SVIFCT57S

Company profile

Sartorius is a leading international partner for the biopharma sector. Our solutions are supporting our customers to develop and produce drugs safely, timely and economically. Founded in 1870, Sartorius currently employs around 15,000 people in more than 30 countries. Its operational business is subdivided into two divisions: Bioprocess Solutions and Lab Products & Services.

Investment Highlights

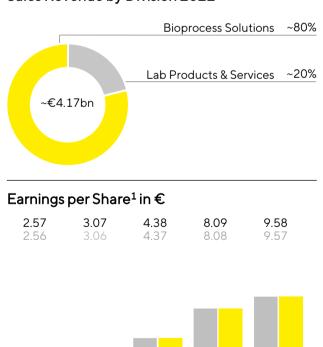
- 1. Clear focus on the attractive biopharma sector
- 2. Long-term growth drivers and significant market entrance barriers
- 3. Market leading position in key technologies and recognized brand
- 4. High share of recurring revenue as well as diversified earnings base
- 5. Proven track record with alliances and acquisitions; strong presence in growth regions

Key Figures	114 0000		111 0000	5,0000	51/0001	=
in millions of $ \in $ unless otherwise specified	H1 2023	∆ in %	H1 2022	FY 2022	FY 2021	FY 2020
Sales revenue (%∆ in const. fx)	1,735.2	-15.4	2,060.3	4,174.7	3,449.2	2,335.7
Order intake (%∆ in const. fx)	1,449.6	-32.7	2,168.8	4,007.3	4,267.9	2,836.3
Underlying EBITDA ²	516.7	-25.9	697.5	1,410.4	1,175.0	692.2
Underlying EBITDA margin ² in %	29.8	-4.1pp	33.9	33.8	34.1	29.6
Relevant net profit ³	202.5	-39.4	333.9	655.4	553.4	299.3
Underlying EPS per ordinary share ³ in \in	2.95	-39.4	4.88	9.57	8.08	4.37
Underlying EPS per preference share ³ in \in	2.96	-39.3	4.89	9.58	8.09	4.38
Equity ratio in %	38.4	+2.5pp	35.9	38.1	30.2	30.8
Net debt to underlying EBITDA ⁴	2.1		1.4	1.7	1.5	2.6

1 Relevant / underlying EBITDA: earnings before interest, taxes, depreciation and amortization and adjusted for extraordinary items

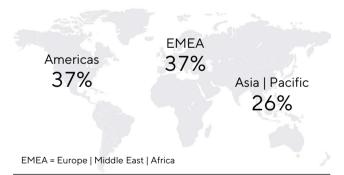
2022

2 After non-controlling interest, adjusted for extraordinary items as well as amortization, and based on a normalized financial result and the normalized tax rate 3 Ratio of net debt to underlying EBITDA of the past twelve months including the pro forma contribution of acquisitions for this period



Sales Revenue by Division 2022

Sales Revenue by Region 2022



2023 Guidance¹

	Sales Revenue Growth (in %)	EBITDA- Margin ²	
Sartorius Group	Decline in the low to mid- teens percentage range	~30%	
Bioprocess Solutions	Decline in the low to mid- teens percentage range	~31%	
Lab Products & Services	low single-digit negative to stable	~26%	
Capex Ratio	~15%		

1 In constant currencies

2 Underlying = excluding extraordinary items

1 After non-controlling interest, adjusted for extraordinary items as well as amortization, and based on a normalized financial result and the normalized tax rate

2020

2021

Preference shares (SRT3)

2018

Ordinary shares (SRT)

2019

SVILCILLS

Division Strategy

Bioprocess Solutions

- Positioned as a total solution provider for the biopharma industry with a product portfolio covering nearly all steps of the customers' production processes
- Global leading positions in key technology platforms; e.g. process filtration, fluid management, fermentation and cell cultivation
- One of the widest portfolios in the industry; clear focus on single-use technologies such as filters, bags and cell culture media
- Continuous expansion through complementary acquisitions and targeted alliances

Lab Products & Services

Sartorius Shares in Comparison (indexed)

July 1, 2022, to June 30, 2023

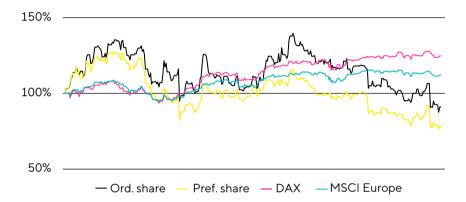
- Premium provider of laboratory instruments consumables and services for sample preparation and bioanalytic
- · Leading international positions in laboratory balances, pipettes and lab consumables
- Outstanding brand reputation, strong product range and global services as the basis for further expansion of market share
- Extension of portfolio with complementary laboratory products through acquisitions and alliances

Facts about Preference Shares

Ticker symbol: SRT3 Ticker symbol Bloomberg: SRT3 GY Ticker symbol Reuters: SATG_pe.DE WKN | ISIN: 716 563 | DE0007165631

Facts about Ordinary Shares

Ticker symbol: SRT Ticker symbol Bloomberg: SRT GY Ticker symbol Reuters: SATG WKN | ISIN: 716 560 | DE0007165607

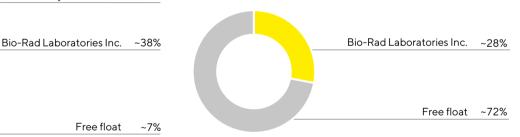


Shareholder Structure: Ordinary Shares

December 31, 2022

Administred by an executor ~55%

Shareholder Structure: Preference Shares December 31, 2022



The ownership relates to outstanding shares and thus excludes treasury shares. Information on shareholdings and shares in free float pursuant to Sections 33 et seq. of the German Securities Trading Act (WpHG) and the shareholders' own disclosures. The legal disclosure requirements refer only to ordinary shares and not to non-voting preference shares.

Financial Calendar

October 19, 2023 Publication of Quarterly Statement Jan.-Sep. 2023

Disclaimer

This fact sheet contains statements concerning the future performance of the Sartorius Group. 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 materialize. This is because our assumptions harbour 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. Throughout this fact sheet, differences may be apparent as a result of rounding during addition.

Petra Müller Head of Investor Relations Tel.: +49.551.308.6035 Email: petra.mueller2@sartorius.com