9.8 million in 2017). Concerning the remaining losses to be carried forward, no deferred tax amounts were recognized because of the lack of visibility of future taxable profits.

Deferred tax assets in the amount of €0.0 million (€2.3 million) relate to companies that reported losses in this year under review or in the earlier reporting year.

Deferred Tax Liabilities

The deferred tax liabilities in connection with intangible assets refer to assets acquired in business combinations and consequently are mainly linked to customer relationships and technologies.

The Group did not record deferred tax liabilities on approx. €658 million (€480 million) in cumulative undistributed earnings of subsidiaries because these earnings are intended to be reinvested in these operations. When the dividends are paid out, an amount of 5% of the dividends will be taxed under the French and German taxation rules and, if applicable, with withholding tax. Furthermore, additional income tax consequences could arise in the case of an intermediate holding company.

In fiscal 2018, as in the previous years, the tax effect from hedging instruments, and the deferred tax assets from the recognition of the remeasurements of defined benefit liabilities (assets) were recognized in other comprehensive income. The income taxes recognized in other comprehensive income are disclosed as follows in the table:

| € in K | 2018 | 2017 |
|---|--------------|---------------|
| Cash flow hedges | 3,463 | -7,889 |
| Remeasurements of the net defined benefit obligations | -160 | 481 |
| Total | 3,303 | -7,408 |

20. Inventories

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|--------------------------------|-------------------------|-------------------------|
| Raw materials and supplies | 68,318 | 60,130 |
| Work in progress | 71,985 | 53,205 |
| Finished goods and merchandise | 108,520 | 68,008 |
| Payments on account | 3,179 | 4,261 |
| Total | 252,002 | 185,604 |

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|-------------------------------|-------------------------|-------------------------|
| Gross amount inventories | 269,013 | 199,824 |
| Write-downs | -17,011 | -14,221 |
| Net Amount Inventories | 252,002 | 185,604 |

Raw materials and supplies, including merchandise, are reported under "Inventories" at average cost. In principle, finished goods and work in progress are reported at cost of conversion. This cost includes direct costs, which can be allocated to these materials, and the appropriate portion of production and materials handling overhead, general administrative expenses and fixed assets at normal depreciation and/or amortization rates, based on the normal production capacity, provided that these expenses are caused by production.

Inventories must be valued at the lower amount of cost and the net realizable value. The net realizable value represents the estimated selling price in the ordinary course of business less the estimated costs of completion and the estimated costs necessary for marketing, sales and distribution. Where inventory risks exist, such as the risk of reduced shelf life as a result of storage periods or limited usability, inventories are marked down accordingly.

21. Issued Capital

Sartorius Stedim Biotech S.A. share capital consists of 92,180,190 shares with a par value of 0.20 €.

As of December 31, 2017, and December 31, 2018, there were no dilutive instruments. Shares registered

in the name of the same owner for at least four years benefit from a double voting right.

| | Dec. 31, 2018 | Dec. 31, 2017 |
|--|-------------------|-------------------|
| Number of shares at the beginning of the period | 92,180,190 | 92,180,190 |
| Number of shares at the end of the period | 92,180,190 | 92,180,190 |
| Nominal value per share (in €) | 0.20 | 0.20 |
| Issued capital amount (€ in K) | 18,436 | 18,436 |

Dividends

The Board of Directors will submit a proposal to the Annual General Shareholders' Meeting for payment of a dividend for the year ended December 31, 2018, as follows: payment of a net dividend of €0.57 per share (2017: €0.46), i.e., a total disbursement of 52,540,761.00€ (excluding treasury shares; 2017: 42,402,887.00€).

22. Non-Controlling Interest

The non-controlling interest recognized in the statement of financial position amounting to €8,476K relate to the subsidiary Sartorius Korea Biotech. The Group's interest in Sartorius Korea Biotech is 69%, the remaining 31% are subject to an exercisable call option.

The purchase price for this non-controlling interest is variable and depends on the future performance of the entity.

Key Figures

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|----------------------------------|-------------------------|-------------------------|
| Sartorius Korea Biotech Co. Ltd. | | |
| Sales revenue | 63,111 | 61,477 |
| Net result | 6,150 | 8,631 |
| Total assets | 36,225 | 34,399 |
| Attributed profit or loss | 1,907 | 2,676 |

There are no significant restrictions on the Group's ability to access or use the assets or settle the liabilities of the mentioned entity.

23. Pension and Employee Benefits Provisions

Pension Obligations

Pension provisions and similar obligations are recognized in the consolidated financial statements of Sartorius Stedim Biotech Group in accordance with actuarial principles. IAS 19, Employee Benefits, stipulates the Projected Unit Credit Method as the method of measurement. In addition to known pensions and life expectancies, this expected cash value method takes into account future salary and pension increases.

All remeasurements of the net defined benefit liability are recognized in other comprehensive income (pension reserves) in accordance with the standard IAS 19.

Defined Contribution Plans

Most of the Sartorius Stedim Biotech Group companies make payments under defined contributions plans, primarily related to government-run pension plans. In 2018, the total expense recognized for the defined contribution plans amounted to €20,807 K (2017: €19,527 K).

Defined Benefit Plans

The remeasurements of defined benefit liabilities (asset) are shown in other comprehensive income according to the standard IAS 19. The actuarial losses, which were transferred to the pension reserves, essentially resulted from a change in the discount rate and totaled €469 K (€-667 K in 2017).

An amount of €24,441 K relates in particular to pension provisions for retirement pension plans in Germany. These provisions totaled €25,068 K in 2017 and primarily relate to direct commitments under defined benefit pension plans. Under these commitments, the employees earn benefits for each year of service rendered to the company. The benefits earned depend on the salary level and the age of the respective employees. The pension benefits are generally not funded with assets.

The assumed discount rates reflect the interest rates payable on the reporting date for high-quality corporate bonds with matching maturities and denominated in the relevant currencies (mainly Euro). If such corporate bonds are not available with matching long-term maturities or are insufficiently

available, their matching interest rates are determined by extrapolation.

Measurement of the post-employment benefit obligations is based on the following actuarial assumptions:

For Germany:

| in % | Dec. 31, 2018 | Dec. 31, 2017 |
|--------------------------|---------------|---------------|
| Discount rate | 1.81 | 1.75 |
| Future salary increases | 3.00 | 3.00 |
| Future pension increases | 2.00 | 2.00 |

With regard to the assumptions for mortality and disability the tables "Richttafeln (RT) 2018 G" by Klaus Heubeck were applied.

For France:

| in % | Dec. 31, 2018 | Dec. 31, 2017 |
|--------------------------|---------------|---------------|
| Discount rate | 1.80 | 1.70 |
| Future salary increases | 2.00 | 2.50 |
| Future pension increases | 2.00 | 2.00 |

The amounts reported in the statement of profit or loss and other comprehensive income consist of the following:

| | 2018 € in K | 2017 € in K |
|---|----------------|----------------|
| Current service cost | -1,978 | -1,856 |
| Past service cost | 485 | 408 |
| Net interest expenses | -531 | -509 |
| Components of defined benefit costs recognized in profit or loss | -2,023 | -1,958 |
| Return on plan assets (excl. interest) | 56 | 57 |
| Remeasurements | 413 | -723 |
| Components of defined benefit costs recognized in other comprehensive income | 469 | -667 |
| Total | -1,554 | -2,624 |

In the statement of profit or loss, the current service cost is disclosed according to the assignment of employees to the respective functions.

The amount included in the consolidated statement of financial position arising from the Group's obligation in respect of defined benefit plans is as follows:

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|----------------------------------|-------------------------|-------------------------|
| Present value of the obligations | 46,459 | 43,752 |
| Fair value of the plan assets | -10,865 | -8,306 |
| Net Liability | 35,595 | 35,446 |

The present value of the defined benefit obligation developed as follows:

| | 2018 € in K | 2017 € in K |
|---|----------------|----------------|
| Present value of the obligations as of Jan. 1 | 43,752 | 40,841 |
| Current service cost | 1,978 | 1,856 |
| Past service cost | -485 | -408 |
| Interest cost | 649 | 585 |
| Remeasurements | -548 | 745 |
| Foreign currency translation differences | 386 | -725 |
| Retirement benefits paid in the reporting year | -1,011 | -852 |
| Employee contributions | 261 | 220 |
| Contributions by plan participants | 1,466 | 1,550 |
| Other changes | 14 | -60 |
| Present value of the obligations as of Dec. 31 | 46,459 | 43,752 |

The remeasurements of defined benefit liability (asset) can be allocated as follows:

| | 2018 € in K | 2017 € in K |
|------------------------------------|----------------|----------------|
| Experience adjustments | 367 | 850 |
| Changes in demographic assumptions | 306 | 632 |
| Changes in financial assumptions | -1,221 | -737 |
| Total | -548 | 745 |

Plan Assets

| | 2018 € in K | 2017 € in K |
|--|----------------|----------------|
| Plan assets as of Jan. 1 | 8,306 | 6,622 |
| Interest income | 118 | 75 |
| Return on plan assets (excl. interest) | 56 | 57 |
| Remeasurements | -136 | 22 |
| Group contribution & payments | -887 | -709 |
| Foreign currency translation differences | 279 | -481 |
| Employee contributions | 261 | 220 |
| Employer contributions | 1,402 | 894 |
| Contributions by plan participants | 1,466 | 1,606 |
| Other changes | 0 | 0 |
| Plan assets as of Dec. 31 | 10,865 | 8,306 |

Composition of Plan Assets

The plan assets do primarily refer to insurance contracts in Germany and Switzerland, there are no major equity or debt investments included. The subsidiary in South Korea has deposited an amount of €2.4 million (€1.8 million in 2017) to local banks as cash and cash equivalents.

Sensitivity Analysis

An increase|decrease of the actuarial assumptions would have the following impacts on the defined benefit obligations (a positive sign (+) means an increase of the obligation):

2017:

| € in K | | |
|--------------------------------|----------|----------|
| Demographic assumptions | | |
| Life expectancy | +1 year | -1 year |
| Effect | 874 | -870 |
| Financial assumptions | | |
| Discount rate | +100 bps | -100 bps |
| Effect | -6,105 | 7,848 |
| Future salary increases | +50 bps | -50 bps |
| Effect | 865 | -803 |
| Future pension increases | +25 bps | -25 bps |
| Effect | 984 | -937 |

2018:

| € in K | | |
|--------------------------------|----------|----------|
| Demographic assumptions | | |
| Life expectancy | +1 year | -1 year |
| Effect | 1,053 | -1,052 |
| Financial assumptions | | |
| Discount rate | +100 bps | -100 bps |
| Effect | -6,389 | 8,171 |
| Future salary increases | +50 bps | -50 bps |
| Effect | 938 | -861 |
| Future pension increases | +25 bps | -25 bps |
| Effect | 1,020 | -972 |

The sensitivity analysis presented above might not be representative of the actual change in the defined benefit obligation as it is unlikely that the change in assumptions would occur in isolation of one another. Furthermore, the present value of the defined benefit obligation has been calculated using the same method that was applied in calculating the defined benefit obligation liability recognized in the statement of financial position (projected unit credit method).

Maturity Analysis

The undiscounted cash flows from defined benefits obligations can be allocated to maturities as follows:

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|--------------|-------------------------|-------------------------|
| <1 year | 1,664 | 1,477 |
| 1-5 years | 8,336 | 7,616 |
| 6-10 years | 12,430 | 11,564 |
| >10 years | 73,605 | 58,288 |
| Total | 96,035 | 78,944 |

The weighted average duration of the defined benefit obligations is 16.6 years (2017: 17.6 years).

24. Other Provisions

A provision is recognized when a present legal or constructive obligation to third parties arising from past events has been incurred, an outflow of resources is probable and the amount of the obligation can be reasonably estimated. The amount recognized as a provision represents the best estimate of the obligation at the closing date.

Restructuring provisions are set up in connection with programs that materially change the scope of business performed by a segment or business unit or the manner in which business is conducted. In most cases, restructuring expenses include termination benefits and compensation payments due to the termination of agreements with suppliers and dealers, including leasing contracts. Restructuring provisions are recognized when the Group has a detailed formal plan that has either commenced implementation or been announced.

Other Non-current Provisions

| | Payments to employees on early retirement plan € in K | Other € in K | Total € in K |
|---------------------------------|---|-----------------|-----------------|
| Balance at Jan. 1, 2017 | 2,331 | 752 | 3,083 |
| Currency translation | 0 | -11 | -11 |
| Consumption | -1,035 | -16 | -1,051 |
| Additions | 877 | 205 | 1,082 |
| Balance at Dec. 31, 2017 | 2,173 | 930 | 3,103 |

| | Payments to employees on early retirement plan € in K | Other € in K | Total € in K |
|---------------------------------|---|-----------------|-----------------|
| Balance at Jan. 1, 2018 | 2,173 | 930 | 3,103 |
| Currency translation | 0 | 9 | 9 |
| Consumption | -1,124 | -41 | -1,165 |
| Additions | 824 | 105 | 929 |
| Balance at Dec. 31, 2018 | 1,873 | 1,004 | 2,877 |

The non-current provisions comprise mainly provisions for partial retirement and employee anniversary bonuses (included in the item "other"). These obligations arise mainly in German Group companies. The partial retirement plans allow employees to work part-time for 3 - 5 years before their actual retirement.

According to IAS 19 these obligations are treated as severance payments to be earned in future periods and

are therefore recognized in profit or loss over the respective period of service. Actuarial gains and losses, as well as past service costs, on these obligations are recognized as income or expense.

Non-current provisions are reported at their present value on the reporting date. The discount rate for employees on the early retirement plan is 0.1% (2017: 0.0%).

Current Provisions

During financial 2017 and 2018, the current provisions developed as follows:

| | Warranties € in K | Other € in K | Total € in K |
|---------------------------------|----------------------|-----------------|-----------------|
| Balance at Jan. 1, 2017 | 4,914 | 4,367 | 9,281 |
| Currency translation | -56 | -73 | -129 |
| Consumption | -1,362 | -1,095 | -2,457 |
| Release | -837 | -1,539 | -2,377 |
| Additions | 2,165 | 3,651 | 5,816 |
| Other changes | 0 | -577 | -577 |
| Balance at Dec. 31, 2017 | 4,824 | 4,735 | 9,558 |

| | Warranties € in K | Other € in K | Total € in K |
|---------------------------------|----------------------|-----------------|-----------------|
| Balance at Jan. 1, 2018 | 4,824 | 4,735 | 9,558 |
| Currency translation | 4 | -14 | -9 |
| Consumption | -934 | -1,045 | -1,980 |
| Release | -2,550 | -649 | -3,199 |
| Additions | 5,021 | 2,892 | 7,913 |
| Other changes | 0 | 0 | 0 |
| Balance at Dec. 31, 2018 | 6,364 | 5,919 | 12,283 |

Warranty provisions contain expenses for replacement deliveries and repairs. Specific risks are recognized when the occurrence is more likely than not. General warranty risks are considered on the basis of experiences in the past. The other provisions contain mainly onerous contracts and uncertain liabilities to employees.

25. Other Financial Obligations | Contingent Assets and Liabilities

As was the case in the previous years there are no significant contingent liabilities or contingent assets to be reported.

Financial Instruments | Financial Risks

A financial instrument is any contract that gives rise to a financial asset of one entity and a financial liability or equity instrument of another entity. The following sections give an overview of the impact of financial instruments on the financial statements of the Sartorius Stedim Biotech Group and provide additional information on the balance sheet items, which contain financial instruments.

26. Initial Application of IFRS 9, Financial Instruments: Transition Effects and Accounting Policies

The initial date of application of IFRS 9, Financial Instruments is January 1, 2018. The Group applies IFRS 9 retrospectively without restating the prior-year figures. The comparative period is presented in line with previous rules (IAS 39). The aggregate effects from the transition to IFRS 9 are accounted for by way of an adjustment to the opening balance of retained earnings. The effect from the initial application of IFRS 9 results from the adjustment of the credit impairment approach for trade receivables and amounts to €0.3 million. As a result, this increased retained earnings from €612.7 million as of December 31, 2017 to €613.0 million as of January 1, 2018.

IFRS 9 replaces the existing guidelines in IAS 39, Financial Instruments: recognition and measurement. IFRS 9 includes revised guidelines for classification and measurement of financial instruments, including a new model of expected credit losses for the calculation of impairments of financial assets, as well as the new general guidelines for hedge accounting. This standard also adopts the guidance of IAS 39 for recognition and derecognition of financial instruments.

Under IFRS 9, the new classification and measurement approach for financial assets reflects both the entity's business model (held-to-collect, held-to-collect-and-sell, other) within the scope of which assets are held and the contractual cash flow characteristics (so-called "SPPI" criterion, solely payments of principal and interest). For each class of the Group's financial assets, the following table provides an overview of the categories according to IAS 39 and IFRS 9, as well as the respective carrying amounts at the date of initial application of IFRS 9.

| Carrying amounts of financial assets at the date of initial application of IFRS 9 | Category acc. to IAS 39 | Carrying amount acc. to IAS 39 Dec. 31, 2017 € in mn | Reclassification € in mn | Remeasure-ment € in mn | Carrying amount acc. to IFRS 9 Jan. 1, 2018 € in mn | Category acc. to IFRS 9 |
|---|------------------------------|---|-----------------------------|---------------------------|--|---|
| Cash and cash equivalents | Loans and receivables | 32.6 | | | 32.6 | Measured at amortized cost |
| Receivables and other assets | Loans and receivables | 12.8 | | | 12.8 | Measured at amortized cost |
| Derivative financial instruments in hedge relationships | n/a | 7.2 | | | 7.2 | n/a |
| Other financial assets (current) | | 20.0 | | | 20.0 | |
| Trade receivables | Loans and receivables | 204.6 | | 0.3 | 205.0 | Measured at amortized cost |
| Amounts due from customers for contract work | n/a | 7.0 | | 0.0 | 7.0 | n/a |
| Trade receivables | | 211.6 | | 0.3 | 211.9 | |
| Financial assets | Loans and receivables | 2.7 | 0.7 | | 3.4 | Measured at amortized cost |
| Financial assets | Available for sale | 1.6 | -1.0 | | 0.6 | Debt instruments at fair value through profit or loss |
| Financial assets | Available for sale | 0.1 | | | 0.1 | Equity instruments at fair value through profit or loss |
| Financial assets (investments in non-consolidated subsidiaries) | n/a | 0.1 | | | 0.1 | n/a (investments in non-consolidated subsidiaries) |
| Derivative financial instruments in hedge relationships | n/a | 7.3 | | | 7.3 | n/a |
| Financial assets (non-current) | | 11.8 | -0.4 | | 11.5 | |

There were no effects on the Group's financial liabilities. The few reclassifications of financial assets were determined on the basis of individual assessments of the financial instruments, especially with regard to the contractual cash flow characteristics. For equity instruments that existed as of the date of initial application of the Standard and that were not held for trading, the Group decided to recognize future changes in the fair value of those instruments in profit or loss. This choice is generally to be made on an instrument-by-instrument basis upon initial recognition of the instrument. Reclassification at the date of initial application did not lead to measurement effects. The measurement effects presented result from the adjustment of the impairment approach for trade receivables. There were no reclassifications of financial instruments during the reporting period.

With regard to the impairment of financial assets, IFRS 9 replaces the incurred-loss model by the expected-loss model. Financial assets are generally regarded as credit-impaired when there are objective

indications that cast doubt about the full collection of the cash flows of the respective financial assets. With regard to the financial assets of the Group, the simplified approach which is applied to trade receivables is especially of relevance.

At the date of initial application of IFRS 9, the allowance for expected credit losses amounted to €0.1 million. This amount includes the expected credit loss in relation to contract assets according to IFRS 15 that do not contain a significant financing component. In the course of the transition to IFRS 9, the Group also analyzed and revised its approach for the recognition of incurred losses, given its low level of historical losses. At the date of initial application of IFRS 9, the total effect from the adjustments to the impairment approach for trade receivables amounted to €0.3 million and increased the Group's retained earnings as of January 1, 2018.

The following table illustrates the effects from the adjusted impairment model for trade receivables:

| Valuation allowances at the date of initial application of IFRS 9 | Category acc. to IAS 39 | Loss allowances acc. to IAS 39 Dec. 31, 2017 € in mn | Remeasure-ment due to implementa-tion of IFRS 9 € in mn | Loss allowances acc. to IFRS 9 Jan. 1, 2018 € in mn | Category acc. to IFRS 9 |
|---|-------------------------|---|--|--|----------------------------|
| Trade receivables | Loans and receivables | -2.4 | 0.3 | -2.1 | Measured at amortized cost |
| Amounts due from customers for contract work | n/a | 0.0 | 0.0 | 0.0 | n/a |
| Trade receivables | | -2.4 | 0.3 | -2.1 | |

Besides trade receivables, cash and cash equivalents are the most material financial assets on the Group's statement of financial position as of the date of initial application of IFRS 9 and as of the reporting date, December 31, 2018. No impairment is recognized for these financial assets due to materiality considerations.

For the remaining financial assets that are measured at amortized cost, no impairment is recognized as of the date of initial application of IFRS 9 and as of December 31, 2018, for the 12-month expected credit losses, given the Group's immaterial historical losses.

Derivatives are measured at fair value determined according to the mark-to-market method in which recognized mathematical methods are used. The fair values are based on the market data available at the time the value of these derivatives is calculated and reflect the estimates of the market conditions at the end of the year. Those instruments that are not designated as hedging instruments and for which no hedge accounting is applied, are categorized as held for trading. Changes of the fair values of derivative financial instruments are either recognized in profit or loss or, in case of hedging relationships, in other comprehensive income.

The Group applies the new hedge accounting rules of IFRS 9 prospectively. The Group uses forward transactions to hedge cash flow risks that result from changes in foreign exchange rates in relation to sales of products and the procurement of materials, and designates only the spot element of the hedging instrument. No transition effects resulted from the application of the new hedge accounting requirements.

Financial Assets

Financial assets of the Group mainly include cash and cash equivalents, trade and loan receivables and derivative financial instruments with a positive fair value.

27. Cash and Cash Equivalents

The Group considers all highly liquid investments with less than three months maturity from the date of acquisition to be cash equivalents. This mainly includes checks, cash on hand and deposits in banks. Cash and cash equivalents are measured at cost. For purposes of the consolidated cash flow statement, cash and cash equivalents include cash and cash equivalents as defined above. As of December 31, 2018, cash and cash equivalents amounted to €23,975 K (2017: €32,552 K).

28. Current Trade Receivables | Other Receivables

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|---|-------------------------|-------------------------|
| Trade receivables from third parties | 204,931 | 192,866 |
| Amounts due from customers for contract work* | 3,902 | 6,967 |
| Receivables from subsidiaries of the Sartorius AG Group | 11,399 | 11,782 |
| Trade receivables | 220,231 | 211,616 |

*Contract assets according to IFRS 15.

The book values of trade receivables and other receivables are representative of their fair value considering the maturity date and the credit risks. The contract assets are recognized in connection with customer-specific construction contracts that meet the requirements for revenue recognition over time according to IFRS 15 (see section 9). The "Receivables from subsidiaries of the Sartorius AG Group" item refers to other companies of the Sartorius Group (please refer to section 41). Impairment losses on trade and other receivables are recognized using separate allowance accounts. For details on the determination of the impairment allowances see section 38.

30. Loans and Borrowings

| | Balance at Dec. 31, 2018 € in K | of which current Dec. 31, 2018 € in K | Balance at Dec. 31, 2017 € in K | of which current Dec. 31, 2017 € in K |
|--|---------------------------------------|--|---------------------------------------|---|
| Liabilities to banks | 38,278 | 35,153 | 29,367 | 23,117 |
| Loans from Sartorius AG | 94,501 | 54,501 | 109,481 | 69,481 |
| Other loans from Sartorius Group companies | 163 | 163 | 3,335 | 3,335 |
| Total loans and borrowings | 132,942 | 89,817 | 142,183 | 95,933 |

Sartorius Stedim Biotech Group has signed a loan agreement with its parent company Sartorius AG which secures the financing of the Sartorius Stedim Biotech Group over the long term. The credit volume of this agreement is approx. 300 million euros and the interest rate is variable with a credit margin based on arms'-length principles.

29. Other Financial Assets

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|----------------------------------|-------------------------|-------------------------|
| Derivative financial instruments | 2,824 | 7,207 |
| Other financial assets | 19,212 | 12,842 |
| Current financial assets | 22,036 | 20,049 |

The amount shown as derivative financial instruments represents the fair value of foreign currency hedging instruments, mainly forward contracts (for details refer to section 35).

Other financial assets are measured at amortized cost using the effective interest method less any impairment losses. The item "Other financial assets" includes loan receivables to other entities of the Sartorius AG Group in the amount of €2,588 K (2017: €1,542 K).

Financial Liabilities

Financial liabilities of the Group mainly comprise loans borrowed from Sartorius AG, trade payables, finance lease liabilities and derivative financial instruments with a negative fair value. The financial liabilities other than derivative financial instruments are measured at amortized cost.

The non-current loans and borrowings do not include the liabilities in connection with acquisitions which are presented in the caption "other non-current liabilities".

31. Other Non-current Liabilities

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|-------------------|--------------------------------|-------------------------|
| Other liabilities | 8,889 | 30,392 |
| Total | 8,889 | 30,392 |

The other non-current liabilities include the liability for phantom units that was incurred in connection with the acquisition of the non-controlling interests in the company AllPure Technologies, LLC amounting to €8,739 K. The purchase price depends on the future sales and is due 2022 at the latest. As a result of the positive development of the AllPure business, the expected payments are determined by considering future revenue at an annual growth rate of about 20% on average. Due to this change in expectations regarding revenue growth, the liability has been increased by about €3 million. The effect is recognized in profit or loss. An increase (decrease) of the sales revenue by 10% in each of the following years would lead to an increase (decrease) of the liability by €0.9 million (€0.9 million).

In the previous reporting period, the other non-current liabilities included the non-current part of the liability for the remaining purchase price for the cell culture media business of the company Lonza. In connection with the change in cooperation between the Group and Lonza, the Group was released from the obligation to pay the remaining purchase price in the reporting period. Therefore, the related liability was released in 2018 and the effects on profit or loss are recognized in other operating income (see note 11).

Trade Payables

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|--|--------------------------------|-------------------------|
| Payments received on account of orders* | 51,000 | 40,749 |
| Trade payables to third parties | 86,187 | 69,900 |
| Payables to participations | 6 | 2 |
| Payables to subsidiaries of the Sartorius AG Group | 17,375 | 7,309 |
| Total | 154,568 | 117,959 |

*Contract liabilities according to IFRS 15.

32. Other Current Financial Liabilities

| | Dec. 31, 2018 € in K | Dec. 31, 2017 € in K |
|----------------------------------|--------------------------------|-------------------------|
| Derivative financial instruments | 1,346 | 1,928 |
| Other liabilities | 12,966 | 33,036 |
| Total | 14,312 | 34,964 |

The derivative financial instruments refer to the fair values of foreign currency hedging transactions such as forward contracts (mainly related to the US\$).

The reduction in other financial liabilities is linked to the release of the liability for the acquisition of the cell culture media business (see note 11).

33. Carrying Amounts and Fair Values of Financial Instruments according to Categories

The following tables show the carrying amounts and fair values of financial assets and liabilities by category of financial instrument as of December 31, 2018, under IFRS 9 and as of December 31, 2017, under IAS 39:

| | Category acc. to IFRS 9 | Carrying amount Dec. 31, 2018 € in K | Fair value Dec. 31, 2018 € in K |
|---|---|--|---------------------------------------|
| Investments in non-consolidated subsidiaries | n/a | 109 | 109 |
| Financial assets | Equity instruments at fair value through profit or loss | 50 | 50 |
| Financial assets | Debt instruments at fair value through profit or loss | 671 | 671 |
| Financial assets | Measured at amortized cost | 4,814 | 4,814 |
| Financial assets (non-current) | | 5,644 | 5,644 |
| Amounts due from customers for contract work | n/a | 3,902 | 3,902 |
| Trade receivables | Measured at amortized cost | 216,330 | 216,330 |
| Trade receivables | | 220,231 | 220,231 |
| Receivables and other assets | Measured at amortized cost | 19,212 | 19,212 |
| Derivative financial instruments designated as hedging instruments* | n/a | 2,824 | 2,824 |
| Other financial assets (current) | | 22,036 | 22,036 |
| Cash and cash equivalents | Measured at amortized cost | 23,975 | 23,975 |
| Loans and borrowings | Financial liabilities at cost | 132,942 | 133,175 |
| Trade payables | Financial liabilities at cost | 103,568 | 103,568 |
| Trade payables payments received for orders | n/a | 51,000 | 51,000 |
| Trade payables | | 154,568 | 154,568 |
| Derivative financial instruments | Held for trading | 1,346 | 1,346 |
| Other financial liabilities | Financial liabilities at cost | 21,855 | 21,326 |
| Other financial liabilities | | 23,201 | 22,672 |

*The amounts include the non-designated part of the contracts.

| | Category acc. to IAS 39 | Carrying amount Dec. 31, 2017 € in K | Fair value Dec. 31, 2017 € in K |
|--|-------------------------------------|---|---------------------------------------|
| Financial Assets | Available for sale | 1,813 | 1,813 |
| Financial Assets | Loans and receivables | 2,734 | 2,734 |
| Derivative financial instruments in hedging relationship | Hedging Instruments* | 7,260 | 7,260 |
| Non-current financial assets | | 11,808 | 11,808 |
| Trade receivables | Loans and receivables | 211,616 | 211,616 |
| Financial Assets | Loans and receivables | 12,842 | 12,842 |
| Derivative financial instruments in hedging relationship | Hedging Instruments* | 7,207 | 7,207 |
| Other financial assets | | 20,049 | 20,049 |
| Cash and cash equivalents | Loans and receivables | 32,552 | 32,552 |
| Loans and borrowings | Financial liabilities at cost | 142,183 | 142,671 |
| Finance lease liabilities | IAS 17 | 17,487 | 28,086 |
| Trade payables | Financial liabilities at cost | 77,211 | 77,211 |
| Trade payables | n/a | 40,749 | 40,749 |
| Trade payables | | 117,959 | 117,959 |
| Derivative financial instruments | Held for trading | 552 | 552 |
| Derivative financial instruments in hedging relationship | Hedging Instruments* | 1,376 | 1,376 |
| Other financial liabilities | Financial liabilities at cost | 63,428 | 67,719 |
| Other financial liabilities | | 65,356 | 69,647 |

*The amounts include the non-designated part of the contracts.

The fair values of the financial instruments were determined on the basis of the market information available on the reporting date and are to be allocated to one of the three levels of the fair value hierarchy in accordance with IFRS 13.

Level 1 financial instruments are measured on the basis of prices quoted on active markets for identical assets and liabilities. In Level 2, financial instruments are measured on the basis of input factors which are derivable from observable market data or on the basis of market prices for similar instruments. Level 3 financial instruments are measured on the basis of input factors that cannot be derived from observable market data.

The financial instruments to be recognized at fair value on the reporting date are mainly derivatives in the form of forward contracts. They were measured on the basis of their quoted exchange rates and market yield curves (Level 2).

The fair values to be disclosed for financial liabilities recognized at amortized cost, especially liabilities to banks and finance leases, were measured on the basis of the market interest rate, taking the current indicative credit spreads into account (Level 2).

The fair values of the remaining financial assets and liabilities to be disclosed approximate the carrying amounts on account of their predominantly short-term maturity. The maximum credit loss risk is reflected by the carrying amounts of the financial assets recognized in the Statement of Financial Position.

The Group recognizes transfers between the levels of the fair value hierarchies at the end of the reporting period during which a change has occurred. In the current reporting period, there were no transfers between the levels.

34. Net Gains and Losses from Financial Instruments

The net gains and losses of the various categories of financial instruments are presented in the following table:

| Categories according to IFRS 9 | 2018 12 months € in K |
|--|-----------------------------|
| Financial assets measured at amortized cost | 2,771 |
| Financial assets and liabilities measured at fair value through profit or loss | 552 |
| Financial liabilities measured at amortized cost | -12,733 |

| Categories according to IAS 39 | 2017 12 months € in K |
|---|-----------------------------|
| Loans and receivables | -4,121 |
| Financial assets and liabilities held for trading | 2,358 |
| Financial liabilities at cost | 1,606 |

The net result from financial assets measured at amortized cost mainly includes the effects of currency translation and changes in allowances.

The net result from financial assets and liabilities held for trading predominantly comprises changes in the fair value of derivative financial instruments as well as interest income and interest expenses for these financial instruments.

The net result from liabilities measured at amortized cost mainly comprises the effects of foreign currency translation.

Total interest income and expenses for financial assets and liabilities that are not measured at fair value through profit or loss were as follows:

| | 2018 12 months € in K | 2017 12 months € in K |
|-------------------|-----------------------------|-----------------------------|
| Interest income | 463 | 282 |
| Interest expenses | -4,563 | -5,454 |

Capital and Financial Risk Management

Capital Risk Management

In the Sartorius Stedim Biotech Group, capital is managed in order to maximize earnings of those participating in the company by optimizing the ratio of equity to liabilities. Furthermore, we ensure that all Group companies operate under the premise of the going-concern principle.

The financial liabilities detailed above are regarded as managed capital and, furthermore, so are the cash and cash equivalents as well as equity capital.

Goals of Financial Risk Management

The Treasury Department of the Sartorius Stedim Biotech Group is centrally focused in Sartorius Corporate Administration GmbH, a subsidiary of Sartorius AG. This centralized Treasury Department performs services for all companies of the Sartorius Group, including the Sartorius Stedim Biotech Group, and coordinates access to national and international financial markets. In addition, the Treasury Department monitors and controls financial risks by internal risk reporting, which analyzes risks according to their degree and scope. Essentially, these risks entail currency, interest rate and liquidity risks as well as credit risks.

The Sartorius Stedim Biotech Group strives to minimize the impact of currency and interest rate risks using appropriate derivative financial instruments. Hedging transactions and their controlling are carried out by different staff members. Moreover, the Group's Internal Auditing Department regularly monitors the use of such financial instruments. Trading with derivative financial instruments is done for hedging purposes only.

35. Management of Exchange Rate Risks and Hedge Accounting

The Group is exposed to currency risks as more than one third of sales revenue is generated in U.S. dollars or currencies linked to the U.S. dollar and, to a lesser extent, in other foreign currencies. At the same time, the Group is able to compensate the major part of the revenues denominated in foreign currencies with costs incurred in the same currencies due its global production network. The share of revenues generated in foreign currencies that exceeds such costs, the so-called net currency exposure, is hedged with derivative

financial instruments to a large extent (generally 70% to 80%). The Group generally follows a rolling hedging strategy of up to 12 months in advance. Also, the hedging measures are reviewed at regular intervals in order to adapt them to currency fluctuations.

For currency hedging, forward contracts are used. Forward contracts secure the right, and simultaneously create the obligation, to sell an established foreign currency amount on the exercise date at a specific exchange rate against the euro, independently of the exchange rate actually valid on this date. The profit or loss resulting from the difference between the current and the previously established exchange rate is

recognized as income or expense in the statement of profit or loss.

At the balance sheet date, forward contracts have been carried out in an amount of \$137 million (2017: \$190 million) to hedge against the risk of fluctuation in the EUR|USD exchange rate. This amount covers roughly half of the expected net exposure for the U.S. dollar within the period of 12 months. Furthermore, other foreign currencies have been hedged in smaller volumes.

The following table shows the forward transactions as of the balance sheet date:

| Dec. 31, 2017 | Currency | Volume | Maturity | Fair value € in K |
|------------------|------------|----------------|----------|-------------------|
| Forward contract | USD | 110,000 | 2018 | 6,988 |
| | USD | 80,000 | 2019 | 7,260 |
| | USD | 190,000 | | 14,248 |
| Forward contract | JPY | 700,000 | 2018 | 72 |
| | JPY | 700,000 | | 72 |
| Forward contract | CHF | 24,500 | 2018 | -1,375 |
| | CHF | 24,500 | | -1,375 |
| Forward contract | CAD | 8,800 | 2018 | 148 |
| | CAD | 8,800 | | 148 |

| Dec. 31, 2018 | Currency | Volume | Maturity | Fair value € in K |
|------------------|------------|------------------|----------|-------------------|
| Forward contract | USD | 50,000 | 2019 | -53 |
| | USD | 50,000 | 2019 | 296 |
| | USD | 21,000 | 2019 | 712 |
| | USD | 16,000 | 2019 | 644 |
| | USD | 137,000 | | 1,599 |
| Forward contract | JPY | 400,000 | 2019 | -138 |
| | JPY | 200,000 | 2019 | -67 |
| | JPY | 200,000 | 2019 | -51 |
| | JPY | 250,000 | 2019 | -52 |
| | JPY | 1,050,000 | | -308 |
| Forward contract | CHF | 3,000 | 2019 | 4 |
| | CHF | 3,000 | | 4 |
| Forward contract | GBP | 3,500 | 2019 | 18 |
| | GBP | 3,500 | | 18 |
| Forward contract | CAD | 2,000 | 2019 | 32 |
| | CAD | 4,000 | 2019 | 90 |
| | CAD | 2,000 | 2019 | 43 |
| | CAD | 8,000 | | 165 |

Derivative financial instruments are measured at the time of acquisition at cost and at fair value on subsequent balance sheet dates. The changes in value of the derivative financial instruments are generally recognized in the statement of profit or loss on the reporting date. If the derivative financial instruments serve to hedge against cash flow risk arising from exchange rate risks and a qualified hedging relationship exists based on the criteria of IFRS 9, the valuation adjustments of the effective portion of the instrument are recognized in other comprehensive income (cumulative amount in 2018: €4.8 million; 2017: €16.4 million). Only the spot element of the forward contracts used to hedge the cash flow risks is designated as hedging instrument. The amounts recognized in equity are included in the profit or loss in the period in which the hedged transactions affect profit or loss. The changes of the hedging reserves are explained below as well as in the statement of changes

in equity. The non-designated or ineffective portion of the hedging instruments is recognized within the financial result in profit or loss.

The economic relationship between hedging instrument and hedged item and the effectiveness of the hedge relationship is determined based on consistency of the significant contractual features of the transactions ("Critical Terms Match"). In this regard, the Group performs a qualitative assessment. Hedge ineffectiveness may possibly arise when the timing of future transactions deviates from the original assumptions or the credit risk of the counterparties of the hedging instrument changes.

The following table presents the effects of the hedging instruments related to exchange rate risks on the financial position and performance of the Group:

| Currency | Carrying amount (asset) as of Dec. 31, 2018 | Carrying amount (liability) as of Dec. 31, 2018 | Hedge ratio | Change in value of hedging instruments | Change in value of hedged item | Nominal amount | Maturity: 1-6 months | Maturity: 7-12 | Average exercise price |
|----------|---|---|-------------|--|--------------------------------|-----------------------------|----------------------|----------------|------------------------|
| | € in K | € in K | | € in K | € in K | in K of respective currency | | | |
| USD | 5,554 | -661 | 100% | 4,893 | 4,893 | 137,000 | 100,000 | 37,000 | 1.14 |
| CHF | 5 | 0 | 100% | 5 | 5 | 3,000 | 3,000 | 0 | 1.13 |
| CAD | 194 | 0 | 100% | 194 | 194 | 8,000 | 6,000 | 2,000 | 1.52 |
| JPY | 0 | -302 | 100% | -302 | -302 | 1,050,000 | 600,000 | 450,000 | 130.33 |
| GBP | 16 | 0 | 100% | 16 | 16 | 3,500 | 3,500 | 0 | 0.90 |

Hedging instruments that have a positive fair value are shown in the line item "financial assets (non-current)" or "other financial assets (current)" in the statement of financial position. Hedging instruments that have a negative fair value are shown in the line item "other financial liabilities (non-current)" or "other financial liabilities (current)" in the statement of financial position.

The amounts that are recognized in the reporting period in connection with the cash flow hedges in other comprehensive income as well as those amounts that were reclassified from other comprehensive income to profit or loss (in the line item "sales revenue") are presented in the statement of other comprehensive income and the statement of changes in equity.

If the U.S. dollar would have depreciated 10% against the euro, the equity would have increased by €10.7 million (2017: increase by €13.2 million), the

impact on the result would have been +€1.6 million (2017: €0.1 million). Vice versa, if the U.S. dollar would have appreciated 10% against the euro, the resulting impact on the result would have been -€2.0 million (2017: -€0.1 million) and on the other comprehensive income -€13.1 million (2017: -€16.1 million).

36. Interest Risk Management

Sartorius Stedim Biotech is mainly financed through its parent company Sartorius AG. This major loan is taken out at variable interest rates; therefore the Group continues to be exposed to interest rate risks. To control the interest risk, an appropriate ratio between fixed and variable loans is maintained. As of December 31, 2018, the Group has no open interest rate derivative contracts to hedge the risk of increasing interest rates.

As of Dec. 31, 2018 the raised loans with variable interest rates amount to approx. €95 million.

If the market interest rate had been 1.0 percentage point higher, the interest expenses in the statement of profit or loss would have been €0.95 million (2017: €0.8 million) higher.

With regard to a decrease in interest rates a base interest rate of 0% has been considered. The impact on the financial result would have been €0.0 million (2017: –€0.2 million).

37. Liquidity Risk Management

The maturity of the financial liabilities excluding derivative financial instruments shows the following pattern:

| | Carrying amount Dec. 31, 2017 € in K | Cash Flow Dec. 31, 2017 € in K | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K |
|---|---|-----------------------------------|--------------------|-----------------------|---------------------|
| Loans and borrowings | 142,183 | 142,723 | 96,224 | 46,499 | 0 |
| Finance Leases | 17,487 | 32,825 | 1,900 | 7,558 | 23,368 |
| Trade payables | 77,211 | 77,211 | 77,211 | 0 | 0 |
| Other liabilities (excluding derivatives) | 63,428 | 68,965 | 33,036 | 35,780 | 150 |
| Financial Liabilities | 300,309 | 321,724 | 208,370 | 89,837 | 23,518 |

| | Carrying amount Dec. 31, 2018 € in K | Cash Flow Dec. 31, 2018 € in K | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K |
|---|---|--------------------------------------|--------------------|-----------------------|---------------------|
| Loans and borrowings | 132,942 | 133,191 | 89,997 | 43,194 | 0 |
| Finance Leases | 16,693 | 33,100 | 1,988 | 9,374 | 21,738 |
| Trade payables | 103,568 | 103,568 | 103,568 | 0 | 0 |
| Other liabilities (excluding derivatives) | 21,855 | 22,208 | 17,515 | 4,542 | 150 |
| Financial Liabilities | 275,059 | 292,068 | 213,069 | 57,111 | 21,888 |

The cash flows shown in the above tables include the undiscounted expected payments in connection with the respective financial liabilities including the associated interest payments based on the interest rates as of the balance sheet date.

The loans and borrowings include the loan raised from the parent company Sartorius AG. The other liabilities include the liability for the phantom units in AllPure.

The following tables illustrate the liquidity analysis for derivative financial instruments based on undiscounted cash flows:

| | Carrying amount Dec. 31, 2017 € in K | Cash Flow Dec. 31, 2017 € in K | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K |
|--------------------------|---|-----------------------------------|--------------------|-----------------------|---------------------|
| Gross fulfilment | | | | | |
| Forward contracts | 1,376 | 1,375 | 1,375 | 0 | 0 |
| Payment obligation | | 22,308 | 22,308 | 0 | 0 |
| Payment claim | | -20,932 | -20,932 | 0 | 0 |
| Net fulfilment | | | | | |
| Interest rate swaps | 552 | 552 | 552 | 0 | 0 |
| Derivatives | 1,928 | 1,927 | 1,927 | 0 | 0 |

| | Carrying amount Dec. 31, 2018 € in K | Cash Flow Dec. 31, 2018 € in K | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K |
|--------------------------|---|--------------------------------------|--------------------|-----------------------|---------------------|
| Gross fulfilment | | | | | |
| Forward contracts | 1,344 | 1,344 | 1,344 | 0 | 0 |
| Payment obligation | | 42,177 | 42,177 | | |
| Payment claim | | -40,833 | -40,833 | | |
| Derivatives | 1,344 | 1,344 | 1,344 | 0 | 0 |

The Group controls liquidity risks by maintaining credit lines and additional facilities with banks, by continuously tracking the forecasted and actual cash flows and by managing the maturity profiles of financial assets and liabilities. It is not expected that

the cash outflows will occur at significantly different times or at significantly different amounts.

The table below provides an overview of the credit lines available on the reporting date:

| | Credit line at Dec. 31, 2017 | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K | Interest rate | Credit line used at Dec. 31, 2017 | Credit line unused as of Dec. 31, 2017 |
|------------------------|---------------------------------|--------------------|-----------------------|---------------------|---------------|---|--|
| Loan from Sartorius AG | 300,000 | 0 | 300,000 | 0 | Variable | 109,481 | 190,519 |
| Bilateral loans | 9,375 | 2,344 | 7,031 | 0 | fix | 9,375 | 0 |
| Bilateral credit line | 31,335 | 31,335 | 0 | 0 | Variable | 19,235 | 12,100 |
| Total | 340,710 | 33,679 | 307,031 | 0 | | 138,091 | 202,619 |

| | Credit line at Dec. 31, 2018 | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K | Interest rate | Credit line used at Dec. 31, 2018 | Credit line unused as of Dec. 31, 2018 |
|------------------------|---------------------------------|--------------------|-----------------------|---------------------|---------------|---|--|
| Loan from Sartorius AG | 309,601 | 0 | 309,601 | 0 | Variable | 94,500 | 215,101 |
| Bilateral loans | 6,250 | 3,125 | 3,125 | 0 | fix | 6,250 | 0 |
| Bilateral credit line | 23,163 | 23,163 | 0 | 0 | Variable | 16,750 | 6,413 |
| Total | 339,014 | 26,288 | 312,726 | 0 | | 117,500 | 221,514 |

38. Credit Risk Management

Credit risk is the risk of financial loss to the Group if a counterparty to a financial instrument fails to meet its contractual obligations. Credit risk arises principally from cash and cash equivalents and trade receivables. In addition to that, the Group is exposed to credit risks arising from derivative financial instruments with positive fair values and, to a low extent, from contractual cash flows from debt securities.

Credit risk is controlled centrally for the Group by the Treasury Management unit. For counterparties such as banks and financial institutions the creditworthiness is continuously monitored in order to recognize increases in credit risks at an early stage. If no new information is obtained, the Group assumes that the related financial assets still have a low credit risk.

Customers are assigned risk limits that principally depend on the business volume, past experience and the financial position of the customer. Compliance with the limits is regularly reviewed by the management responsible. In some cases, the Group receives advance payments in order to avoid credit risks. There are no significant concentrations of credit risks from individual customers or regions.

For some trade receivables the Group possibly has collateral such as guarantees that can be used within the contractual agreements in case the counterparty does not meet its contractual payment obligations.

Impairment of Trade Receivables and Contract Assets

The new impairment model of IFRS 9, recognition of expected credit losses, is of particular relevance for the Group's trade receivables and contract assets according to IFRS 15. The Group applies the simplified approach according to IFRS 9 to trade receivables and contract assets. Accordingly, lifetime expected credit losses are recognized for these assets. The starting point of the new impairment model is an analysis of the actual historical credit loss rates. These are adjusted, taking into consideration forward-looking information and the effects of current changes in the macroeconomic environment, if significant. Due to the immaterial level of historical credit losses the Group currently determines the expected credit losses for the Group's portfolio of trade receivables as a whole. However, historical loss rates are analyzed regularly in more detail in order to apply different loss rates to different portfolios, where appropriate.

The contract assets are related to projects for typical customers of the Group. Therefore, it is assumed that the loss rates applied to trade receivables are appropriate approximations for the loss rates of the contract assets. Accordingly, there is no further differentiation between trade receivables and contract assets.

On this basis, the allowances for trade receivables and contract assets were determined as follows as of December 31, 2018 and as of the date of initial application of IFRS 9:

| December 31, 2018 € in K | Not due | 1 – 30 days overdue | 31 – 60 days overdue | 61 – 90 days overdue | More than 90 days overdue | Total |
|--|---------|------------------------|-------------------------|-------------------------|------------------------------|--------------|
| Gross carrying amount of trade receivables | 133,707 | 39,781 | 17,502 | 6,115 | 22,843 | 219,947 |
| Gross carrying amount of contract assets | 3,902 | 0 | 0 | 0 | 0 | 3,902 |
| Impairment loss allowance | 77 | 24 | 49 | 280 | 3,188 | 3,617 |
| January 1, 2018 € in K | Not due | 1 – 30 days overdue | 31 – 60 days overdue | 61 – 90 days overdue | More than 90 days overdue | Total |
| Gross carrying amount of trade receivables | 149,266 | 27,460 | 10,757 | 4,180 | 15,419 | 207,083 |
| Gross carrying amount of contract assets | 6,967 | 0 | 0 | 0 | 0 | 6,967 |
| Impairment loss allowance | 78 | 18 | 108 | 4 | 1,903 | 2,111 |

The expected credit losses are determined based on a loss rate of 0.05%. In addition, impairments are determined on the basis of individual assessments. Days overdue are one essential criterion in this context. A default is generally presumed when there is no

reasonable expectation of recovering a financial asset. In such a case, the respective receivables are derecognized.

The movements in the allowance for impairment in respect of trade receivables and contract assets in the reporting period are presented below:

| | € in K |
|--|---------------|
| Balance at January 1, 2018 under IAS 39 | -2,433 |
| Adjustment on initial application of IFRS 9 | 321 |
| Balance at January 1, 2018 under IFRS 9 | -2,111 |
| Net remeasurement of loss allowance recognized in profit or loss in the reporting period | -2,041 |
| Derecognition and consumption | 39 |
| Recoveries of amounts previously impaired | 474 |
| Foreign currency translation differences | 23 |
| Balance at December 31, 2018 under IFRS 9 | -3,617 |

Impairment of Other Financial Assets

Besides trade receivables, cash and cash equivalents are the most material financial assets on the Group's statement of financial position as of the reporting date, December, 2018. The expected credit losses are monitored at regular intervals. Due to the high creditworthiness of the counterparties and the short maturities, the impairment which would have to be recognized for these financial assets is immaterial. Therefore, no impairment is recognized for cash and cash equivalents.

For the other financial assets measured at amortized cost no impairment is recognized as of December 31, 2018 for the twelve months of expected credit losses due to immaterial historical credit losses. In the event of a significant increase in credit risk, which is generally presumed when a payment is more than 30 days past due, the lifetime expected credit losses are

recognized for the respective financial asset. A default is generally presumed when there is no reasonable expectation of recovering a financial asset. This is generally presumed when payments are more than 90 days past due. As of the reporting date, there are no indications of increases in credit risk to a material extent. The carrying amounts of the financial assets reflect the maximum credit loss for these assets at the end of the reporting period.

39. Other Risks Associated with Financial Instruments

As of the reporting date, the Sartorius Stedim Biotech Group has not been exposed to the risk of volatility in share prices.

40. Share-based Payments

On the level of Sartorius Stedim Biotech's majority shareholder Sartorius AG, share-based payments exist in the form of so-called phantom stock units. Under this plan the respective board member is granted a certain number of phantom stocks each year that represent an agreed amount of money. The exercise of these stocks is not possible before four years and is depending on certain requirements regarding the performance of the Sartorius AG shares.

When the stocks are paid out the amount is based on the share price at the exercise date. The payment is capped at an amount of 2.5 times the share price at the time these virtual options were granted. For further details please refer to the Remuneration Report.

The fair value of the phantom stock units is disclosed as follows:

| | Number of phantom stock units | Subscription price in € | Fair value when granted on Jan. 1 of the particular year € in K | Fair value at year-end on Dec. 31, 2018 € in K | Paid out € in K | Exercisable |
|---|-------------------------------|-------------------------|---|--|-----------------|-------------|
| Tranche of phantom stock units for 2014 | 8,032 | 21.01 | 169 | 0 | 422 | |
| Tranche of phantom stock units for 2015 | 7,360 | 24.70 | 182 | 454 | 0 | no |
| Tranche of phantom stock units for 2016 | 3,484 | 57.41 | 200 | 359 | 0 | no |
| Tranche of phantom stock units for 2017 | 2,950 | 70.51 | 208 | 281 | 0 | no |
| Tranche of phantom stock units for 2018 | 2,685 | 80.32 | 216 | 239 | 0 | no |
| Total | 24,511 | | 974 | 1,333 | 422 | |

Other Disclosures

The consolidated financial statements were prepared on a going concern basis.

Material Events after the Reporting Date

No material events occurred after the reporting date.

Number of Employees

The average workforce employed during the year 2018 was 5,412 (4,991 in 2017).

41. Related Parties

General

The majority shareholder of Sartorius Stedim Biotech S.A. is Sartorius AG, which holds a controlling stake in the company of 74.3% in equity capital – and 84.5% of the voting rights. The Sartorius Group itself is organized in two divisions: Bioprocess Solutions (mainly run by the Sartorius Stedim Biotech Group) and Lab Products & Services (mainly run by the other companies of Sartorius Group). This structure leads to the fact that the Group holds two subsidiaries in most of the countries and these companies partially share space, staff and other resources. Furthermore, the German group companies carry out various central functions and accordingly deliver services to the worldwide entities (e.g. IT support). The company Sartorius Corporate Administration GmbH, a 100% subsidiary of Sartorius AG has incorporated numerous Group functions like Group Finance, HR, IT, Investor Relations and Legal. These services are charged within the Group and to a significant extent also to Sartorius Stedim Biotech.

The described structures lead to various relations and transactions with related parties. Transactions between Sartorius Stedim Biotech S.A. and its subsidiaries (presented in Note 7), which are related parties of the company, have been eliminated on consolidation and are not disclosed under this Note. Details of transactions between the Group and other related parties, belonging to the Sartorius Group, are disclosed below.

Sales, Purchases and Commissions

In certain business areas members of the Sartorius Group act as contract manufacturers for the Sartorius Stedim Biotech Group and vice versa. The respective transactions are carried out at arms' length principles and are disclosed in the table below as "sales revenue" and "purchases".

| | Sales revenue 2018 € in K | Purchases 2018 € in K |
|------------------------------------|---------------------------------|-----------------------------|
| Related parties of Sartorius Group | 72,221 | 7,061 |
| | Sales revenue 2017 € in K | Purchases 2017 € in K |
| Related parties of Sartorius Group | 60,292 | 5,085 |

Certain product groups of the Sartorius Stedim Biotech portfolio are sold through the sales force of other Sartorius entities. For the arranging of the sale the Sartorius Stedim Biotech Group has paid commissions in the amount of €0.4 million (€1.1 million in 2017). These commissions are typically calculated as a percentage of the generated sales revenue.

Management Fees and Other Shareholder Costs

One of Sartorius Stedim Biotech S.A.'s board members is also member of the Sartorius AG Executive Board and is paid by the German parent company. For his services for Sartorius Stedim Biotech a portion of his remuneration is charged to Sartorius Stedim Biotech S.A. (€0.7 million in 2018; €0.8 million in 2017) and charged to Sartorius Stedim Biotech GmbH (€0.7 million in 2018; €0.8 million in 2017).

Other shareholder functions like Group Financial Reporting, Compliance and Investor Relations are performed by the above mentioned Sartorius Corporate Administration GmbH in Germany. These services have been charged to Sartorius Stedim Biotech S.A. in the amount of €0.8 million (2017: €0.8 million).

Shareholder Loan

As described in note 30 the Sartorius Stedim Biotech Group has raised a loan from its parent company Sartorius AG with a credit volume of approx. €300 million and a current utilization of approx. €95 million (2017: €109 million). The interest charged is based on a variable interest rate plus an arms'-length credit margin.

Administration Charges and Shared Costs

As described above the companies in most of the countries share certain functions and costs. The underlying contracts include mainly subleases for office space and central administrative functions, such as accounting and controlling, human resource

management and IT. In this respect, the relevant companies charge rent, salaries, social security costs and other expenses for such services, as well as a pro-rated profit margin for the services they provide.

The most significant contract in this context is the one between Sartorius Stedim Biotech GmbH, Germany, and Sartorius Corporate Administration GmbH. This company provides all central service and administrative functions to Sartorius Stedim Biotech GmbH and other Group companies. The calculation for services fees typically includes a surcharge of 3% on total costs. 3% is a surcharge compliant with arm's length principles for routine tasks, following OECD and EU guidelines. In 2018, services for approx. €53.4 million were provided to Sartorius Stedim Biotech GmbH (€40.1 million in 2017). This amount covers the following functions:

- Marketing Communication, e-Business, Business Development
- Environment, Health & Security, Factory Maintenance
- Finance, Human Resources, Information Technology
- Central Services & General Organization.

During the reporting period, the Sartorius Stedim Biotech Group acquired a perpetual right to use the Sartorius global process template for some of its entities for a total amount of €19.9 million. This transaction enhances the independence of the Sartorius Stedim Biotech Group from Sartorius Corporate Administration GmbH and will lead to a corresponding reduction in future IT service charges.

Compensation of Key Management Personnel

In 2017 and 2018, the Executive Board Management received the following remuneration:

| | Total € in K | Short-term benefits € in K | Post-employment benefits € in K | Other long-term benefits € in K | Termination benefits € in K | Share-based payments € in K |
|--------------------------|-----------------|----------------------------------|---------------------------------------|---------------------------------------|-----------------------------------|-----------------------------------|
| 2018¹⁾ | 2,522 | 1,333 | 257 | 189 | 0 | 743 |
| 2017 ¹⁾ | 3,056 | 1,210 | 258 | 184 | 0 | 1,404 |

¹⁾ For more information please refer to the chapter Corporate Governance (See pages 57 to 87)

Statutory Auditors' Report on the Consolidated Financial Statements

This is a translation into English of the statutory auditors' report on the consolidated financial statements of the Company issued in French and it is provided solely for the convenience of English speaking users.

This statutory auditors' report includes information required by European regulation and French law, such as information about the appointment of the statutory auditors or verification of the information concerning the Group presented in the management report.

This report should be read in conjunction with, and construed in accordance with, French law and professional auditing standards applicable in France

For the year ended 31 December 2018

To the Shareholders of SARTORIUS STEDIM BIOTECH S.A.,

Opinion

In compliance with the assignment entrusted to us by your shareholders' meetings, we have audited the accompanying consolidated financial statements of SARTORIUS STEDIM BIOTECH S.A. for the year ended 31 December 2018.

In our opinion, the consolidated financial statements give a true and fair view of the assets and liabilities and of the financial position of the Group as at 31 December 2018 and of the results of its operations for the year then ended in accordance with International Financial Reporting Standards as adopted by the European Union.

The audit opinion expressed above is consistent with our report to the Audit Committee.

Basis for Opinion

Audit Framework

We conducted our audit in accordance with professional standards applicable in France. We believe that the audit evidence we have obtained is sufficient and appropriate to provide a basis for our opinion.

Our responsibilities under those standards are further described in the Statutory Auditors' Responsibilities for the Audit of the Consolidated Financial Statements section of our report.

Independence

We conducted our audit engagement in compliance with independence rules applicable to us, for the period from 1 January 2018 to the date of our report and specifically we did not provide any prohibited non-audit services referred to in Article 5(1) of Regulation (EU) No 537/2014 or in the French Code of ethics (code de déontologie) for statutory auditors.

Justification of Assessments – Key Audit Matters

In accordance with the requirements of Articles L.823 - 9 and R.823 - 7 of the French Commercial Code (code de commerce) relating to the justification of our assessments, we inform you of the key audit matters relating to risks of material misstatement that, in our professional judgment, were of most significance in our audit of the consolidated financial statements of the current period, as well as how we addressed those risks.

These matters were addressed in the context of our audit of the consolidated financial statements as a whole, and in forming our opinion thereon, and we do not provide a separate opinion on specific items of the consolidated financial statements.

Goodwill valuation – Impairment test

Identified risk

As of 31 December 31 2018, goodwill amounts to 384.7 M€, or 25% of total consolidated assets.

As described in Note 5 to the consolidated financial statements, SARTORIUS STEDIM BIOTECH is an "integrated solution provider" for its customers, and as a result there is only one operating segment from a product and customer perspective: "Biopharma ". In addition, as indicated in Note 16 to the consolidated financial statements, because of the interdependence of the market in which your group operates, the

lowest level at which goodwill can be allocated is the Biopharma segment. The goodwill has therefore been fully allocated to the Cash Generating Units (C.G.U.) corresponding to the Biopharma segment.

Goodwill is subject to an annual impairment test and whenever there are indicators of impairment in accordance with the methods and assumptions described in Notes 3 and 16 to the consolidated financial statements. In particular, in view of what has been described above, the impairment test is carried out at the level of the Biopharma segment.

We considered that the determination of the value of goodwill is a key point of our audit given its significance in the consolidated financial statements of your group, and because the determination of the recoverable amount taken into account in the impairment test on the basis of the value in use of the C.G.U. requires the use of estimates and assumptions (including future cash flows, perpetual revenue growth rates and the discount rate) that require significant management judgment.

Responses obtained during our audit

We obtained the impairment test from the C.G.U. group corresponding to the Biopharma segment as well as the forecasts underlying the calculation (4-year plan).

We reviewed the compliance of the company's methodology in accordance with applicable accounting standards.

We also performed a critical review of how the company has implemented this methodology, including the following procedures:

- Assess the reasonableness of the key assumptions used to determine the cash flow of the Biopharma segment as well as that used for the perpetual growth rate;
- Assess, with the support of our valuation specialists, the discount rate used by management. We compared this rate with our own estimates and analysed its various constituent elements;
- Check the arithmetical accuracy of the impairment test performed by your group.

We have also obtained and assessed the sensitivity analyses carried out by management, as shown in Note 16 of the notes to the consolidated financial statements. As a result, we were able to verify that

only an extremely large change in the main assumptions could lead to the recognition of an impairment of goodwill.

Lastly, we verified the appropriateness of the information provided in Notes 3 and 16 to the consolidated financial statements.

Specific verifications

As required by French law and regulations, we have also verified in accordance with professional standards applicable in France the information pertaining to the Group presented in the management report of the Board of Directors.

We have no matters to report as to its fair presentation and its consistency with the consolidated financial statements

Report on Other Legal and Regulatory Requirements

Appointment of the Statutory Auditors

We were appointed as statutory auditors of SARTORIUS STEDIM BIOTECH S.A. by the annual general meeting held on 7 April 2015 for KPMG S.A. and on 19 May 2006 for Deloitte & Associés.

As at 31 December 2018, KPMG S.A. was in its 4th year of the audit mandate without interruption and Deloitte & Associés was in its 13th year.

Responsibilities of Management and Those Charged with Governance for the Consolidated Financial Statements

Management is responsible for the preparation and fair presentation of the consolidated financial statements in accordance with International Financial Reporting Standards as adopted by the European Union and for such internal control as management determines is necessary to enable the preparation of consolidated financial statements that are free from material misstatement, whether due to fraud or error.

In preparing the consolidated financial statements, management is responsible for assessing the Company's ability to continue as a going concern, disclosing, as applicable, matters related to going concern and using the going concern basis of

accounting unless it is expected to liquidate the Company or to cease operations.

The Audit Committee is responsible for monitoring the financial reporting process and the effectiveness of internal control and risks management systems and where applicable, its internal audit, regarding the accounting and financial reporting procedures.

The consolidated financial statements were approved by the Board of Directors.

Statutory Auditors' Responsibilities for the Audit of the Consolidated Financial Statements

Objectives and audit approach

Our role is to issue a report on the consolidated financial statements. Our objective is to obtain reasonable assurance about whether the consolidated financial statements as a whole are free from material misstatement. Reasonable assurance is a high level of assurance, but is not a guarantee that an audit conducted in accordance with professional standards will always detect a material misstatement when it exists. Misstatements can arise from fraud or error and are considered material if, individually or in the aggregate, they could reasonably be expected to influence the economic decisions of users taken on the basis of these consolidated financial statements.

As specified in Article L.823-10-1 of the French Commercial Code (code de commerce), our statutory audit does not include assurance on the viability of the Company or the quality of management of the affairs of the Company.

As part of an audit conducted in accordance with professional standards applicable in France, the statutory auditor exercises professional judgment throughout the audit and furthermore:

- Identifies and assesses the risks of material misstatement of the consolidated financial statements, whether due to fraud or error, designs and performs audit procedures responsive to those risks, and obtains audit evidence considered to be sufficient and appropriate to provide a basis for his opinion. The risk of not detecting a material misstatement resulting from fraud is higher than for one resulting from error, as fraud may involve collusion, forgery, intentional omissions, misrepresentations, or the override of internal control. Obtains an understanding of internal control relevant to the audit in order to design audit

procedures that are appropriate in the circumstances, but not for the purpose of expressing an opinion on the effectiveness of the internal control.

- Obtains an understanding of internal control relevant to the audit in order to design audit procedures that are appropriate in the circumstances, but not for the purpose of expressing an opinion on the effectiveness of the internal control.
- Evaluates the appropriateness of accounting policies used and the reasonableness of accounting estimates and related disclosures made by management in the consolidated financial statements.
- Assesses the appropriateness of management's use of the going concern basis of accounting and, based on the audit evidence obtained, whether a material uncertainty exists related to events or conditions that may cast significant doubt on the Company's ability to continue as a going concern. This assessment is based on the audit evidence obtained up to the date of his audit report. However, future events or conditions may cause the Company to cease to continue as a going concern. If the statutory auditor concludes that a material uncertainty exists, there is a requirement to draw attention in the audit report to the related disclosures in the consolidated financial statements or, if such disclosures are not provided or inadequate, to modify the opinion expressed therein.
- Evaluates the overall presentation of the consolidated financial statements and assesses whether these statements represent the underlying transactions and events in a manner that achieves fair presentation.
- Obtains sufficient appropriate audit evidence regarding the financial information of the entities or business activities within the Group to express an opinion on the consolidated financial statements. The statutory auditor is responsible for the direction, supervision and performance of the audit of the consolidated financial statements and for the opinion expressed on these consolidated financial statements.

Report to the Audit Committee

We submit a report to the Audit Committee which includes in particular a description of the scope of the audit and the audit program implemented, as well as the results of our audit. We also report, if any, significant deficiencies in internal control regarding

the accounting and financial reporting procedures that we have identified.

Our report to the Audit Committee includes the risks of material misstatement that, in our professional judgment, were of most significance in the audit of the consolidated financial statements of the current period and which are therefore the key audit matters, that we are required to describe in this audit report.

We also provide the Audit Committee with the declaration provided for in Article 6 of Regulation (EU) N° 537/2014, confirming our independence within the meaning of the rules applicable in France such as they are set in particular by Articles L.822-10 to L.822-14 of the French Commercial Code (code de commerce) and in the French Code of Ethics (code de déontologie) for statutory auditors. Where appropriate, we discuss with the Audit Committee the risks that may reasonably be thought to bear on our independence, and the related safeguards.

Marseille, 8 February 2019

The Statutory Auditors

French original signed by

KPMG Audit
A division of KPMG S.A

John Evans

Deloitte & Associés

Vincent Gros

Annual Financial Statements of
Sartorius Stedim Biotech S.A. and Notes

05

Annual Financial Statements

Parent Company Balance Sheet: Assets (in thousands of €)

| | Gross at Dec. 31, 2018 | Depreciation, amortization and provisions Dec. 31, 2018 | Net at Dec. 31, 2018 | Net at Dec. 31, 2017 |
|------------------------------------|---------------------------|--|-------------------------|-------------------------|
| Intangible assets | 552 | -117 | 435 | 463 |
| Property, plant and equipment | 19,176 | -12,022 | 7,155 | 12,384 |
| Financial investments | 128,749 | 0 | 128,749 | 128,138 |
| Total non-current assets | 148,478 | -12,139 | 136,339 | 140,985 |
| Inventories and work in progress | 0 | 0 | 0 | 0 |
| Trade receivables to third parties | 0 | 0 | 0 | 0 |
| Other receivables | 66,066 | 0 | 66,066 | 23,812 |
| Deposits and cash equivalents | 62 | | 62 | 96 |
| Total current assets | 66,128 | 0 | 66,128 | 23,908 |
| Prepaid expenses | 231 | | 231 | 49 |
| Currency translation adjustment | 482 | | 482 | 0 |
| Total assets | 215,319 | -12,139 | 203,180 | 164,942 |

Parent Company Balance Sheet: Equity and Liabilities (in thousands of €)

| | At Dec. 31, 2018 | At Dec. 31, 2017 |
|---|------------------|------------------|
| Share capital | 18,436 | 18,436 |
| Share premium | 12,609 | 12,609 |
| Reserves | 2,434 | 2,434 |
| Retained earnings carried forward | 34,346 | 27,286 |
| Profit for the period | 49,521 | 49,463 |
| Regulated provisions | 4,088 | 4,088 |
| Total equity | 121,434 | 114,316 |
| Provisions for liabilities and charges | 482 | 0 |
| Total provisions for liabilities and charges | 482 | 0 |
| Loans and borrowings | 0 | 0 |
| Trade payables | 881 | 455 |
| Tax and social charges payable | 63 | 96 |
| Liabilities for non-current assets | 313 | 0 |
| Other liabilities | 79,705 | 50,075 |
| Total liabilities | 80,962 | 50,626 |
| Currency translation adjustment | 301 | 0 |
| Total equity and liabilities | 203,180 | 164,942 |

Parent Company: Income Statement (in thousands of €)

| | At Dec. 31, 2018 | At Dec. 31, 2017 |
|---|------------------|------------------|
| Sales revenue | 1,999 | 2,198 |
| Inventory movements | 0 | 0 |
| Capitalized production costs | 0 | 0 |
| Depreciation or amortization reversals | 0 | 0 |
| Other operating income and expense reallocation | 1,490 | 130 |
| Purchases consumed | 0 | 0 |
| External charges for services | -3,775 | -3,453 |
| Tax and duties | -411 | -830 |
| Personnel costs | 0 | 0 |
| Additions to amortization, depreciation and provision | -816 | -825 |
| Other operating expenses | -857 | -417 |
| Operating profit (EBIT) | (2,371) | (3,197) |
| Net financing income (expense) | 48,576 | 47,108 |
| Profit (loss) from ordinary activities | 46,205 | 43,911 |
| Exceptional income (expense) | 0 | 0 |
| Income tax | 3,316 | 5,552 |
| Net profit (loss) | 49,521 | 49,463 |

1. Materiel Events during the Year

None

2. Materiel Events after the Reporting date

None

3. Accounting Principles and Methods

The parent company's financial statements for the year ended December 31, 2018, were prepared and presented in accordance with French accounting rules in compliance with the principles of prudence, reporting on distinct financial years and the presumption of going concern.

The annual financial statements have been prepared in accordance with the clauses of the CRC Regulation 2014-03 of September 8, 2014 on the French chart of accounts.

Sartorius Stedim Biotech S.A. is listed in Compartment A of the Euronext Paris Stock Exchange (ISIN FR code 0000053266) and also prepares consolidated financial statements in accordance with IFRS standards, as adopted by the European Union on December 31, 2018. Sartorius Stedim Biotech S.A. is consolidated by Sartorius A.G.

3.1. Non-current Assets

Non-current intangible and tangible assets are valued at their acquisition costs, excluding costs incurred for their acquisition.

For intangible assets and property, plant and equipment, the Company applied the French Regulation CRC No. 2002-10, recodified by Article 2-4 of Regulation CRC No. 2004-06 relative to the amortization, depreciation and impairment of assets according to the "Component approach."

3.1.1. Intangible Assets

The following is thus valued under this heading: incorporation costs, patents and software.

All these assets are amortized on a straight-line basis using the following indicative useful lives:

- Incorporation costs: One to five years
- Software: One to three years
- Patents: Twenty years
- Leasehold: Eighteen years (Based on the period of use).

As part of the implementation of integrated software, the direct labor costs concerned are included in the

amount capitalized as cost, as a function of the time elapsed.

Intangible assets are valued at acquisition cost less amortization and impairments reported, on an ongoing basis.

3.1.2. Property, Plant and Equipment

Property, plant and equipment (PPE) are recognized at their acquisition value, including the installation cost of these assets.

Depreciation is calculated over the standard and economic life of the assets using the straight-line method.

All these non-current assets are depreciated on a straight-line basis using the following indicative periods of use:

- Buildings: Twenty to forty years
- Improvements, fixtures and fittings:
Ten to fifteen years
- Plant and equipment: Four to ten years
- Office and IT equipment: Three to five years
- Motor vehicles: Four to five years

Property, plant and equipment are valued at acquisition cost less depreciation and impairments reported, on an ongoing basis.

3.1.3. Financial Investments

Investments relate mainly to shareholdings in subsidiaries and other treasury shares held within the scope of the share buyback program; they are recorded at their acquisition cost, including fees linked to their acquisition.

An impairment provision may be recorded to take into account, in particular, either the stock exchange price or the underlying assets of these subsidiaries, their financial position and their prospects.

Shareholdings in subsidiaries are subject to impairment tests.

3.2. Receivables and Payables

Receivables and payables are recorded at their nominal value.

Receivables whose collection is doubtful are subject to a provision for doubtful debts.

4. Non-Current Assets (in thousands of €)

4.1. Intangible Assets

| | At Dec. 31, 2017 | Increase in 2018 | Decrease in 2018 | At Dec. 31, 2018 |
|-------------------------------|------------------|------------------|------------------|------------------|
| Gross values | | | | |
| Incorporation costs | 4 | 0 | 0 | 4 |
| Patents | 0 | 0 | 0 | 0 |
| Software, licenses | 0 | 0 | 0 | 0 |
| Business goodwill | 548 | 0 | 0 | 548 |
| Intangible assets in progress | 0 | 0 | 0 | 0 |
| Total | 552 | 0 | 0 | 552 |
| Amortization and depreciation | 89 | 28 | 0 | 117 |
| Net amount | 463 | -28 | 0 | 435 |

4.2. Property, Plant and Equipment

| | At Dec. 31, 2017 | Increase in 2018 | Decrease in 2018 | At Dec. 31, 2018 |
|---|------------------|------------------|------------------|------------------|
| Gross values | | | | |
| Land | 496 | 0 | 0 | 496 |
| Buildings | 15,561 | 154 | 0 | 15,715 |
| Plant and equipment | 0 | 0 | 0 | 0 |
| Other | 2,127 | 145 | 0 | 2,272 |
| Property, plant and equipment in progress | 5,434 | 559 | -5,300 | 693 |
| Total | 23,618 | 858 | -5,300 | 19,176 |
| Amortization and depreciation | At Dec. 31, 2017 | Addition | Release | At Dec. 31, 2018 |
| Buildings | 10,388 | 516 | 0 | 10,904 |
| Plant and equipment | 0 | 0 | 0 | 0 |
| Other | 846 | 271 | 0 | 1,117 |
| Total | 11,234 | 788 | 0 | 12,022 |
| Property, plant and equipment, net | 12,384 | 70 | -5,300 | 7,154 |

The increase in tangible assets includes fixtures and fittings for a net amount of €299 K and assets under construction for an amount of €559 K.

4.3. Financial Investments

| Investments | At Dec. 31, 2017 | Increase in 2018 | Decrease in 2018 | At Dec. 31, 2018 |
|-------------------------------|------------------|------------------|------------------|------------------|
| Shareholdings | 127,977 | 0 | 0 | 127,977 |
| Write-down of shareholdings | 0 | 0 | 0 | 0 |
| Deposits and guarantees | 160 | 0 | -25 | 135 |
| Treasury shares | 0 | 637 | 0 | 637 |
| Write-down of treasury shares | 0 | 0 | 0 | 0 |
| Other non-current assets | 0 | 0 | 0 | 0 |
| Total | 128,137 | 637 | -25 | 128,749 |

The following is included under "Financial investments":

- 99.99% of the share capital of Sartorius Stedim Bioprocess SARL, a Tunisian company;
- 100% of the share capital of Sartorius Stedim Biotech GmbH, a company governed by German law, following the merger of the Sartorius and the Stedim Groups in June 2007;
- 100% of the share capital of Sartorius Stedim Aseptics S.A., a French company acquired in 2004;

- 100% of the share capital of Sartorius Stedim FMT S.A.S., a French company created in connection with the Contribution Assets transfer in 2013;

- Other investments: €1.0 K.

The amount now corresponds to the share of Sartorius Stedim Biotech in the Russian company Sartorius Stedim RUS.

A new liquidity contract between the entity Sartorius Stedim Biotech S.A. and the brokerage company Gilbert Dupont began the 20th of April 2018. Therefore, Sartorius Stedim Biotech holds 3,416 shares of SSB S.A. in portfolio at the closing.

5. Trade Receivables (in thousands of €)

Maturity of Receivables at Year-end (in thousands of €)

| Type of receivable | Net amount | Less than 1 year | More than 1 year |
|-----------------------------|---------------|------------------|------------------|
| Deposits and guarantees | 772 | 772 | |
| Non-current assets | 772 | 772 | 0 |
| Advance payments on account | 0 | 0 | 0 |
| Trade receivables | 0 | 0 | 0 |
| Personnel | 0 | 0 | 0 |
| Social security | 0 | 0 | 0 |
| Taxes and duties | 1,784 | 1,784 | 0 |
| Group | 64,282 | 64,282 | 0 |
| Other receivables | 0 | 0 | 0 |
| Current assets | 66,066 | 66,066 | 0 |
| Prepaid expenses | 231 | 231 | 0 |
| Total receivables | 67,069 | 67,069 | 0 |

The "Group" item for receivables from Group subsidiaries (€64,282 K) relates to current account cash advances provided to Sartorius Stedim France, Sartorius Stedim Aseptics, Sartorius Stedim FMT SAS and Sartorius Stedim Bioprocess Tunisia.

The "Taxes and duties" (€1,784 K) captures primarily includes the net tax receivable relating to the tax grouping system.

6. Maturity of Liabilities at Year-end (in thousands of €)

| Type of liability | Net amount | Less than 1 year | Between 1 and 5 years | More than 5 years |
|---|---------------|------------------|-----------------------|-------------------|
| Loans and borrowings from credit institutions | | | | |
| Originally less than 2 years | 0 | 0 | 0 | 0 |
| Originally more than 2 years | 0 | 0 | 0 | 0 |
| Current bank overdrafts and accrued interest | 0 | 0 | 0 | 0 |
| Trade payables | 881 | 881 | 0 | 0 |
| - including bills of exchange | 0 | 0 | 0 | 0 |
| Advances and payments on account for orders | 0 | 0 | 0 | 0 |
| Tax and social security payable | 63 | 63 | 0 | 0 |
| Liabilities for non-current assets | 313 | 313 | 0 | 0 |
| Group and associates | 79,322 | 79,322 | 0 | 0 |
| Other | 384 | 384 | 0 | 0 |
| Total liabilities | 80,962 | 80,962 | 0 | 0 |

The "Group" item for liabilities from Group subsidiaries (€79,322 K) relates to cash-pooling liabilities and current account cash advances provided by Sartorius AG, Sartorius Stedim Biotech GmbH, Sartorius Stedim France SAS, Sartorius Stedim FMT SAS and Sartorius Stedim Aseptics SA.

Accrued expenses included in these accounts represented €411 K and concerned the following items:

| Type of expense | At Dec. 31, 2018 |
|--|------------------|
| Accrued banking charges | 0 |
| Suppliers' invoices to be received | 411 |
| Paid vacation including social charges | 0 |
| Bonuses, including social charges and profit sharing | 0 |
| Social security payable | 0 |
| Taxes payable | 0 |
| Employee profit sharing | 0 |
| Total charges payable | 411 |

7. Parent Company Statement of Changes in Equity (in thousands of €)

7.1. Equity

At December 31, 2017, the share capital was €18,436 K, comprising 92,180,190 shares of a €0.20 par value.

At December 31, 2018, the share capital is €18,436 K, comprising 92,180,190 shares of a €0.20 par value.

The Annual General Shareholders' Meeting on April 3, 2018, approved the appropriation of the net profit for the year of €49,463 K, as follows:

– Allocation to the retained earnings carried forward:
– €7,060 K

– Paid into the legal reserves: None

A dividend total of €42,403 K, or a net dividend per share of €0.46, was paid.

| | Appropriation of profit in 2017 | | | Movements 2018 | | Equity before appropriation of profit in 2018 |
|-------------------------------|---------------------------------|----------|----------------|----------------|----------------|---|
| | Before | Changes | After | Increases | Decreases | Total |
| Number of shares: | 92,180,190 | | 92,180,190 | | | 92,180,190 |
| Share capital | 18,436 | | 18,436 | | | 18,436 |
| Share premium | 0 | | 0 | | | 0 |
| Merger premium | 12,609 | | 12,609 | | | 12,609 |
| Legal reserve | 1,844 | | 1,844 | | | 1,844 |
| Other reserves | 591 | | 591 | | | 591 |
| Balance carried forward | 27,286 | 7,060 | 34,346 | | | 34,346 |
| Dividends paid | 0 | 42,403 | 42,403 | | (42,403) | 0 |
| Net profit to be appropriated | 49,463 | (49,463) | 0 | | | 0 |
| Profit for the reporting year | | | 0 | 49,521 | | 49,521 |
| Regulated provisions | 4,088 | | 4,088 | | | 4,088 |
| Total | 114,317 | 0 | 114,317 | 49,521 | -42,403 | 121,435 |

7.2. Stock Options

None

8. Risks and Provisions (in thousands of €)

8.1. Provisions

| Type of provision | Provisions at Dec. 31, 2017 | Additions 2018 | Releases 2018 | Provisions at Dec. 31, 2018 |
|---|-----------------------------|----------------|---------------|-----------------------------|
| Regulated provisions | | | | |
| Accelerated amortization and depreciation | 4,088 | 0 | 0 | 4,088 |
| Subtotal (1) | 4,088 | 0 | 0 | 4,088 |
| Provisions for liabilities and charges | | | | |
| Exchange risk | 0 | 482 | 0 | 482 |
| Other costs | 0 | 0 | 0 | 0 |
| Taxation | 0 | 0 | 0 | 0 |
| Subtotal (2) | 0 | 482 | 0 | 482 |
| Grand total | 4,088 | 482 | 0 | 4,570 |

8.2. Market Risk Exposure

Operating Cash Flow risks

At December 31, 2018, the impacts on net amount in foreign currency in current assets and liabilities are respectively of €482 K and €301 K.

Current and Future Tax Position (in thousands of €)

As of January 1, 2008, the company chose to adopt the French tax integration regime within the framework of a tax group. The lead company of this group is Sartorius Stedim Biotech S.A. The other member companies of this tax integration group for

tax relief are Sartorius Stedim Aseptics S.A., Sartorius Stedim France S.A.S. and Sartorius Stedim FMT S.A.S.

The member companies report income tax as if there were no integration tax regime. The parent corporation benefits from tax relief related to consolidating the gains and losses of the other members companies.

For 2018, the net impact according to the consolidation rules of the French tax integration regime for tax relief is an income of €3,316 K. Taking into account the tax credits not yet compensated, the company SSB holds a receivable from the State of €1,633 K.

9. Operating Income (in thousands of €)

9.1. Sales Revenue by Operating Segment

| Operating segment | At Dec. 31, 2018 | | At Dec. 31, 2017 | |
|-------------------|---------------------|-------------|---------------------|-------------|
| Services | 1,999 | 100% | 2,198 | 100% |
| Total | 1,999 | 100% | 2,198 | 100% |

9.2. Sales Revenue by Geographical Region

| Geographical region | At Dec. 31, 2018 | | At Dec. 31, 2017 | |
|--------------------------|---------------------|-------------|---------------------|-------------|
| France | 1,999 | 100% | 2,198 | 100% |
| Export | 0 | | 0 | 0% |
| EU and other countries | 0 | | 0 | |
| North American continent | 0 | | 0 | |
| Total | 1,999 | 100% | 2,198 | 100% |

The Sale revenue corresponds to the rent invoiced to the entity Sartorius Stedim FMT S.A.S. for the use of premises located in Aubagne within its operational activity.

10. Breakdown of Income Tax (in thousands of €)

| | At Dec. 31, 2018 | | | At Dec. 31, 2017 | | |
|--------------------------------|----------------------|----------------------|---------------------|----------------------|----------------------|------------------|
| | Profit before tax | Income tax charge | Profit after tax | Profit before tax | Income tax charge | Profit after tax |
| Gross taxable income | 46,205 | 0 | 46,205 | 43,911 | 2,206 | 46,117 |
| Exceptional income (expense) | 0 | 0 | 0 | 0 | 0 | 0 |
| French tax integration relief | 0 | 3,316 | 3,316 | 0 | 3,346 | 3,346 |
| Net taxable income | 46,205 | 3,316 | 49,521 | 43,911 | 5,552 | 49,463 |

11. Information on Directors' Remuneration

Remuneration allocated and paid to members of the Board of Directors as directors' meeting attendance fees amounted to €313.0K. These fees related to the 2017 fiscal year and were paid in 2018.

No meeting attendance fees were paid by Sartorius Stedim Biotech S.A. to the general management of the company in fiscal 2018. A Part of the Executive Board's remuneration has been recharged by Sartorius AG to Sartorius Stedim Biotech S.A. for an amount of €1,312 K.

12. Off-Balance Sheet Commitments (in thousands of €)

| Type of commitment | Comment | At Dec. 31, 2018 | At Dec. 31, 2017 |
|--|---------|---------------------|---------------------|
| Commitments given | | | |
| Guarantees for bilateral credit lines | | 0 | 0 |
| Guarantees for currency hedging contracts | | 0 | 0 |
| Commitments from renting / leasing | | 0 | 0 |
| Commitments received | | | |
| Contractual loan capacity from credit institutions | | 0 | 0 |

The commitments in connection with the lease are summarized in the following table:

| Leasing | < 1 year € in K | 1 - 5 years € in K | > 5 years € in K | Total | Buy-back value |
|----------------------------|--------------------|--------------------------|-------------------------|------------------------------|----------------------------|
| Tangible Assets | | | | | |
| Buildings and Improvements | 267 | 770 | 0 | 1,037 | 0 |
| Total | 267 | 770 | 0 | 1,037 | |
| Leasing | Historical value | Payments for the Year | Cumulatives Payments | Depreciation for the Year | Cumulative Depreciation |
| Tangible Assets | | | | | |
| Buildings and Improvements | 2,391 | 272 | 1,500 | 245 | 679 |
| Total | 2,391 | 272 | 1,500 | 245 | 679 |

The building will be operational from the 1st of January 2015.

13. Information on Related Parties (in thousands of €)

Affiliates are its parent company, Sartorius AG, and the companies owned by Sartorius Stedim Biotech S.A., and are Sartorius Stedim FMT S.A.S., Sartorius Stedim Bioprocess SARL, Sartorius Stedim Aseptics S.A. and Sartorius Stedim Biotech GmbH.

The company Sartorius Stedim Biotech S.A. is consolidated in the financial statements of Sartorius AG, Otto-Brenner-Strasse 20, 37079 Goettingen (Germany).

In the following, you will find the table of the main amounts with the related parties:

| Items | At Dec. 31, 2018 | At Dec. 31, 2017 |
|-------------------------|------------------|------------------|
| Investments | 127,977 | 127,977 |
| Trade receivables | 0 | 896 |
| Other receivables | 64,282 | 14,372 |
| Trade payables | 0 | 0 |
| Other liabilities | 79,322 | 49,960 |
| Income from investments | 49,007 | 48,409 |
| Other financial income | 336 | 15 |
| Finance expense | 939 | 1,727 |

In the following, you will find the table of subsidiaries and shareholdings:

| At Dec. 31, 2018 | Share capital | Reserves, share premium and retained earnings before appropriation | Ownership in % | Book value of shares held | | Loans outstanding and advances granted | Changes in deposits and pledges | Sales (ex-VAT) for the financial year | Net profit | Dividends received |
|---|---------------|--|----------------|---------------------------|--------|--|---------------------------------|---------------------------------------|------------|--------------------|
| | | | | Gross | Net | | | | | |
| | | | | | | | | | | |
| Sartorius Stedim Biotech GmbH | | | 100.00% | | | | | | | |
| (Euros) | 6,000 | 370,416 | | 79,949 | 79,949 | -46,029 | 0 | 559,070 | 121,563 | 45,000 |
| Sartorius Stedim FMT S.A.S. | | | 100.00% | | | | | | | |
| (Euros) | 42,940 | -7,597 | | 42,940 | 42,940 | -2,885 | 0 | 173,918 | 859 | 0 |
| Sartorius Stedim Bioprocess SARL | | | 99.99% | | | | | | | |
| (Dinars) | 5,950 | 779 | | | | | | 78,366 | 2,534 | 3,462 |
| (Euros) | | | | 3,132 | 3,132 | 5,832 | 0 | 25,173 | 814 | 1,007 |
| Sartorius Stedim RUS | | | 100.00% | | | | | | | |
| (Rubles) | 8,000 | 38,292 | | | | | | 321,051 | 32,738 | 0 |
| (Euros) | 100 | 480 | | 109 | 109 | 0 | 0 | 4,027 | 411 | 0 |
| Sartorius Stedim Aseptics S.A. | | | 100.00% | | | | | | | |
| (Euros) | 448 | 5,145 | | 1,848 | 1,848 | -6,999 | 0 | 13,515 | 3,971 | 3,000 |
| | | | | | | | | | | |
| At Dec. 31, 2017 | Share capital | Reserves, share premium and retained earnings before appropriation | Ownership in % | Book value of shares held | | Loans outstanding and advances granted | Changes in deposits and pledges | Sales (ex-VAT) for the financial year | Net profit | Dividends received |
| | | | | Gross | Net | | | | | |
| | | | | | | | | | | |
| Sartorius Stedim Biotech GmbH | | | 100.00% | | | | | | | |
| (Euros) | 6,000 | 299,853 | | 79,949 | 79,949 | -9,806 | 0 | 496,937 | 82,637 | 45,000 |
| Sartorius Stedim FMT S.A.S. | | | 100.00% | | | | | | | |
| (Euros) | 42,940 | -7,597 | | 42,940 | 42,940 | -2,885 | 0 | 148,938 | -7,061 | 0 |
| Sartorius Stedim Bioprocess SARL | | | 99.99% | | | | | | | |
| (Dinars) | 5,950 | 779 | | | | | | 66,983 | 3,342 | 0 |
| (Euros) | | | | 3,132 | 3,132 | 5,832 | 0 | 22,493 | 1,198 | 0 |
| Sartorius Stedim RUS | | | 100.00% | | | | | | | |
| (Rubles) | 8,000 | 5,554 | | | | | | 154,151 | -5,851 | 0 |
| (Euros) | 115 | 80 | | 109 | 109 | 0 | 0 | 2,221 | -84 | 0 |
| Sartorius Stedim Aseptics S.A. | | | 100.00% | | | | | | | |
| (Euros) | 448 | 5,145 | | 1,848 | 1,848 | -6,999 | 0 | 10,181 | 2,852 | 2,200 |

The previous list contains only information on transactions in Company shares received in accordance with the Article 19 MAR (Operations realized by

Executive Directors). Therefore, we are not aware of all transactions whose cumulative trade volumes have remained below the notification threshold of €20,000 per calendar year.

Statutory Auditors' Report on the Financial Statements

This is a translation into English of the statutory auditors' report on the financial statements of the Company issued in French and it is provided solely for the convenience of English speaking users.

This statutory auditors' report includes information required by European regulation and French law, such as information about the appointment of the statutory auditors or verification of the management report and other documents provided to shareholders.

This report should be read in conjunction with, and construed in accordance with, French law and professional auditing standards applicable in France.

For the year ended 31 December 2018

To the Shareholders of SARTORIUS STEDIM BIOTECH S.A.,

Opinion

In compliance with the assignment entrusted to us by your shareholders' meetings, we have audited the accompanying financial statements of SARTORIUS STEDIM BIOTECH S.A. for the year ended 31 December 2018.

In our opinion, the financial statements give a true and fair view of the assets and liabilities and of the financial position of the Company as at 31 December 2018 and of the results of its operations for the year then ended in accordance with French accounting principles.

The audit opinion expressed above is consistent with our report to the Audit Committee.

Basis for Opinion

Audit Framework

We conducted our audit in accordance with professional standards applicable in France. We believe that the audit evidence we have obtained is sufficient and appropriate to provide a basis for our opinion.

Our responsibilities under those standards are further described in the Statutory Auditors' Responsibilities for the Audit of the Financial Statements section of our report.

Independence

We conducted our audit engagement in compliance with independence rules applicable to us, for the period from 1 January 2018 to the date of our report and specifically we did not provide any prohibited non-audit services referred to in Article 5(1) of Regulation (EU) No 537/2014 or in the French Code of ethics (code de déontologie) for statutory auditors

Justification of Assessments – Key Audit Matters

In accordance with the requirements of Articles L.823-9 and R.823-7 of the French Commercial Code (code de commerce) relating to the justification of our assessments, we inform you of the key audit matters relating to risks of material misstatement that, in our professional judgment, concerned the appropriateness of the accounting principles applied, the reasonableness of the significant estimates used and the overall presentation of the accounts.

These assessments were addressed in the context of our audit of the financial statements as a whole, and in forming our opinion thereon, and we do not provide a separate opinion on specific items of the financial statements.

Specific verifications

We have also performed, in accordance with professional standards applicable in France, the specific verifications required by French law and regulations.

Information given in the management report and in the other documents with respect to the financial position and the financial statements provided to the Shareholders

We have no matters to report as to the fair presentation and the consistency with the financial statements of the information given in the management report of the Board of Directors and in the other documents with respect to the financial position and the financial statements provided to Shareholders.

We attest the fair presentation and the consistency with the financial statements of the information relating to payment terms, required under Article

D.441-4 of the French Commercial Code (Code de commerce).

Report on corporate governance

We attest that the Board of Directors' report on corporate governance sets out the information required by Articles L.225-37-3 and L.225-37-4 of the French Commercial Code.

Concerning the information given in accordance with the requirements of Article L.225-37-3 of the French Commercial Code (code de commerce) relating to remunerations and benefits received by the directors and any other commitments made in their favour, we have verified its consistency with the financial statements, or with the underlying information used to prepare these financial statements and, where applicable, with the information obtained by your company from controlling and controlled companies. Based on this work, we attest the accuracy and fair presentation of this information.

Other informations

In accordance with French law, we have verified that the required information concerning the purchase of investments and controlling interests and the identity of the shareholders and holders of the voting rights has been properly disclosed in the management report.

Report on Other Legal and Regulatory Requirements

Appointment of the Statutory Auditors

We were appointed as statutory auditors of Sartorus Stedim Biotech S.A. by the annual general meeting held on 7 April 2015 for KPMG S.A. and on 19 May 2006 for Deloitte & Associés.

As at 31 December 2018, KPMG S.A. was in its 4th year of the audit mandate without interruption and Deloitte & Associés was in its 13th year.

Responsibilities of Management and Those Charged with Governance for the Financial Statements

Management is responsible for the preparation and fair presentation of the financial statements in accordance with French accounting principles and for

such internal control as management determines is necessary to enable the preparation of financial statements that are free from material misstatement, whether due to fraud or error.

In preparing the financial statements, management is responsible for assessing the Company's ability to continue as a going concern, disclosing, as applicable, matters related to going concern and using the going concern basis of accounting unless it is expected to liquidate the Company or to cease operations.

The Audit Committee is responsible for monitoring the financial reporting process and the effectiveness of internal control and risks management systems and where applicable, its internal audit, regarding the accounting and financial reporting procedures.

The financial statements were approved by the Board of Directors.

Statutory Auditors' Responsibilities for the Audit of the Financial Statements

Objectives and audit approach

Our role is to issue a report on the financial statements. Our objective is to obtain reasonable assurance about whether the financial statements as a whole are free from material misstatement. Reasonable assurance is a high level of assurance, but is not a guarantee that an audit conducted in accordance with professional standards will always detect a material misstatement when it exists. Misstatements can arise from fraud or error and are considered material if, individually or in the aggregate, they could reasonably be expected to influence the economic decisions of users taken on the basis of these financial statements.

As specified in Article L.823-10-1 of the French Commercial Code (code de commerce), our statutory audit does not include assurance on the viability of the Company or the quality of management of the affairs of the Company.

As part of an audit conducted in accordance with professional standards applicable in France, the statutory auditor exercises professional judgment throughout the audit and furthermore:

- Identifies and assesses the risks of material misstatement of the financial statements, whether due to fraud or error, designs and performs audit procedures responsive to those risks, and obtains audit evidence considered to be sufficient and appropriate to provide a basis for his opinion. The risk of not detecting a material misstatement resulting from fraud is higher than for one resulting from error, as fraud may involve collusion, forgery, intentional omissions, misrepresentations, or the override of internal control.
- Obtains an understanding of internal control relevant to the audit in order to design audit procedures that are appropriate in the circumstances, but not for the purpose of expressing an opinion on the effectiveness of the internal control.
- Evaluates the appropriateness of accounting policies used and the reasonableness of accounting estimates and related disclosures made by management in the financial statements.
- Assesses the appropriateness of management's use of the going concern basis of accounting and, based on the audit evidence obtained, whether a material uncertainty exists related to events or conditions that may cast significant doubt on the Company's ability to continue as a going concern. This assessment is based on the audit evidence obtained up to the date of his audit report. However, future events or conditions may cause the Company to cease to continue as a going concern. If the statutory auditor concludes that a material uncertainty exists, there is a requirement to draw attention in the audit report to the related disclosures in the financial statements or, if such disclosures are not provided or inadequate, to modify the opinion expressed therein.
- Evaluates the overall presentation of the financial statements and assesses whether these statements represent the underlying transactions and events in a manner that achieves fair presentation.

Report to the Audit Committee

We submit a report to the Audit Committee which includes in particular a description of the scope of the audit and the audit program implemented, as well as the results of our audit. We also report, if any,

significant deficiencies in internal control regarding the accounting and financial reporting procedures that we have identified.

Our report to the Audit Committee includes the risks of material misstatement that, in our professional judgment, were of most significance in the audit of the financial statements of the current period and which are therefore the key audit matters that we are required to describe in this report.

We also provide the Audit Committee with the declaration provided for in Article 6 of Regulation (EU) N° 537/2014, confirming our independence within the meaning of the rules applicable in France such as they are set in particular by Articles L.822 - 10 to L.822 - 14 of the French Commercial Code (code de commerce) and in the French Code of Ethics (code de déontologie) for statutory auditors. Where appropriate, we discuss with the Audit Committee [ou autre terminologie retenue par la société pour désigner le comité spécialisé visé à l'article L.823 - 19 du code de commerce] the risks that may reasonably be thought to bear on our independence, and the related safeguards.

Marseille, 8 February 2019

The Statutory Auditors

French original signed by

John Evans

Vincent Gros

Supplementary Information

06

Other Information of a Legal Nature

General Information on the Issuer

Corporate Name

The corporate name of the company is: "Sartorius Stedim Biotech".

In all legal deeds and documents issued by the company, this is always preceded or followed by the words "société anonyme" or the abbreviation "S.A." and a statement of the share capital (Company bylaws, Heading 1, Article 1).

Registered Office

The registered office is in Aubagne (13400), France, Z.I. Les Paluds, avenue de Jouques.
Phone number: +33 (0)4 42 84 56 00.

This office may be transferred to another location in the same "département" [French county or state] or an adjacent county or state by simple decision of the Board of Directors subject to ratification by the next Annual General Shareholders' Meeting and anywhere else in France by a decision taken by an Extraordinary General Shareholders' Meeting.

If the Board of Directors decides to transfer the registered office, it is authorized to revise the bylaws as a result (Company bylaws, Heading 1, Article 2).

Legal Form and Applicable Law

The company is a public limited liability company or joint stock company [société anonyme], subject to the French legislation particularly to the French Commercial Code.

Date of Incorporation – Duration

The company was incorporated on September 28, 1978, as a "société anonyme." The company's duration is for 99 years, effective upon registration in the French trade and commercial register ("registre du commerce et des sociétés"), unless subject to dissolution or extension provided by the present company bylaws (Heading 1, Article 3).

Corporate Purpose

In France and abroad, the company's purpose is:

- to purchase, develop, administrate and manage a portfolio of equity security, securities, voting rights and other social rights in all companies regardless of their activity and this, by all means including by way of setting up of new companies, contribution in kind of any types of social rights, subscription rights, mergers, purchases of other social rights or incorporation of companies;
- to manage, conduct and coordinate the activities of its subsidiaries and affiliates; when applicable, to provide to said companies all services of an administrative, financial, accounting and legal nature and any opinion and advise or to order any studies or researches that are necessary for their development or growth;
- and more generally, all financial, commercial, industrial, personal and real property operations linked, directly or indirectly, to the above-mentioned corporate purpose or to all other complementary, related or similar purposes, which may promote the development or accomplishment thereof (Company bylaws, Heading 1, Article 4).

Trade and Commercial Register – APE Code

The company is registered with the "registre du commerce et des sociétés" de Marseille, under the number RCS B 314 093 352. Its economic activity code (APE) is 6420Z (Holding company activity).

Inspection of Legal Documents at the Registered Office of the Company

The reference document may be viewed at the registered office of the company, on its website and on the website of the AMF. During the validity of the present Reference Document, the bylaws, the Statutory Auditors' reports and the financial statements of the last three fiscal years, although with reports, mails and other documents, historical financial information of the company and its subsidiaries of the last three fiscal year, evaluation and declarations made by an expert, when these documents are statutory and any other statutory document, can be found at the registered office.

Financial Year

The financial year, also referred to as fiscal year, covers a period of twelve months, beginning on January 1 and ending on December 31 of each year (Company bylaws, Heading 1, Article 6).

Share capital

As of 31 December 2018, the share capital of the Company amounts to €18,436,038, divided in 92,180,190 shares of €0.20 fully paid; 74.3% of which are held by Sartorius AG.

Specific Clauses in the Company Bylaws

Form of Shares

Shares may be in nominative or bearer form according to the shareholder's choice. These shares are entitled to be recorded in an account in accordance with French law (Company bylaws, Heading 1, excerpt of Article 1).

Appropriation of Profits

The income statement that summarizes the income and expenses of the reporting year discloses by difference, after deduction of amortization, depreciation and provisions, the profit for said reporting year. At least 5% must be deducted from the annual profit reduced, where appropriate, by prior losses, to set up the legal reserve. This deduction ceases to be obligatory when the legal reserve amounts to one tenth of the share capital. This obligatory deduction resumes when, for whatever reason, the legal reserve falls below this one tenth. The distributable profit comprises the profit for the reporting year less prior losses and amounts transferred to reserves, pursuant to French laws and the company bylaws, and increased by profit brought forward. This profit is distributed among all shareholders in proportion to the number of shares each one holds. The Annual General Shareholders' Meeting may decide to distribute amounts taken from reserves available to it by expressly indicating the reserve from which the transfers are made. However, dividends are disbursed by way of priority from the annual profit for the reporting year. Except for a reduction in capital, no distribution may be made to shareholders when the equity falls below, or would consequently fall below, the amount of the capital together with the reserves that French laws or the

company bylaws do not permit to distribute. Revaluation surplus is not distributable. It may be incorporated in full or part into the company's capital. However, after transferring the amounts to the reserves, pursuant to French law, the Annual General Shareholders' Meeting may transfer any amount it considers necessary to all available reserves, ordinary or extraordinary reserves, or carry it forward.

Shareholders' Meetings

Convening

Annual (or Ordinary) General Shareholders' Meetings are those convened to take all decisions that do not result in a revision of the bylaws. Extraordinary General Shareholders' Meetings are those called to decide or authorize direct or indirect revisions to the bylaws. Special Meetings bring together the holders of a specific class of share to consider revisions to the rights of this class of share. Decisions made at the General Meetings are binding for all shareholders, even those who are absent, dissenting or legally incapable or incapacitated (Company bylaws, Heading 3, Article 13). General Meetings are convened by the Board of Directors or, by default, the independent auditors or a person thus empowered. General Meetings are held at the registered office or any other place stated in the notice of convocation (Company bylaws, Heading 3, excerpt of Article 14). The forms and timescale of the notice of convocation are governed by French laws.

Agenda

The notices and letters of call shall indicate the indications required by the law, particularly the agenda, the company electronic address where written questions of Shareholders may be sent and, eventually the mention of the obligation to collect the opinion or the prior approval of the mass of securities Shareholders giving access to the share capital.

The meeting may only deliberate on the matters placed on the agenda. It may, however, remove one or more directors at any time.

One or more shareholders representing the percentage of share capital required by law may, under the conditions and time limits set forth by law, require the inclusion on the agenda of draft resolutions.

In accordance to the Articles R 225 - 71 to R 225 - 74 of the Commercial Code, requests made by the Shareholders to register draft resolutions on the agenda and written questions are sent to the Headquarters by

registered letter with recorded delivery beginning on the publication of the Meeting announcement and until 25 days before the General Meeting, or in a delay of 20 days beginning on the publication of the Meeting announcement, when this one is published more than 45 days before the General Meeting (date of reception of the request by the company will be taken into account).

The request of a new item on the agenda must be motivated. The request to register draft resolutions is provided with the text of draft resolutions, which may have a short explanation of reasons. These requests are subject to justification of possession or representation of required Share capital, in accordance to regulatory rules (Company bylaws, Heading 3, Article 14, excerpt of point 2).

If the meeting has been unable to make a valid decision due to a lack of the required quorum, the second meeting and, where appropriate, the second meeting adjourned are called at least ten days in advance in the same form as the first meeting (Company bylaws, Heading 3, Article 14, excerpt of point 1).

Admission to Meetings – Powers

Every shareholder has the right to attend General Meetings and to participate in the discussions, in person or by proxy, regardless of the number of shares held, on simple proof of identity and the ownership of shares. The right to participate in a General Meeting is subject to the condition that the shares must be recorded, in the name of the shareholder or the shareholder's appointed broker, either in the nominative share accounts held by the company or in the bearer share accounts held by the authorized broker, by zero hours, Paris time, on the second working day prior to the meeting. The recording or registration of the shares in the bearer share accounts held by the authorized broker must be confirmed by a share certificate provided by the broker. This share certificate must be attached to the postal voting form, the proxy form or the application for an admission pass, issued in the name of the shareholder or on behalf of the shareholder represented by the appointed broker. A certificate must also be supplied to shareholders who wish to attend the General Meeting in person but who have not received an admission pass by zero hours, Paris time, on the second working day prior to the meeting.

A Shareholder may be represented by another Shareholder, his or her spouse or by the partner with who he or she signed a Civil Partnership. Furthermore, he or she may be represented by any other moral or

physical person of his choice in accordance to the Articles L. 225-106 to -106-3 of the Commercial Code; in that aim, the representative must present valid proof of proxy.

The legal representatives of shareholders who are legally incapable or incapacitated and individuals representing corporate shareholders take part in meetings, whether or not they are shareholders (Company bylaws, Heading 3, Article 14, point 3).

All Shareholders may also have a postal voting, using a registration form and sent to the company according to the law and regulations; to be acceptable this registration must be received by the company three days before the date of the Meeting.

In case of remote voting using an electronic vote, or a proxy vote given by electronic signature, this vote is made according to the conditions of the current regulations (Company bylaws, Heading 3, Article 14, point 4).

All legal documents relative to legal information for shareholders are made available to them at the registered office of the company.

Board of Directors

(Company bylaws, Heading 3, Point 3, Articles 6 to 9.)

1. Subject to legal exemptions, the Company is directed by a Board of Directors composed of a minimum of three members and a maximum of eighteen.

The composition of the Board of Directors is made with a balance number of men and women.

2. During the duration of the company's existence, directors shall be appointed or renewed in office by the ordinary general meeting. However, in case of merger, directors may be appointed by the extraordinary general meeting deciding on the transaction.

3. Each director must, during his entire term of office, own at least one share.

4. Directors have a term of office of three years.

Directors' duties shall cease at the end of the ordinary general meeting deciding on the accounts of the financial year elapsed, held in the year when the term of office of the director concerned expires.

Directors may be renewed in office. They may be removed from office at any time by the ordinary general meeting.

5. No person may be appointed director if, having reached the age of 75, his appointment would result in more than one third of the members of the board of directors exceeding that age. If that proportion is exceeded, the oldest director shall automatically be deemed to have resigned at the end of the ordinary general meeting approving the accounts of the financial year when exceeded.

6. Directors may be individuals or legal entities. Directors who are legal entities are required, upon their appointment, to appoint a permanent representative who is subject to the same conditions and obligations and who incurs the same liability as though personally a director, without prejudice to the several liability of the legal entity represented.

When the legal entity who is a director terminates the mandate given to its permanent representative, it shall promptly notify the Company, by registered letter, of its decision as well as the identity of its new permanent representative. The same applies in the event of death or resignation of the permanent representative.

7. If one or more directors' seats become vacant between two general meetings due to death or resignation, the board of directors may proceed to make appointments on an interim basis so as to fill the seats on the Board. These appointments must be made within three months of the vacancy, when the number of directors has fallen below the minimum under the articles of association but without falling below the statutory minimum.

Interim appointments made in this manner by the Board are subject to ratification by the next ordinary general meeting. Failing ratification, the decisions taken or the acts accomplished shall nonetheless remain valid.

When the number of directors falls below the statutory minimum, the directors remaining in office are required to immediately call an ordinary meeting so as to fill the vacant seats on the Board.

A director appointed in replacement of another shall only remain in office for the remaining term of office of his predecessor.

8. Directors who are individuals cannot concomitantly hold more than three seats on the board of directors or supervisory boards of sociétés anonymes having their registered office in metropolitan France, subject to the exceptions provided by law.

9. A Company employee may not be appointed a director unless his employment agreement corresponds to effective employment. He shall not lose the benefit of his employment agreement. The number of directors bound to the Company by an employment agreement may not exceed one third of the directors in office.

Organization and management of the Board of Directors

1. The Board of Directors elects a Chairman from among its members who are individuals and determines his remuneration. It sets the duration of the Chairman's term of office, which may not exceed his office as director.

2. No person may be appointed Chairman of the Board of Directors if over the age of 75. If the Chairman in office exceeds that age, he shall be deemed to have automatically resigned.

3. The Chairman represents the Board of Directors. He organizes and directs its work, and reports on it to the general meeting. He ensures the proper operation of the Company's decision-making bodies and ensures, in particular, that the directors are themselves in a position to fulfill their duties.

4. In case of absence or impediment affecting the Chairman, the Board of Directors appoints an acting Chairman of the meeting.

5. The Board of Directors appoints a secretary who may be chosen, either from among the directors or outside them. The secretary shall be replaced by simple decision of the Board.

Meetings and decisions of the Board

1. The Board of Directors meets, upon the call of its Chairman, as often as required by the interest of the Company. However, directors representing at least one third of the members of the Board of Directors may, by precisely indicating the meeting's agenda, call a Board if it has not met within the last two months.

The CEO, if not chairing the Board of Directors, may request the Chairman to call a Board meeting with a specified agenda.

2. The meeting shall take place at the registered office or in any other location indicated in the notice of call. The call to meeting, indicating the agenda, should be sent at least 7 days beforehand by letter, telegram, telex or fax. The call may be verbal and the meeting may be held immediately if all of the directors are in agreement.

3. For the Board of Directors to validly deliberate, at least one half of the directors are required to be present or represented.

The Board's decisions are taken at a majority of the members present or represented.

The acting Chairman has a casting vote.

4. An attendance sheet shall be held and signed by directors participating in the Board meeting.

5. The internal regulations established by the Board of Directors may provide that directors participating in a Board meeting by videoconference in accordance with the applicable regulations are deemed present for the purposes of calculating quorum and majority.

This provision shall not apply for the adoption of the following decisions:

- appointment, remuneration, removal of the Chairman, CEO and Executive Vice Presidents;
- closing of annual accounts, consolidated accounts and preparation of management report and report on the management of the group.

6. The Board of Directors' deliberations are recorded in minutes held in accordance with the applicable laws. The minutes are signed by the acting Chairman and by one or two directors.

Copies or excerpts of the minutes of the Board of Directors' deliberations shall be validly certified by the Chairman or by the CEO.

Powers of the Board of Directors

1. The Board of Directors determines the Company's business guidelines and ensures that they are implemented. Subject to the powers expressly granted by law to shareholders' meetings and within the limit of its corporate objects, it deals with any matter relating to the proper running of the Company and by its deliberations governs the affairs of the company.

In its dealings with third parties, the Company is bound even by acts of the Board of Directors that are outside its corporate purpose, unless it can prove that the third party knew that that act was ultra vires or could not reasonably have been unaware thereof in view of the circumstances, it being specified that mere publication of the articles of association does not suffice to establish proof thereof.

2. The Board of Directors shall carry out any controls and verifications it deems appropriate.

Each director shall receive the information necessary to the performance of his duties and may obtain all documents he considers useful from the General Management.

3. The Board of Directors may give all delegations of authority to the representatives of its choice within the limit of its authority under the law and under these articles of association.

The Board may decide on the creation of review committees in charge of studying the issues that the Board or its Chairman submits to it.

General Management (Company bylaws, Heading 3, Article 10)

Mode of operation

In accordance with Article L. 225-51-1 of the Commercial Code, the Company's General Management is ensured, under his responsibility, either by the Chairman of the Board of Directors or by any other individual appointed by the Board of Directors with the title of CEO.

The choice between these two modes of operation of General Management is made by the Board of Directors. The Board's decision concerning the choice of mode of operation of General Management is taken by majority vote of the directors present or represented. Shareholders and third parties are informed of the choice made by the Board of Directors under the conditions set forth by the applicable regulations.

The Board of Directors may modify the option chosen at any time.

A change in the mode of operation of General Management shall not entail any modification of the articles of association.

Depending on the mode of exercise chosen by the Board of Directors, the Chairman or a CEO shall ensure, under his responsibility, the General Management of the Company.

The CEO is appointed by the Board of Directors, which sets the duration of his term of office, determines his remuneration and, as applicable, the restrictions on his powers.

For the performance of his duties, the CEO must be under the age of 75. When this age limit is exceeded during the course of his term of office, the CEO shall be deemed to have automatically resigned and a new CEO shall be appointed.

The CEO may be removed from office at any time by the Board of Directors. Removal of a CEO who is not also the chairman may give rise to damages if decided without valid cause.

Powers of the CEO

The CEO is vested with the broadest powers to act in all circumstances in the name of the Company. The CEO shall exercise these powers within the limit of the corporate objects, and subject to the powers expressly granted by law to shareholders' meetings and to the Board of Directors.

The CEO represents the Company in its dealings with third parties. The Company is bound even by those acts of the CEO that are outside its corporate objects, unless it can prove that the third party knew that that act was ultra vires or could not reasonably have been unaware thereof in view of the circumstances, it being specified that mere publication of the articles of association does not suffice to establish proof thereof.

Executive Vice Presidents

Upon the motion of the CEO, whether this position is filled by the Chairman of the Board of Directors or by another person, the Board of Directors may name one or more individuals with responsibility for assisting the CEO with the title of Executive Vice Presidents.

The maximum number of Executive Vice Presidents may not exceed five.

In agreement with the CEO, the Board of Directors shall determine the scope and the extent of the powers granted to the Executive Vice Presidents and set their remuneration.

As regards third parties, the Executive Vice Presidents or the Executive Vice Presidents have the same powers as the CEO.

Upon the cessation of his duties or in case of impediment affecting the CEO, the Executive Vice Presidents shall retain, unless otherwise decided by the Board of Directors, their office and authority until the appointment of a new CEO.

The CEO may be removed from office at any time by the Board of Directors. Removal of a CEO who is not

also the chairman may give rise to damages if decided without valid cause.

Conditions for the Exercise of Voting Rights – Majority Quorum (Company bylaws, Heading 3, Article 15)

At Annual and Extraordinary General Meetings, the quorum is calculated on the basis of the shares comprising the share capital and, in Special Meetings, on the basis of all the shares of the class concerned, net of shares not entitled to voting rights by virtue of the law.

In the event of postal voting, only the forms received by the company prior to the meeting will be considered when calculating the quorum, under the conditions and timeframe set by the decree.

The right to vote conferred to shares is proportional to the capital they represent. With an equal par value, every share in capital or income right carries the right to one vote.

In the event that the shares are pledged, the voting right is exercised by the holder of the securities. The issuing company may not validly vote with shares subscribed, acquired or taken in pledge by it; these shares are not taken into account to calculate the quorum.

The voting takes place and the votes are cast by show of hands, or by those sitting and standing, or by roll call, as decided by the officers of the meeting.

Further Information on Voting Rights

There is no limit in the bylaws on voting rights.

A double voting right is conferred to the holders of registered shares that are fully paid up and that have been registered in the name of the same holder for at least four years.

In the event of conversion to bearer form, the converted share immediately forfeits its double voting right. In the event of a capital increase by incorporation of reserves, profits or share premium, this double voting right applies to new shares issued and allocated free of charge to a shareholder on the basis of existing shares that already carry this right (Heading 2, Article 3, of the company bylaws). This revision to the bylaws was unanimously passed by the General Shareholders' Meeting in an extra-ordinary session on August 24, 1994. It may be cancelled by a General Shareholders' Meeting convened in an

extraordinary session and after ratification by a Special Meeting of the beneficiary shareholders.

As of December 31, 2018, there were 68,844,424 shares with a double voting right out of a total of 92,180,190 shares. Thus, the total voting rights are 161,024,614.

The Annual General Shareholders' Meeting is held at least once a year, within six months of the year end, to consider the financial statements of that year, subject to an extension of this timeframe by a legal decision. The Annual General Shareholders' Meeting may only validly deliberate, upon the first convocation, if the shareholders present – represented or voting by post – hold at least one quarter of the shares with a right to vote. No quorum is required upon the second convocation. The meeting decides on the basis of the majority of votes held by shareholders present or represented, including shareholders voting by post (Company bylaws, Heading 3, Article 16).

Shareholders' agreement

None

Crossing Legal Thresholds

Any shareholder whose shareholdings cross the legal thresholds defined by French law, either upwards or downwards, must declare said crossing by notification of the Autorité des Marchés Financiers, pursuant to the law in force. The bylaws of the company do not provide for any additional threshold declarations.

Identification of Shareholders

Within the legal and regulatory framework, the company is authorized to seek the identity of bearer shareholders.

Payment of Dividends

The Annual General Shareholders' Meeting has the power to give every shareholder, for all or part of a dividend payable, the option of receiving this dividend in shares, as provided by French law, or in cash.

The terms of the payment of the dividend in cash are set by the General Meeting or, by default, the Board of Directors. Cash dividends must be paid within a maximum of nine months after the end of the reporting year, unless this timeframe is extended by legal authorization. However, this profit may be distributed as an interim dividend prior to the approval of the annual financial statements when a balance sheet prepared during or at the end of a financial year and certified by the independent auditors discloses that the company has realized a profit since the close of the previous financial year, after recognition of the necessary amortization, depreciation and provisions, as well as after deduction, where relevant, of prior losses and amounts to be transferred to the reserves, as required by French laws or the company bylaws. These interim dividends may not exceed the profit thus defined. No reimbursement of dividends may be required from shareholders unless the distribution was made in violation of legal provisions and the company determines that the beneficiaries were aware of the illegality of this distribution at the time it occurred or could not ignore this nature of the dividends. Where this occurs, the shares in reimbursement are time-barred three years after the payment of these dividends. Dividends not collected within five years of their payment are time-barred (Company bylaws, Heading 3, Article 22).

Financial score

None

Other Information on the Assets, Financial Position and Results for the Group

Major Contracts

Several service agreements were entered into between entities of the divisions of the Sartorius Group and Sartorius Stedim Biotech Group, in order to enable the entities from both divisions to benefit from certain general administrative services under the same terms.

Among these service agreements, the service agreement with the highest volume and importance is in place between Sartorius Stedim Biotech GmbH and Sartorius Corporate Administration GmbH, a 100% subsidiary of Sartorius AG. Sartorius Corporate Administration GmbH provides general administrative services to Sartorius Stedim Biotech and the other entities of the Sartorius Group. Such services include, among others, accounting, treasury management, payroll accounting for human resources, IT systems and legal services. Sartorius Corporate Administration GmbH invoices its services on the basis of the internal and external costs incurred plus a margin of 3%. The services invoiced by Sartorius Corporate Administration GmbH to Sartorius Stedim Biotech GmbH in 2018 totalled million €53.8 against million €40.1 in 2017.

Apart from the above-mentioned service agreements, there are no other contracts with material obligations or commitments that have been concluded outside the ordinary course of the company's business or to which a member of the Sartorius Stedim Biotech Group is a party.

The strategy of the Sales and Marketing organization within the Sartorius Stedim Biotech Group towards customers is to create valuable long-term relationships. Therefore, for example, key account management endeavors to conclude long-term framework contracts with customers. As a total solution provider, Sartorius Stedim Biotech strives to use such contracts to cover the entire product portfolio of Sartorius Stedim Biotech that fits into the validated processes of the customer.

Registered Trademarks and Trademark Applications

| Name | EU | Germany | France | International registration in the countries designated | USA | Australia | Brazil | Mexico | UK | Canada |
|--------------------------------|---|--|--|--|--|--|---|--|--|---|
| SARTORIUS STEDIM BIOTECH | 13/08/2007 No. 006228019 13/08/2027 | | | 16/11/2007 No. 962279 16/11/2027 + AU CH KR RU SG TR VN | 17/08/2007 No. 3709002 10/11/2019 | | 14/01/2008 9 Trademarks 3 Trademark Applications | | | 09/11/2007 No. TMA844101 21/02/2028 |
| BIOSTAT | 23/10/2014 No. 013398722 23/10/2024 | 04/10/1968 No. 873661 31/10/2028 | | 26/06/1985 No. 494574 26/06/2025 + AT BX CH DE ES FR IT PT | 22/07/1988 No. 1572999 26/12/2019 | | 16/12/2014 4 Trademark Applications | | 16/07/1988 No. 1246230 16/07/2026 | |
| HYDROSART | 12/11/2001 No. 002458461 12/11/2021 | 07/04/1983 No. 1065357 07/04/2023 | | | 10/12/2001 No. 2677224 21/01/2023 | | | | | 28/11/2001 No. 609610 06/05/2019 |
| MAXICAPS | 04/10/1999 No. 001330885 04/10/2019 | | | | 15/11/1999 No. 2450203 08/05/2021 | | | | | |
| MIDICAPS | 15/02/2005 No. 004289724 15/02/2025 | | | | 16/02/2005 No. 3195052 02/01/2027 | | | | | |
| MINISART | | 09/08/1978 No. 980370 09/08/2028 | 26/10/1988 No. 1495753 26/10/2028 | | 07/02/1979 No. 1144895 30/12/2020 | | | | 18/01/1979 No. 1107904 09/08/2019 18/01/1979 No. 1107903 18/01/2020 | |
| SARTOCHECK | | 13/06/1979 No. 987883 13/06/2019 | 17/10/1989 No. 1555685 17/10/2019 | | 05/12/1979 No. 1200237 06/07/2022 | | 18/11/2014 No. 908615248 30/05/2027 | | 20/12/1986 No. 1125952 20/12/2020 | |
| SARTOCON | | 06/06/1979 No. 988000 06/06/2019 | 17/10/1989 No. 1555684 17/10/2019 | | 15/06/1982 No. 1197792 15/06/2022 | | | | 20/12/1986 No. 1125951 20/12/2020 | |
| VIROSART | 02/11/2004 No. 004103701 02/11/2024 | 28/07/2004 No. 30443764 31/07/2024 | | | 08/02/2016 No. 86900738 24/01/2027 | | | | | |
| SARTOFLOW | | 03/06/1983 No. 1057870 30/06/2023 | | 06/03/1985 No. 494396 06/03/2025 + AT BX CH DE DZ EG ES FR HU IT KP LI MA MC PT RO RS RU SD VN | 08/08/2007 No. 3689721 09/29/2019 | | | | 25/10/1984 No. 1228900 25/10/2025 | |
| SARTOPORE | 10/01/2000 No. 001454461 10/01/2020 | | | | 15/02/2000 No. 2429825 20/02/2021 | | 18/11/2014 2 Trademark Applications | | | |
| FLEXBOY | 31/08/2005 No. 004614038 31/08/2025 | | 19/04/1993 No. 93465632 19/04/2023 | 24/01/1995 No. 630378 24/01/2025 + CH CN GB KR SE SG 27/02/2006 No. 879252 27/02/2026 + JP | 31/08/1993 No. 2041550 04/03/2027 | 31/01/1995 No. 651778 31/01/2025 | | 03/09/2003 No. 810249 03/09/2023 | 31/01/1995 No. 2009384 31/01/2025 | |
| FLEXEL | 20/02/1998 No. 000753202 20/02/2028 | | 02/09/1997 No. 97693975 02/09/2027 | | 27/02/1998 No. 2414947 26/12/2020 | | | 03/09/2003 No. 810250 03/09/2023 | | |
| PALLETANK | 01/07/1998 No. 000865865 01/07/2028 | | | 11/07/2016 No. 1314189 11/07/2026 + CH IN US | | | | | | |
| RAFT | 31/08/2005 No. 004614046 31/08/2025 | | | | | | | | | |
| EVAM | 15/10/1999 No. 001344266 15/10/2019 | | | | | | | | | |
| NUTRIKIT | | | 05/06/1989 No. 1535354 05/06/2019 | | | | | | | |
| NUTRIPOCHE | | | 05/06/1989 No. 1535352 05/06/2019 | | | | | | | |
| BIOSAFE | | | 01/02/1995 No. 95556118 01/02/2025 | 22/02/2001 No. 758706 22/02/2021 + DE DK GB CH | | | | | | |
| FLEXACT | 07/05/2009 No. 008285173 07/05/2019 | | | 16/10/2009 No. 1028463 16/10/2019 + AU CN JP KR US TR MX SG | | | 06/11/2009 4 Trademarks | | | 26/10/2009 No. 793270 18/11/2026 |
| FLEXSAFE | 22/04/2014 No. 012807996 22/04/2024 | | | 22/10/2014 No. 1226740 22/10/2024 + CN IN JP KR MX SG TR US | | | 21/10/2014 No. 9084706060 25/04/2027 | | | |

Apart from the trademarks mentioned above, the Sartorius Stedim Biotech Group is the owner | applicant of 335 different trademarks in various countries [the dates are indicated as day/month/year].

Registered Trademarks and Trademark Applications

| Name | Japan | Denmark | Finland | Ireland | Malaysia | Norway | Sweden | China | Argentina | India | Taiwan |
|--------------------------------|--|--|--|--|---|--|--|--|---|--|-----------------------------|
| SARTORIUS STEDIM BIOTECH | 08/11/2007 No. 5170560 03/10/2028 | | | | 28/11/2007 12 Trademarks | | | 14/01/2008 13 Trademarks | | 19/11/2007 13 Trademarks | 18/01/2008 13 Trademarks |
| BIOSTAT | 22/02/1988 No. 2021770 22/02/2028 27/08/1986 No. 1880889 27/08/2026 | 28/06/1985 No. 233586 29/08/2026 | 05/01/1988 No. 100350 05/01/2028 | 01/07/1985 No. 116688 30/06/2026 | 11/07/1985 No. 8502982 11/07/2022 | 27/05/1987 No. 128877 27/05/2027 | 31/03/1988 No. 209760 31/03/2028 | 26/04/2012 No. 10830519 14/03/2025 | 17/12/2014 3 Trademarks 1 Trademark Applications | 04/05/2012 No. 2326343 04/05/2022 | |
| HYDROSART | 21/11/2001 No. 4663672 18/04/2023 | | | | | | | | | | |
| MAXICAPS | 15/10/1999 No. 4535058 11/01/2022 | | | | | | | | | | |
| MIDICAPS | 25/02/2005 No. 4906540 04/11/2025 | | | | | | | | | | |
| MINISART | 09/02/1979 No. 1583197 26/04/2023 | | | | | | | | | | |
| SARTOCHECK | 29/09/1983 No. 1618759 29/09/2023 | | | | | | | | 14/11/2014 No. 3367508 16/10/2025 | | |
| SARTOCON | | | | | | | | | | | |
| VIROSART | 28/01/2005 No. 5040228 13/04/2027 | | | | | | | 24/11/2004 No. 4379959 21/06/2028 | | | |
| SARTOFLOW | | | | | | | | | | | |
| SARTOPORE | 02/02/2000 No. 4495393 03/08/2021 | | | | | | | | 12/11/2014 2 Trademarks | | |
| FLEXBOY | | | | | | | | | | | |
| FLEXEL | 02/03/1998 No. 4470133 27/04/2021 | | | | | | | | | | |
| PALLETANK | 28/02/2006 No. 5005301 24/11/2026 | | | | | | | | | | |
| RAFT | | | | | | | | | | | |
| EVAM | | | | | | | | | | | |
| NUTRIKIT | | | | | | | | | | | |
| NUTRIPOCHE | | | | | | | | | | | |
| BIOSAFE | | | | | | | | | | 10/08/2016 1 Trademark Application | |
| FLEXACT | | | | | | | | | 12/11/2014 4 Trademarks | 30/10/2009 4 Trademarks | |
| FLEXSAFE | | | | | | | | | 21/10/2014 No. 2857836 12/12/2026 | | |

Apart from the trademarks mentioned above, the Sartorius Stedim Biotech Group is the owner | applicant of 321 different trademarks in various countries [the dates are indicated as day/month/year].

Special Report of the Statutory Auditors on Related Party Agreements and Commitments

This is a translation into English of the statutory auditors' Special report on related party agreements and commitments of the Company issued in French and it is provided solely for the convenience of English speaking users.

This report should be read in conjunction with, and construed in accordance with, French law and professional auditing standards applicable in France.

Special report of the Statutory Auditors on related party agreements and commitments

For the year ended 31 December 2018

To the Shareholders of SARTORIUS STEDIM BIOTECH S.A.,

We are required to inform you, on the basis of the information provided to us, of the terms and conditions of those agreements and commitments indicated to us, or that we may have identified in the performance of our engagement. We are not required to comment as to whether they are beneficial or appropriate or to ascertain the existence of any such agreements and commitments. It is your responsibility, in accordance with article R. 225-31 of the French Commercial Code ('Code de Commerce'), to evaluate the benefits resulting from these agreements and commitments prior to their approval.

In addition, we are required, where applicable, to inform you in accordance with article R. 225-31 of the French Commercial Code concerning the implementation, during the year, of the agreements and commitments already approved by the General meeting of shareholders.

We performed those procedures which we considered necessary to comply with professional guidance issued by the French national auditing body ('Compagnie nationale des commissaires aux comptes') relating to this type of engagement. These procedures consisted in verifying that the information provided to us is consistent with the documentation from which it has been extracted.

Agreements and commitments submitted for approval by the general meeting of shareholders

Agreements and commitments authorized during the previous accounting period

We hereby inform you that we have not been advised of any agreements or commitments authorized and concluded during the previous accounting period to be submitted to the General meeting of shareholders for their approval in accordance with article L. 225-38 of the French Commercial Code.

Agreements and commitments from prior years not approved by the General meeting of the shareholders

We hereby inform you that the following agreements and commitments, authorized during the year ended 31 December 2017, which were mentioned in our special report on related party agreements and commitments for the year ended 31 December 2017 and which were not approved by the General meeting of shareholders approving the financial statements for the year ended 31 December 2017.

- General Assistance and Administrative Services Agreement
 - With the company, Sartorius AG (SAG) 74.3% shareholder of the company Sartorius Stedim Biotech S.A. (SSB S.A.)
 - Person concerned: Mr Joachim Kreuzburg (Chairman and Chief Executive Officer of SSB SA and Chief Executive Officer of the Executive Committee of SAG).
 - Nature and purpose: general assistance and administrative services agreement signed on 16 February 2017 with retrospective effect commencing 1 January 2015 for an indefinite duration. This agreement covers the recharging by SAG to SSB S.A. of a part of the remuneration of Mr Joachim Kreuzburg in respect of the services he performs and provides within the company.

- Details: the recharge of the said services of the person concerned is calculated using an allocation based on work performed and time spent by the executive for the benefit of SSB S.A.

The amounts invoiced (excluding tax) by SAG to SSB S.A. for the years ended 31 December 2017 and 2018 are detailed below:

- Year end 2018 : €. 674 216

- Year end 2017 : €. 759 996

- Regulated commitments concerning Mr Joachim Kreuzburg

- With the company Sartorius AG (SAG 74. 3% shareholder of Sartorius Stedim Biotech S.A. (SSB S.A.).

- Person concerned: Mr Joachim Kreuzburg (Chairman and Chief Executive Officer of SSB S.A. and Chief Executive Officer of the Executive Committee of SAG)

- Nature and purpose: commitments relating to early departure indemnities, a non-competition clause and supplementary retirement commitments were taken out by SAG for the benefit of Mr Joachim Kreuzburg

- Details: the details of these commitments are as follows:

Early departure indemnity cap

In the event of an early departure, at the initiative of the company, of Mr Joachim Kreuzburg from his executive function on the Executive Committee of SAG, the amount of the departure indemnity that falls due will be limited to a maximum amount corresponding to two years of remuneration.

Non-competition clause

For two years following the complete termination of his functions within the group SAG, Mr Joachim Kreuzburg will be obliged to comply with a non-competition clause compensated by an indemnity equal to one half of his last annual remuneration, if not waived or terminated.

Supplementary retirement commitments

Mr Joachim Kreuzburg benefits from a supplementary retirement scheme in compliance with German law requirements.

In accordance with the Sartorius group overall remuneration policy, these commitments will be recharged to Sartorius Stedim Biotech S.A. upon their occurrence for 20% of their amount.

Agreements and commitments already approved by the general meeting of shareholders

Agreements and commitments approved in previous financial years, which continued during the past year

We hereby inform you that we have not been advised of any agreements or commitments already approved by the General meeting of shareholders and continued during the previous financial year.

Marseille, 8 February 2019

The Statutory Auditors

French original signed by
KPMG Audit
A division of KPMG S.A.

John Evans

Deloitte & Associés

Vincent Gros

Resolutions Submitted to the Annual Combined Shareholders' Meeting on March 26, 2019

Dear Sir/Madam Shareholder,

We have summoned you in a combined Annual General Shareholders' Meeting in order to submit for your approval the eighteen resolutions whose purpose is described and commented below.

Please note that the description of the Company's activity required by the law is included in the management report related to 2018 financial year.

RESOLUTIONS SUBMITTED TO THE ORDINARY SHAREHOLDERS' MEETING

First resolution

(Approval of financial statements for the year ended 31 december 2018 and discharge to all Directors)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, after having considered the annual financial statements for the year ended 31 December 2018, the report of the Board of Directors and the Report of the statutory auditors concerning these financial statements, approved the financial statements for the year ended 31 December 2018, which disclosed a net profit of €49 521 306 as presented, and the transactions reflected in these financial statements or summarized in these reports.

As a result, the Shareholders' Meeting grants full and unreserved discharge to the Directors for the execution of their management duties for said reporting year.

The Shareholder's Meeting asserts that no overall expenses referred to in article 39, 4° of the general tax code were noted.

Second resolution

(Approval of the consolidated financial statements for the year ended 31 December 2018)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings has, after having considered the corporate consolidated accounts for

the year ended 31 December 2018, the report of the Board of Directors and the report of statutory auditors concerning these consolidated accounts, approved the consolidated financial statements for the year ended 31 December 2018, which disclosed a net profit of 209 959 400 as presented, and the transactions reflected in these financial statements or summarized in these reports.

Third resolution

(Assignment of the financial result for the financial year ended 31 December 2018)

The Annual Shareholders' meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, has decided to assign as follows, the income for the year ended 31 December 2018:

- Income of the year : €49 521 306
- Year-earlier profit carried forward: €34 345 883
- Distributable profit: €83 867 189
- Total amount of dividends to be disbursed to shareholders(*): €52 540 761
- Balance resulting from disbursement: €31 326 428

(*) The amount of dividends was calculated on the basis of the total number of shares as of December 31, 2018 (92,180,190 shares).

Each share of the company with a nominal value of €0,20 will entitle its holder to a payment of a net dividend valued at €0,57.

The dividend will be paid as from 2 April 2019.

The Shareholders' Meeting notes that for individual shareholders domiciled for tax purposes in France, dividends received are subject, pursuant to Article 200 A, 1 A 1° of the French General Tax Code, to a single flat-rate withholding tax of 12.8%, at the shareholder's option, such income may be taxed at the progressive income tax rate. In the latter case, dividends are eligible for the 40% allowance referred to in Articles 158 3 2° and 243 bis of the French General Tax Code. In both cases, when dividends are

paid, they are subject to a non-taxable withholding tax at the rate of 12.8% as an advance payment of income tax, which is deducted from the final tax due.

However, in accordance with the third paragraph of Article 117 quater of the French General Tax Code, individuals belonging to a tax household whose reference tax income is less than 50,000 euros for single, divorced or widowed taxpayers or €75,000 for taxpayers subject to joint taxation, may request exemption from this 12.8% withholding tax under the conditions provided for in Article 242 quater of the French General Tax Code.

In addition, for individual shareholders domiciled in France for tax purposes, social security contributions are applied in all cases to dividends paid of 17.2%.

The Shareholders' meeting acknowledges, pursuant to Article 243bis of the French general tax code, that the dividends paid for the last three financial years are the following:

| Exercise | Dividend ¹⁾ | Amount eligible for the 40% abatement | Amount not eligible for the 40% abatement | Dividend per shares ¹⁾ |
|----------|------------------------|---------------------------------------|---|-----------------------------------|
| 2017 | 42,402,887 | 42,402,887 | 0 | 0.46 € |
| 2016 | 38,713,209 | 38,713,209 | 0 | 0.42 € |
| 2015 | 30,734,476 | 30,734,476 | 0 | 2.00 € |

¹⁾ Prior deduction of social contribution on the dividend paid to physical person.

Fourth resolution

(Approval of regulated agreements covered by Article L.225-38 and subsequent of the French Commercial Code)

The Shareholder's Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, after having considered the report of the Board of Directors and the special report of the Statutory Auditors concerning regulated agreements as referred in articles L.225-38 and subsequent of the French Commercial Code, approves the conclusion of said reports and approves the regulated agreements which are mentioned in said special report.

The Shareholders' Meeting takes note, pursuant to the provisions of the Article L. 225-40 of the French commercial code, that the Shareholders who are parties to the regulated agreements mentioned in the special report cannot vote this resolution.

Fifth resolution

(Approval of regulated commitments covered by Article L.225 - 42-1 of the French Commercial Code related to Mr; Joachim Kreuzburg, Chief Executive Officer)

The Shareholder's Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, after having considered the report of the Board of Directors and the special report of the Statutory Auditors concerning regulated commitments as referred in articles L.225 - 42-1 of the French Commercial Code, takes notice of the conclusions of said report and approves the regulated commitments which are mentioned in such a special report, taken by Sartorius AG to the benefit of Mr. Joachim Kreuzburg, relating to a non-competition clause, an earlier departure severance and a supplementary pension scheme.

Shareholders who are parties to the regulated commitments mentioned in the special report cannot vote this resolution.

Sixth resolution

(Setting of the annual Directors' fees for the members of the Board of Directors)

The Shareholder's Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, after having considered the Board of Directors' report, approves the overall annual amount of the attendance fees allocated for the 2018 financial year and the followings years to come, until the Shareholders' Meeting decides otherwise, amounting to €268,800.

The Shareholders' Meeting grants full powers to the Board of Directors for allowing such attendance fees among its members, in whole or in part, and on such terms as it may determine.

Seventh resolution

(Approval of the fixe, variable and extraordinary components of the remuneration and the benefits of all kinds due or awarded to the Chairman and Chief Executive Officer for the 2018 financial year)

The Shareholder's Meeting, pursuant to the article L. 225-100 II of the French commercial code, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the corporate governance report of the Board of Directors, approves the fixe, variable and extraordinary components of the remuneration and the benefits of all kinds due or awarded to Mr. Joachim Kreuzburg, Chairman of the Board and Chief Executive Officer, for the 2018 financial year.

Eighth resolution

(Approval of the criteria and principles for determining, allocating and awarding components of the compensation and benefits of the Chairman and Chief Executive Officer for the 2019 financial year)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, after having considered the report of the Board of Directors established pursuant to Article L.225-37-2 of the French commercial code, approves the criteria and principles for determining, allocating and awarding components of the compensation and benefits of Mr. Joachim Kreuzburg, Chairman of the Board and Chief Executive Officer, for the 2019 financial year.

Ninth resolution

(Renewal of the term of M. Joachim Kreuzburg as Director)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the Board of Directors' report, takes note that the term of M. Joachim Kreuzburg, as a Director of the Company, expired at the end of the current general meeting, decides the renewal of this term of duty for a new three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

Tenth resolution

(Renewal of the term of M. Lothar Kappich as Director)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the Board of Directors' report, takes note that the term of M. Lothar Kappich, as a Director of the Company, expired at the end of the current general meeting, decides the renewal of this term of duty for a new three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

Eleventh resolution

(Renewal of the term of M. Henri Riey as Director)

The Shareholders' Meeting, in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the Board of Directors' report, takes note that the term of M. Henri Riey, as a Director of the Company, expired at the end of the current general meeting, decides to approve the renewal of this term of duty for a new three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

Twelfth resolution

(Nomination of Mrs. Pascale Boissel as new member of the Board of Directors of the Company)

The Shareholders' Meeting in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the Board of Directors' report, decides to appoint, as from today, Mrs. Pascale Boissel, born on October 15th, 1966, of French nationality, as a new member of the Board of Directors of the Company for a three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

Thirteenth resolution

(Nomination of Mr. René Fáber as new member of the Board of Directors of the Company)

The Shareholders' Meeting in accordance with the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, and after having considered the Board of Directors' report, decides to appoint, as from today, Mr. René Fáber, born on July 18th, 1975, of Slovakian nationality, as a new member of the Board of Directors of the Company for a three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

Fourteenth resolution

(Authorization granted to the Board of Directors to enable the Company to trade in its own shares)

The Shareholders' Meeting, having fulfilled the quorum and majority requirements applicable to Ordinary Shareholders' Meetings, having considered the report of the Board of Directors, in compliance with the provisions of articles L. 225-209 et seq. of the French Commercial Code, the directly applicable provisions of the European Commission regulation no. 2273/2003 of 22nd December 2003, the General regulation of the Autorité des marchés financiers (AMF – Financial Market Authority), and the market practices accepted by the AMF:

1) authorizes the Board of Directors, having the right to sub-delegate in compliance with applicable laws and regulations, to make the Company acquire, hold, or transfer, on one or more occasions, shares of the Company in connection with the implementation of a share buy back program subject to the provisions of Articles L. 225-209 et seq. of the French Commercial Code;

2) decides that the acquisition, sale or transfer of such shares may be effected by any means on the market or over-the-counter, including through the acquisition of blocks of shares; these means include the use of any derivative financial instrument traded on a regulated market or over-the-counter or the delivery of shares as a result of the issuance of securities giving access to the Company's capital through conversion, exchange, redemption, exercise of a warrant or in any other manner either directly or through an investment service provider; the maximum share of the capital acquired or transferred in blocks may reach the entire program; these transactions may be carried out at any time, including during periods of public offer on the

capital of the Company, in compliance with the regulations in force;

3) decides that the share buy back program will have, as objective to promote liquidity and stimulate the market price of the Company's shares under a liquidity contract in accordance with the AMAFI Code of Ethics recognized by the AMF;

4) decides that the terms and conditions of the share buy back program are the followings:

duration of the program: a maximum of 18 months, starting from the date of the present Shareholders' Meeting and expiring on the date when any Shareholders' Meeting of the Company adopts a new share buy back program or, alternatively, on September 26, 2020;

maximum redemption percentage: 0.10% of the share capital, ie 92 180 shares on the basis of 92 180 190 shares making up the share capital at the date of this Shareholders' Meeting; being specified that this limit applies to an amount of the share capital of the Company, which may be adjusted by the Board of Directors to take account of transactions affecting the share capital after the date of the present Shareholders' Meeting, the acquisitions made by the Company can not in any case cause it to hold, directly or indirectly through its subsidiaries, more than 10% of its share capital;

when the shares are acquired in order to encourage the liquidity of the Company's shares under the conditions defined by the AMF's general regulations, the number of shares taken into account for the calculation of this limit will correspond to the number shares purchased net of the number of shares resold during the term of the authorization;

maximum unit purchase price (excluding fees and commissions): 150 euros, ie a maximum theoretical amount allocated to the share buyback program of 13 827 000 euros on the basis of the maximum percentage of 0.10%, excluding trading costs the maximum theoretical amount will be adjusted by the Board of Directors to take into account transactions affecting the share capital after the date of this Shareholders' Meeting.

5) The dividends attached to the treasury shares of the Company shall be affected to the retained earnings account;

6) The Shareholders' Meeting grants to the Board of Directors, with right to sub-delegate in compliance with applicable laws and regulations, all necessary

powers to implement this authorization and in particular to establish the terms and conditions of the share buy-back program in compliance with applicable laws and with the present resolution, and notably to proceed, as the case may be, with any adjustment required by transactions on the share capital; to place any purchase order on the stock market; to enter any agreement, notably for the keeping of registers of sale and purchase of shares, to make any and all declarations to the AMF and any other organization, to carry out all formalities, and more generally, to take all appropriate measures;

7) This authorization invalids any prior authorization given on an identical subject matter.

Fifteenth resolution

(Proxy to carry out formalities)

The Shareholders' Meeting gives full authority to the bearer of an original, a copy or an extract of the minutes from the present Annual Shareholders' Meeting to accomplish each necessary procedure.

RESOLUTIONS SUBMITTED TO THE EXTRAORDINARY SHAREHOLDERS' MEETING

Sixteenth resolution

(Update of Bylaws; adoption of the new Bylaws)

The Shareholders' Meeting, having fulfilled the quorum and majority requirements applicable to Extraordinary Shareholders' Meetings, having considered the report of the Board of directors:

decides to update the Company's Bylaws in order to be compliant with the applicable law,

decides, therefore, to update the Company's bylaws as described in the report of the Board of Directors,

adopts, article by article, and as a whole, the new Company's Bylaws

Seventeenth resolution

(Update of Bylaws for determine the procedures for appointing directors and directors representing employees; update of the article 15 of Bylaws)

The Shareholders' Meeting, having fulfilled the quorum and majority requirements applicable to Extraordinary Shareholders' Meetings, having considered the report of the Board of directors and advice of the works committee:

takes note, that the company exceeds, at the end of two consecutive financial years, the thresholds provided in Article L. 225-27-1 I of the French Commercial Code relating to the obligation to appoint one or more director(s) representing employees

decides, in accordance with the provisions of articles L.225-27-1 et seq. of the French commercial Code, to update the Company's bylaws to determine the procedures of appointment of the representing employees

decides, consequently, to update the article 15 of the Bylaws, as updated in the sixteenth resolution above, as follows :

« Article 15 : Board of directors

[Unchanged]

15.9 Directors representing employees

15.9.1 In accordance with the applicable law, there shall be one director representing employees when the number of directors is equal to or less than 12. The director representing employees is:

- elected by the employees of the company and its direct or indirect subsidiaries which have their registered office located in France under the conditions provided in this article, or
- appointed by the trade union organisation that obtained the most votes during the first round of the elections mentioned in Articles L. 2122-1 et L. 2122-4 of the French Labour Code in the Company and its direct or indirect subsidiaries which have their registered office located on France; or
- appointed by the works council.

When the number of directors is more than 12, a second director representing employees is:

- elected by the employees of the company and its direct or indirect subsidiaries which have their registered office located in France, or
- appointed by the trade union organisation that obtained the most votes during the first round of the elections mentioned in Articles L2122-1 et L2122-4 of the French Labour Code in the Company and its direct or indirect subsidiaries, of which the registered offices are located in France, or
- appointed by the works council, or
- appointed by the European works committee.

The absence of the appointment of one or more directors representing employees in application of the applicable law and the present constitution shall not entail the invalidity of the deliberations of the board of directors.

15.9.2 – Directors representing employees are not included in the minimum number and maximum number of directors specified in Articles L.225-17 and L.225-18-1 of the French Commercial Code..

15.9.3 – Directors representing employees must have an employment contract with the Company or with one of its direct or indirect subsidiaries which have their registered office located in France predating their appointment by at least two years and relating to an actual employment.

15.9.4 – For each directors representing employees, voting procedures according to the applicable law, in particular the Article L.225-28 of the French commercial code.

All the company's employees who comply with the conditions provided by the applicable law shall be electors.

Should there be one directorship to fill, a majority vote with two rounds of voting must be held. In addition to the name of the candidate, each candidacy must include the name of a potential substitute. The candidate and the substitute must not be of the same gender.

Should there be two directorships to fill in an electing body, the election shall be by proportional representation based on the list according to the highest vote and without vote splitting. Each list must have four candidates and must be composed alternatively of a candidate of each gender.

In the event of a tied vote, the candidates with the earliest-dated employment contracts shall be declared elected.

Elections shall be held in such a way that a second round of voting may be held no later than eight days before the end of the term of office of the outgoing directors representing the employees.

At each election, the board of directors shall draw up the list of subsidiaries and fix the date of the elections that would meet the following deadlines:

- The date of the election is held at least [8] weeks before the election,
- The list of the electors is held, at least [6] weeks before the election,
- The submission of candidatures is held, at least [5] weeks before the election,
- The list of the candidates is held at least [four] weeks before the election,
- Sending the documents for the vote by post, at least three weeks before the election.

In case of absence of candidatures, the corresponding seat (s) remain vacant until the next election, which renew the directors representing employees office.

The vote is held electronically and/or in paper form.

In the case of a paper vote, the voting takes place on a single day, at the place of work and during working hours. However, employees can vote by correspondence:

- Staff members who are in a predictable absence situation on the election date,
- Staff members who, because of the nature or conditions of their work, are away from the voting place to which they are assigned,
- Staff members working on sites that do not have a voting place.

In the case of voting by electronic and/or in paper form, the procedures related to the organization and conduct of the election of the directors representing employees not specified by the laws or regulations in force or by the present constitution are adopted by the board of directors, or by delegation by its chairman, where appropriate, by implementing any group agreement that would have been entered into

with respect to the terms of such election, in the companies of the perimeter referred to above.

15.9.5 – Directors representing employees are elected for 3 years. The term of office of the director thus appointed shall end during the ordinary shareholders Meeting of the closing of the accounts, held the year of the end of the term of the office.

Directors representing employees shall be entitled for re-election.

15.9.6 – The termination of the employment contract shall end the office of the directors representing employees.

Directors representing employees may not be dismissed other than for fault in the performance of their office by order of the judge of the Tribunal de Grande Instance territorially competent, ruling by way of summary proceedings at the request of the majority of the members of the board of directors.

15.9.7 – In the event of vacancy of an office of a director representing employees due to death, resignation, dismissal, breach of employment contract or for any reason whatsoever, the vacant office shall be filled pursuant to Article L.225-34 of the French commercial code.»

Eighteenth resolution

(Proxy to carry out formalities)

The Shareholders' Meeting gives full authority to the bearer of an original, a copy or an extract of the minutes from the present Annual Shareholders' Meeting to accomplish each necessary procedure.

Report of the Board of Directors

BOARD OF DIRECTORS' REPORT ON RESOLUTIONS SUBMITTED TO THE ANNUAL GENERAL SHAREHOLDERS' MEETING ON 26 MARCH 2019.

Dear Shareholders,

We have convened you to an Ordinary and Extraordinary General Shareholders' Meeting to be held on 26 March 2019, at 13.30pm, at the Company's headquarters Z.I. Les Paluds, Avenue de Jouques, 13400 Aubagne, in order to deliberate on the eighteen resolutions whose purpose is described and commented below.

The purpose of this report is to present a general explanation of the draft resolutions proposed by the Board of Directors.

Please note that the description of the Company's activity required by the law is included in the management report and the registered document related to 2018 financial year.

For further information, the management report and the auditors reports will be presented at the Shareholders' meeting.

The documentation of the Shareholder's meeting, in particular the management report, the registered document and the auditors reports will be provided under the conditions set forth by the applicable law.

RESOLUTIONS SUBMITTED TO THE ORDINARY SHAREHOLDERS MEETING

Approval of the annual financial statements and the consolidated financial statements for the year ended 31 December 2018 (Resolutions n°1 and n°2)

In the first resolution, we submit to your approval the Sartorius Stedim Biotech SA's financial statements for the year ended on 31 December 2018 which resulted in a net profit of €49 521 306 and to discharge all directors.

Please note the absence of expenditures referred to in article 39.4 of the general tax code.

In the second resolution, we submit to your approval the consolidated financial statements for the year ended 31 December 2018 which resulted in a net profit of €209 959 400 .

The detailed financial statements are included in the Reference Document available on the Company's website www.sartorius-france.fr.

Allocation of the financial result for the year ended 31 December 2018 (Resolution n°3)

The net profit resulting from the 2018 financial statements amounts to €49 521 306 to add to the Year-earlier profit carried forward of €34 345 883. This would yield a distributable profit of €83 867 189.

We propose to allocate the net profit as follows:

– Total amount of dividends to be disbursed to shareholders €52 540 761

– The remaining amount of €31 326 428 to be carried forward to the next year now amounting €31 326 428.

It is proposed to set the 2018 net dividend to €0.57 per share.

The dividend will be paid as from 2 April 2019.

It is stated that the distributed amount of €0.57 per share will be eligible to an allowance of 40% applied to physical people residing in France, as referred in article 158.3-2 of the French general code tax

Pursuant to Article 243 bis of the French general tax code, it is also stated that distributed amounts under the three last financial years have amounted as follows:

| Exercise | Dividend | Amount eligible for the 40% abatement | Amount not eligible for the 40% abatement | Dividend per shares ¹⁾ |
|----------|------------|---------------------------------------|---|-----------------------------------|
| 2017 | 42,402,887 | 42,402,887 | 0 € | 0.46€ |
| 2016 | 38,713,209 | 38,713,209 | 0 € | 0.42€ |
| 2015 | 30,734,476 | 30,734,476 | 0 € | 2.00€ |

1) Prior deduction of social contribution on the dividend paid to physical person.

mentions their financial terms and the amounts invoiced in 2018.

The auditor's special report on the regulated agreements and commitments would be provided under the conditions set forth by the applicable law.

We draw your attention on the fact that shareholders having an interest in said regulated agreements shall not vote the corresponding resolutions.

Approval of the attendance fees (Resolution n°6)

We submit to your approval the overall annual amount of attendance fees allocated to the Board of Directors amounting to €268,800 for the financial year ending 31 December 2018 as well as for the following financial years and until an adverse decision.

Full powers would be granted to the Board of Directors for allowing such attendance fees among its members, in whole or in part, and on such terms as it may determine.

Approval of the principles and the elements of compensation granted to the Chief Executive Officer (Resolutions n°7 and 8)

We submit to your the approval of the following information:

- the fixed, variable and exceptional components of the total remuneration and benefits of all kinds due or granted to the Chief Executive Officer for the financial year ending 31 December 2018, as presented in the Reference Document section "Remuneration of the Executive and Non-executive Members of the Board".
- the remuneration policy of the Chief Executive Officer and, more precisely, the principles and criteria for determining, allocating and awarding the fixed, variable and exceptional components of the total remuneration and benefits of all kinds awarded to the Chairman and Chief Executive Officer, as presented in the corporate governance report included in the Reference Document (section 3). These principles and criteria were determined by the Board of Directors after an opinion of the Remuneration Committee. The amounts resulting from the application of these principles and criteria will be submitted to the approval of the Shareholders Meeting convened to approve the 2019 financial statements.

Expiry of the term of five Directors; renewal of term (Resolutions n°9 to 13)

The term of Directors of Mr. Joachim Kreuzburg, Mr. Lothar Kappich, Mr. Bernard Lemaître, Mr. Henri Riey and Mrs. Liliane de Lassus, will expire at the end of the ordinary part of the Shareholders Meeting as of 26 March 2019.

Following proposal of the Remunerations and Nominations Committee, we submit to your approval the renewal of the term of duty of Mr. Joachim Kreuzburg, Mr. Lothar Kappich and Mr. Henri Riey for a new three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021. As a consequence, we submit to your approval the non-renewal of the mandates of Directors of Mr. Lemaître and Mrs de Lassus.

In addition, following proposal of the Remunerations and Nominations Committee, we submit to your approval the appointment of Mrs. Pascale Boissel and Mr. René Fáber as new members of the Board of Directors of the Company, starting as of this day for a three-year term to expire at the end of the Annual Shareholders' Meeting of 2022 convened to approve the financial statements of the financial year ending 31 December 2021.

In the case of approval of the ninth to thirteenth resolutions, the Board of Directors will be composed of the 7 following Directors:

- Mr. Joachim Kreuzburg
- Mr. Lothar Kappich
- Mr. Henri Riey
- Mrs. Susan Dexter
- Mrs. Anne-Marie Graffin
- Mrs. Pascale Boissel
- Mr. René Fáber

Authorization for the Company to buy back its own shares (Resolutions 14)

The Shareholders' Meeting held on April 3rd, 2018, in its thirteenth resolution, had approved a share buyback program for 18 months, up to 0.10% of the share capital with a maximum purchase price of €150 per share. The objective of the buy back program would be to promote liquidity and stimulate the market price of the Company's shares under a liquidity agreement.

We propose to renew this share buyback program and to allow the Board of Directors' to buy back shares of the Company, for an eighteen month period as from the next Shareholders meeting, up to a maximum of 0.10% of the share capital.

The objective of the buy back program would be to promote liquidity and stimulate the market price of the Company's shares under a liquidity contract in accordance with the AMAFI Code of Ethics recognized by the AMF.

The authorization is granted within the following limits:

- 0.10% of the share capital;
- maximum purchase price: €150.00 per share;
- maximum budget: €13,827,000.

In accordance with the law, the transactions may be carried out at any time, including during the period of a public offer as the company's shares. It is important that the Company is able, where applicable, and even during a public offer, to buy back its own shares with a view to achieving the objectives of the buy back program.

Proxy to carry out formalities (Resolution 15)

The Shareholders' Meeting gives full authority to the bearer of an original, a copy or an extract of the minutes from the present Annual Shareholders' Meeting to accomplish each necessary formalities.

RESOLUTIONS SUBMITTED TO THE EXTRAORDINARY SHAREHOLDERS MEETING

Update of Bylaws; adoption of the new Bylaws (Resolution 16)

We hereby ask for an update of the Company's bylaws in order to be compliant with applicable law and reglementation.

Therefore, we hereby ask for an update of the Company's bylaws as follow:

- Re-numbering of Bylaws' articles;
- Restructuing of Bylaws' articles;
- Updating Article 4 "Registered Office" in order to specify that the registered office may be transferred to any other location in France by decision of the Board of Directors, subject to ratification by the Shareholders' Meeting;
- Updating of Article 11 "Transfer and conveyance of shares" in order to insert the information to be provide in case of crossing of thresholds;
- Updating of article 15 "Board of Directors" in order to remove the rules for minimum shareholding by the Directors;
- Updating of Articles 22 "Shareholders' meeting" and 27 "Dissolution – Liquidation" in a view of simplification and harmonization with the applicable law.

Therefore, we submit to your approval, article by article, and as a whole, the new Company's Bylaws.

Update of Bylaws for determine the procedures for appointing directors and directors representing employees; update of the article 15 of Bylaws (Resolution 17)

The Company exceeds, at the end of two consecutive financial years 2017 and 2018, the thresholds provided in Article L. 225-27-1 I of the French Commercial Code relating to the obligation to appoint one or more director(s) representing employees

Therefore in accordance with the provisions of articles L.225-27-1 et seq. of the French commercial Code, we hereby ask for an update of the Company's bylaws to determine the procedures of appointment of the

employees representative. Consequently, Article 15 of the Bylaws will be updated as follows :

« Article 15 : Board of directors

[Unchanged]

15.8 - Directors representing employees

15.8.1 - In accordance with the applicable law, there shall be one director representing employees when the number of directors is equal to or less than 12. The director representing employees is:

- elected by the employees of the company and its direct or indirect subsidiaries which have their registered office located in France under the conditions provided in this article, or
- appointed by the trade union organisation that obtained the most votes during the first round of the elections mentioned in Articles L. 2122-1 et L. 2122-4 of the French Labour Code in the Company and its direct or indirect subsidiaries which have their registered office located on France, or
- appointed by the works council.

When the number of directors is more than 12, a second director representing employees is:

- elected by the employees of the company and its direct or indirect subsidiaries which have their registered office located in France, or
- appointed by the trade union organisation that obtained the most votes during the first round of the elections mentioned in Articles L2122-1 et L2122-4 of the French Labour Code in the Company and its direct or indirect subsidiaries, of which the registered offices are located in France, or
- appointed by the works council, or
- appointed by the European works committee.

The absence of the appointment of one or more directors representing employees in application of the applicable law and the present constitution shall not entail the invalidity of the deliberations of the board of directors.

15.8.2 - Directors representing employees are not included in the minimum number and maximum number of directors specified in Articles L.225-17 and L.225-18-1 of the French Commercial Code..

15.8.3 - Directors representing employees must have an employment contract with the Company or with one of its direct or indirect subsidiaries which have their registered office located in France predating their appointment by at least two years and relating to an actual employment.

15.8.4 - For each directors representing employees, voting procedures according to the applicable law, in particular the Article L.225-28 of the French commercial code.

All the company's employees who comply with the conditions provided by the applicable law shall be electors.

Should there be one directorship to fill, a majority vote with two rounds of voting must be held. In addition to the name of the candidate, each candidacy must include the name of a potential substitute. The candidate and the substitute must not be of the same gender.

Should there be two directorships to fill in an electing body, the election shall be by proportional representation based on the list according to the highest vote and without vote splitting. Each list must have four candidates and must be composed alternatively of a candidate of each gender.

In the event of a tied vote, the candidates with the earliest-dated employment contracts shall be declared elected.

Elections shall be held in such a way that a second round of voting may be held no later than [eight] days before the end of the term of office of the outgoing directors representing the employees.

At each election, the board of directors shall draw up the list of subsidiaries and fix the date of the elections that would meet the following deadlines:

- The date of the election is held at least [8] weeks before the election,
- The list of the electors is held, at least [6] weeks before the election,
- The submission of candidatures is held, at least [5] weeks before the election,
- The list of the candidates is held at least [four] weeks before the election,
- Sending the documents for the vote by post, at least three weeks before the election.

In case of absence of candidatures, the corresponding seat (s) remain vacant until the next election, which renew the directors representing employees office.

The vote is held electronically and/or in paper form.

In the case of a paper vote, the voting takes place on a single day, at the place of work and during working hours. However, employees can vote by correspondence:

- Staff members who are in a predictable absence situation on the election date,
- Staff members who, because of the nature or conditions of their work, are away from the voting place to which they are assigned,
- Staff members working on sites that do not have a voting place.

In the case of voting by electronic and/or in paper form, the procedures related to the organization and conduct of the election of the directors representing employees not specified by the laws or regulations in force or by the present constitution are adopted by the board of directors, or by delegation by its chairman, where appropriate, by implementing any group agreement that would have been entered into with respect to the terms of such election, in the companies of the perimeter referred to above.]

15.8.5 – Directors representing employees are elected for 3 years. The term of office of the director thus appointed shall end during the ordinary shareholders' Meeting of the closing of the accounts, held the year of the end of the term of the office.

Directors representing employees shall be entitled for re-election.

15.8.6 – The termination of the employment contract shall end the office of the directors representing employees.

Directors representing employees may not be dismissed other than for fault in the performance of their office by order of the judge of the Tribunal de Grande Instance territorially competent, ruling by way of summary proceedings at the request of the majority of the members of the board of directors.

15.8.7 – In the event of vacancy of an office of a director representing employees due to death, resignation, dismissal, breach of employment contract or for any reason whatsoever, the vacant office shall be filled pursuant to Article L.225-34 of the French commercial code.»

Proxy to carry out formalities (Resolution 18)

The Shareholders' Meeting gives full authority to the bearer of an original, a copy or an extract of the minutes from the present Annual Shareholders' Meeting to accomplish each necessary procedure.

We hope that the different proposals made in this report will meet your approval and that you will agree to vote in favor of the corresponding resolutions.

The Board of Directors
represented by its Chairman
Mr Joachim Kreuzburg

Information on the Reference Document and the Annual Financial Report

Declaration of Responsibility for the Reference Document and the 2018 Annual Financial Report

I hereby certify, after having taken all reasonable measures to this effect, that the information contained in the present Reference Document is, to the best of my knowledge, in accordance with the facts and makes no omission likely to affect its import.

I certify, to the best of my knowledge, that the financial statements have been prepared in accordance with applicable accounting standards and give a fair view of the assets, liabilities and financial position and profit or loss of the company and all the activities included in the consolidation, and that the management report enclosed presents a fair review of the development and performance of the business and financial position of the company and of all the activities included in the consolidation as well as a description of the main risks and uncertainties to which they are exposed.

I have received a completion letter from the auditors stating that they have audited the information contained in this Reference Document about the financial position and financial statements and that they have read this document in its entirety.

February 18, 2019



Joachim Kreuzburg
Chairman of the Board and CEO

Table of Reconciliation

In order to facilitate understanding of the present document concerning the presentation of Sartorius Stedim Biotech S.A., the table below has, on the left,

the headings from Note 1 of European Regulation No. 809/2004 of April 29, 2004, of the European Commission and in the column on the right, the corresponding pages of the present document.

| Headings of Part 1 of European Regulation N°809/2004 of April 29, 2004 | | Pages |
|--|--|---------------------------|
| 1. | Persons responsible | |
| 1.1. | Persons responsible for information | 178 |
| 1.2. | Certification of persons responsible for registering document | 178 |
| 2. | Independent auditors | |
| 2.1. | Name and address of Independent Auditors of the issuer | 86-87 |
| 3. | Selected financial information | |
| 3.1. | Presentation of selected historical financial information for every year of the period covered by this financial information | 24-33 |
| 4. | Risk factors | 35-46 |
| 5. | Information on the issuer | |
| 5.1. | History and development of the company | 12 - 15 |
| 5.1.1. | Corporate name and commercial name of issuer | 154 |
| 5.1.2. | Place and registration number of issuer | 5, 154 |
| 5.1.3. | Date of establishment and life of issuer | 154 |
| 5.1.4. | Registered office and legal form of issuer, legislation governing its operation, country of origin, address | 154 |
| 5.2. | Investments | 28 |
| 5.2.1. | Principal investments (including their amounts) carried out | 32, 107 - 111 |
| 6. | Overview of operations | |
| 6.1. | Principal operations | 19, 24 - 26 |
| 6.2. | Principal markets | 22 - 23 |
| 6.3. | Dependence on patents, licenses and contracts | 39, 162 - 163 |
| 6.4. | Competition | 22-23 |
| 7. | Organigram (organizational charts) | |
| 7.1. | Description of group | 18, 100- 101 |
| 7.2. | List of subsidiaries | 101 |
| 8. | Property, plant and equipment | |
| 8.1. | Significant existing or planned property, plant and equipment | 28, 33 |
| 8.2. | Environmental issues | 34 |
| 9. | Analysis of financial situation and results | |
| 9.1. | Financial position | 29 - 31, 91 - 92 |
| 9.2. | Operating profit | 24 - 26, 90, 102 - 106 |
| 10. | Cash position and capital | |
| 10.1. | Issuer's capital (short and long-term) | 46, 93, 113, 140, 145-146 |
| 10.2. | Cash flow | 29, 92 |
| 10.3. | Borrowing conditions and financial structure | 29 - 31, 121, 145 |
| 10.4. | Expected sources of financing | |
| 11. | Research and development, patents and licenses | 27 - 28, 162 - 163 |
| 12. | Information on trends | 47-49 |
| 13. | Profit forecasts or estimates | 47-49 |
| 14. | Governing, management, supervisory and executive bodies | |
| 14.1. | Composition of governing and management bodies Nature of all family links amongst these persons | 66 |
| 14.1.1. | Conviction for fraud within the last five years at least | 66 |
| 14.1.2. | Bankruptcy, sequestration or liquidation of a member of governing bodies | 66 |
| 14.1.3. | Indictment and/or official public sanction against a member of governing bodies | 66 |
| 14.2. | Conflict of interest at the level of governing and management bodies | 66 |
| 15. | Remunerations and benefits | |
| 15.1. | Remuneration paid and benefits in kind | 55, 79 -85 |
| 15.2. | Pensions, retirement or other benefits | 82-85 |

| Headings of Part 1 of European Regulation N°809/2004 of April 29, 2004 | | Pages |
|---|--|-----------------------|
| 16. | Operation of governing and management bodies | |
| 16.1. | Expiration date of current mandates and terms of office | 58 |
| 16.2. | Service agreements with the members | 66,67, 71 |
| 16.3. | Audit and Remuneration Committees of issuer | 69-70 |
| 16.4. | Corporate governance | 72-75 |
| 17. | Employees | 2 |
| 17.1. | Workforce at end of period covered by historical financial information | 2, 132 |
| 17.2. | Shareholding in capital | 54, 146 |
| 17.3. | Employee shareholding in capital | 52 |
| 18. | Principal shareholders | |
| 18.1. | Crossing thresholds | 52, 160 |
| 18.2. | Double voting rights | 52, 159-160 |
| 18.3. | Control of the business | 15, 132 |
| 19. | Transactions with related parties | 132-133 |
| 20. | Financial information on the Issuer's assets, financial situation and profit | |
| 20.1. | Historical financial information (results of the last five years) | 56 |
| 20.2. | 2016/2017 Consolidated Financial Results | 24 - 26, 90 |
| 20.3. | 2017 Statement of Profit or Loss | 90 |
| 20.4. | 2016/2017 Consolidated Financial Statements (Statement of Financial Position, Statement of Profit or Loss, Statement of Cash flows, Statement of Changes in Equity, Notes to the Consolidated Financial Statements) | 90-133 |
| 20.5. | Verification of annual historical information (Independent Auditors' Reports) | 134-137, 150-152 |
| 20.6. | Last financial information | 4 |
| 20.7. | Dividend distribution policy | 15, 50, 113, 145 |
| 20.8. | Legal and arbitration procedures | 42 |
| 20.9. | Significant change in financial or commercial situation | 24-33, 98-100 |
| 21. | Additional information | |
| 21.1. | Share capital | 51, 155 |
| 21.1.1. | Amount of issued capital, number of shares authorized, number of shares issued and fully paid, number of shares issued but not fully paid, par value per share and reconciliation of the number of shares outstanding at the beginning and end of the year | 51-52, 113, 145-146 |
| 21.1.2. | Shares not representing capital | 52 |
| 21.1.3. | Number, book value and face value of shares held by or on behalf of the issuer itself or by subsidiaries of the issuer | 51-52 |
| 21.1.4. | Amount of convertible securities, exchangeable securities or securities with warrants | 52 |
| 21.1.5. | Information about and terms of any acquisition rights or obligations over authorized but unissued capital, or an undertaking to increase the capital | 75-76, 113 |
| 21.1.6. | Information about any capital of any member of the group which is under option or agreed conditionally or unconditionally to be put under option | 76 |
| 21.1.7. | History of share capital for the period covered by the historical financial information | 51-52 |
| 21.2. | Memorandum and articles of association | 154 |
| 21.2.1. | Objects and purposes of the issuer | 154 |
| 21.2.2. | Member of administrative, management and supervisory bodies | 58-64 |
| 21.2.3. | Rights, preferences and restrictions attaching to each class of the existing shares | 155,159 |
| 21.2.4. | Actions required to change the rights of shareholders | 159-160 |
| 21.2.5. | Conditions governing the manner in which annual general meetings and extraordinary general meetings of shareholders are called including the conditions of admission | 155-156 |
| 21.2.6. | Provisions in the issuer's articles of association, statutes, charter or bylaws that would have an effect of delaying, deferring or preventing a change in control of the issuer | non applicable |
| 21.2.7. | Provisions in the articles of association, statutes, charter or bylaws governing the ownership threshold above which shareholder ownership must be disclosed | 159-160 |
| 21.2.8. | Conditions imposed by the memorandum and articles of association, statutes, charter or bylaws governing changes in the capital, where such conditions are more stringent than is required by law | 160 |
| 22. | Major contracts | 160 |
| 23. | Information provided by third parties, declaration by experts and declaration of interests | non applicable |
| 24. | Documents accessible to the public | 4 |
| 25. | Information on shareholdings | 101 |

Glossary

Industrial | Product-specific Terms

Biopharmaceuticals

Biopharmaceuticals, or biologics, are pharmaceutical drugs manufactured in or extracted from biological sources.

Bioreactor

In English-speaking countries, a bioreactor is used as a vessel for cultivating animal or human cells in a culture medium. In non-English-speaking countries, this term is also used synonymously with "fermentor" that is a system in which microorganisms (bacteria, yeast, fungi) multiply. In any case, these vessels are used to obtain cells, parts of these or one of their metabolites.

Downstream processing

Collective term for the various steps that follow fermentation or cell cultivation in the production of biopharmaceuticals; for example separation, purification and concentration

EMA – European Medicines Agency

European Union agency for the evaluation of medicinal products.

FDA – Food and Drug Administration

U.S. governmental agency responsible for monitoring foods and biotechnological, medical, veterinary and pharmaceutical products.

Fermentation

Technical process used to produce or transform intra- or extracellular substances with the help of microorganisms

Fluid management technologies

Technologies and systems for use in handling sensitive biological liquids; for example single-use bags for the preparation, storage and transport of biopharmaceutical solutions, intermediates and final bulk products

Membrane chromatography

Selective separation of mixtures of substances by adsorption to specifically modified membranes (membrane adsorbers) in a flowing system

Membrane

Thin film or foil made of polymers; because of its porous structure, this film is used as core component for all filtration applications.

Monoclonal antibodies

Synthetic antibodies that are increasingly being used in medical diagnosis and treatment

Purification

An important step in downstream processing

Single-use technologies

Technologies and products for a single use, providing significant time and cost savings; for example disposable filters or bags

Upstream processing

Upstream processing is defined as the entire process from early cell isolation and cultivation, to cell banking and culture expansion of cells until final harvesting. It refers to the part of the bioprocess in which cells or cell lines are grown in bioreactors (see bioreactor).

Validation

Systematic checking of essential steps and facilities in research and development and in production, including testing pharmaceuticals, to ensure that the products manufactured can be made reliably and reproducibly in the desired quality

Business | Economic Terms**Amortization**

Amortization relates exclusively to potential reductions in the value of goodwill and the allocation of the purchase price to intangible assets acquired as carried out in accordance with IFRS 3.

Cash flow

Cash balance of inflows and outflows of funds, representing the operating activities of an organization. Alternativ: Difference between the available cash at the beginning of an accounting period and that at the end of the period

Derivative financial instruments

Instruments for hedging against the risks of changes in market prices in foreign currencies

EBIT

Earnings before interest and taxes

EBIT margin

Ratio of EBIT (see EBIT) to sales revenue

EBITDA

Earnings before interest, taxes, depreciation and amortization.

EBITDA margin

Ratio of EBITDA (see EBITDA) to sales revenue

Extraordinary items

Extraordinary items essentially cover one-time expenses for corporate projects and integration and acquisition related items.

Fixed assets

Sum of intangible assets, property, plant and equipment and financial assets

Free float

Shares of a public company that are freely available to the investing public

Goodwill

Difference between the price paid for a company or business and its net assets. Goodwill is a form of intangible asset.

Investment rate

Ratio of capital expenditures to sales revenue

Normalized financial result

Financial result excluding fair value adjustments of hedging instruments, as well as currency effects from foreign currency loans

Normalized income tax

Underlying income tax, based on the underlying profit before taxes and non-cash amortization

Order intake

All customer orders which were legally concluded during the respective reporting period

Supply chain management

Setup and coordinated control of integrated flows of materials, information and finances (supply chains) over the entire value-added process

Treasury

Short- and medium-term liquidity management

Underlying EBITDA

EBITDA (see EBITDA) adjusted for extraordinary items (see extraordinary items)

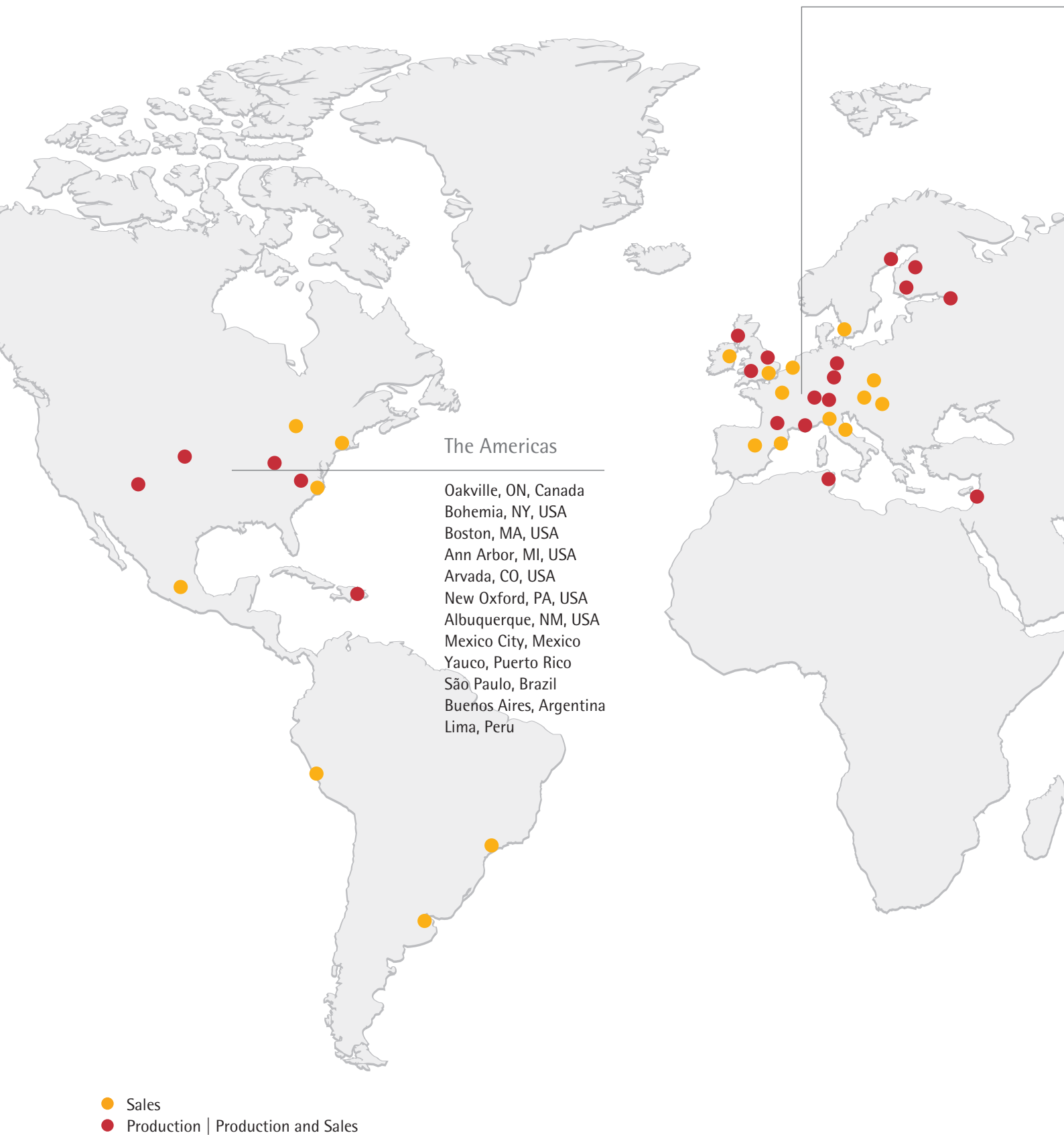
Underlying EBITDA margin

Ratio of operating EBITDA (see underlying EBITDA) to sales revenue

Underlying (consolidated) net profit

Profit adjusted for extraordinary items, non-cash amortization and based on the normalized financial result (see normalized financial result) as well as the corresponding tax effects for each of these items.

A Local Presence Worldwide



Europe | Africa

Dublin, Ireland
Royston, UK
Stonehouse, UK
Epsom, UK
Glasgow, UK
Brussels, Belgium
Dourdan, France
Aubagne, France
 Lourdes, France
Florence, Italy
Milan, Italy
Madrid, Spain
Barcelona, Spain

Kajaani, Finland
Helsinki, Finland
Malmo, Sweden
Umeå, Sweden
St. Petersburg, Russia
Goettingen, Germany
Guxhagen, Germany
Laupheim, Germany
Poznan, Poland
Budapest, Hungary
Vienna, Austria
Tagelswangen, Switzerland
Mohamdia, Tunisia
Beit Haemek, Israel

Asia | Pacific

Beijing, China
Shanghai, China
Hong Kong, China
Seoul, South Korea
Tokyo, Japan
Hanoi, Vietnam
Bangkok, Thailand
Bangalore, India
Kuala Lumpur, Malaysia
Singapore, Singapore
Melbourne, Australia

Financial Schedule

| | |
|--|------------------|
| Annual Shareholders' Meeting in Aubagne, France | March 26, 2019 |
| Payment of dividends ¹⁾ | April 2, 2019 |
| Publication of first-quarter figures for 2019 | April 18, 2019 |
| Publication of first-half figures for 2019 | July 19, 2019 |
| Publication of nine-month figures for 2019 | October 22, 2019 |
| Publication of preliminary figures for fiscal 2019 | January 2020 |
| Annual Shareholders' Meeting in Aubagne, France | March 24, 2020 |
| Publication of first-quarter figures for 2020 | April 2020 |

¹⁾ Subject to approval by the Annual Shareholders' Meeting

Contacts

Petra Kirchhoff
Head of
Corporate Communications & IR

Phone: 0551.308.1686
petra.kirchhoff@sartorius.com

Ben Orzelek
Head of Investor Relations

Phone: 0551.308.1668
ben.orzelek@sartorius.com

About This Publication

Published by
Sartorius AG
Corporate Communications
37070 Goettingen, Germany

Editorial Deadline
February 15, 2019

Published on
February 19, 2019

Financial Reporting System firesys
firesys GmbH
Frankfurt | Main, Germany

Photography
Peter Ginter
Lohmar, Germany
Frank Stefan Kimmel
Goettingen

This is a translation of the original
French-language
"Document de Référence 2018".

Sartorius Stedim Biotech S.A.
Zone Industrielle Les Paluds
Avenue de Jouques – CS 91051
13781 Aubagne Cedex, France

Phone: +33.4.42.84.56.00

Fax: +33.4.42.84.56.19

info@sartorius-stedim.com

www.sartorius-stedim.com

turning science **into solutions**