

Simplifying Progress

SRI Investor Presentation

April 2024



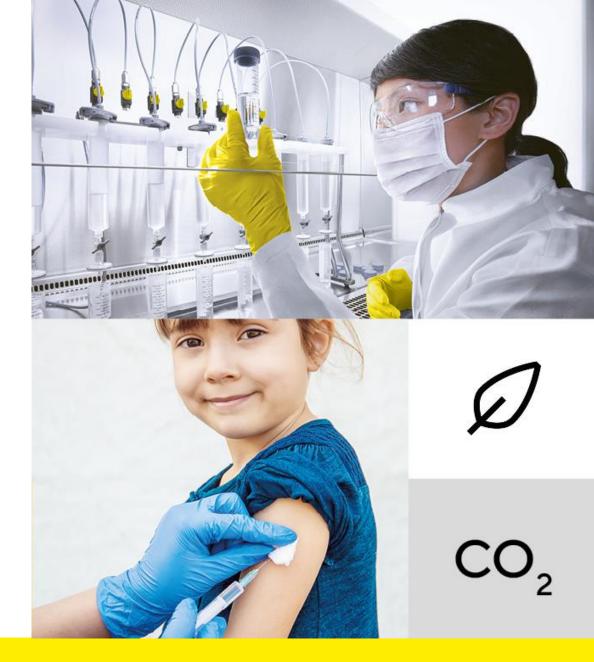
Agenda

Overview | Mission

Green sustainability

Social responsibility

Governance



Sartorius at a glance



~€3.4bn

Sales revenue 2023



~15%

Sales CAGR¹2013-2023



28.3%

EBITDA margin² 2022



60+

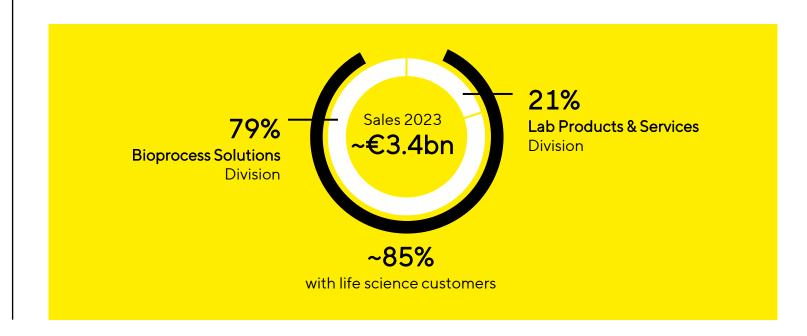
Locations worldwide, HQ in Göttingen, Germany



~14,600

Employees 12/2023

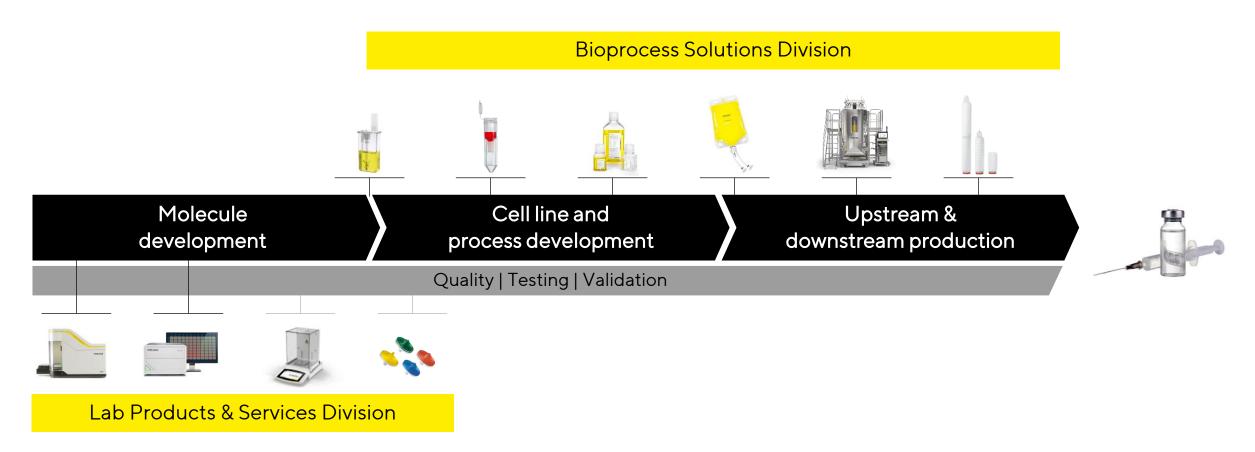
Two divisions with clear focus on attractive biopharma and life science industry



1 In constant currencies 2 Excluding extraordinary items



Supporting customers from drug discovery to commercial-scale production of biologics





Long-term value creation by focusing on four material topics





At the core of our business: "Good health and well-being"



Our mission

We empower scientists and engineers to simplify and accelerate progress in life science and bioprocessing, enabling the development of new and better therapies and more affordable medicine.



Health is one of the most relevant global topics

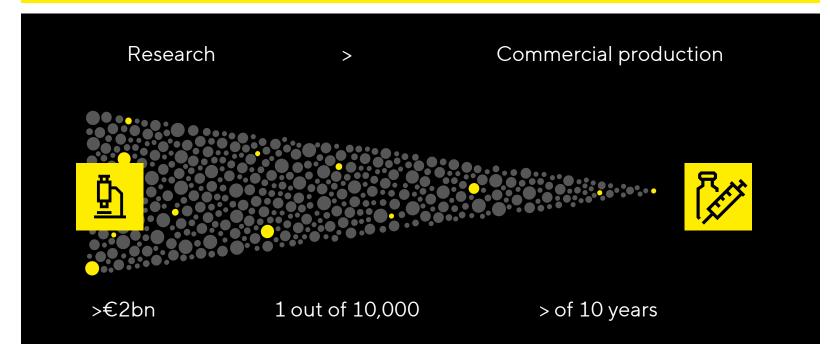
Growing and aging population

World population over 65 years¹
~ 1.6bn

~0.8bn

Q
2022 2050

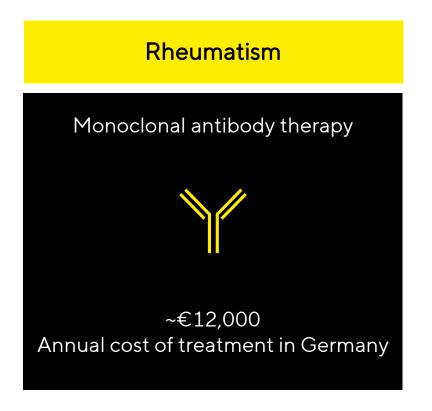
Development and manufacturing of drugs are time-consuming and costly²

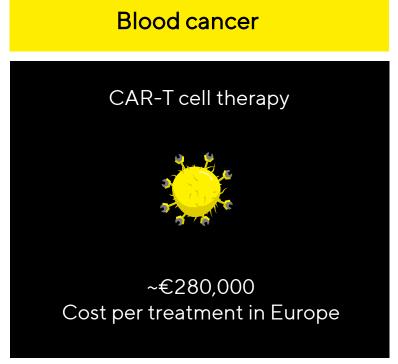


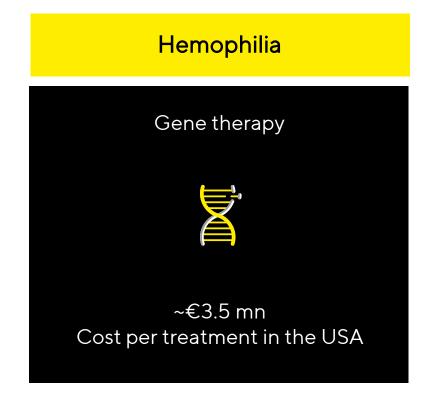
1 United Nations: World Population Prospects, 2022 2 Based on data from the Tufts Center for the Study of Drug Development and the Association of the British Pharmaceutical Industry



The challenge: Many innovative biotech medications are very expensive





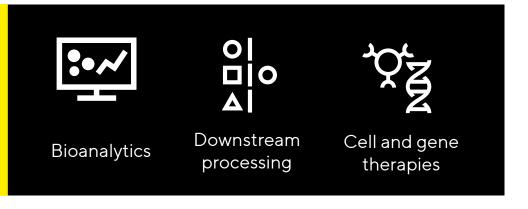


Three drivers for innovation: The key to achieve ESG targets



Acquisitions

of complementary and differentiating technologies



Cooperations

with partners that are leading in areas of future relevance

Advanced materials

Process automation & digitization

3D cell printing technology

Bioanalytics



Sartorius' technologies enable researchers to achieve results faster, ...

Example: digitalization and automation of laboratory processes







iQue®3



- More experiments with lower error rate
- Data and analysis for better decisions
- Shortening time-to-clinic



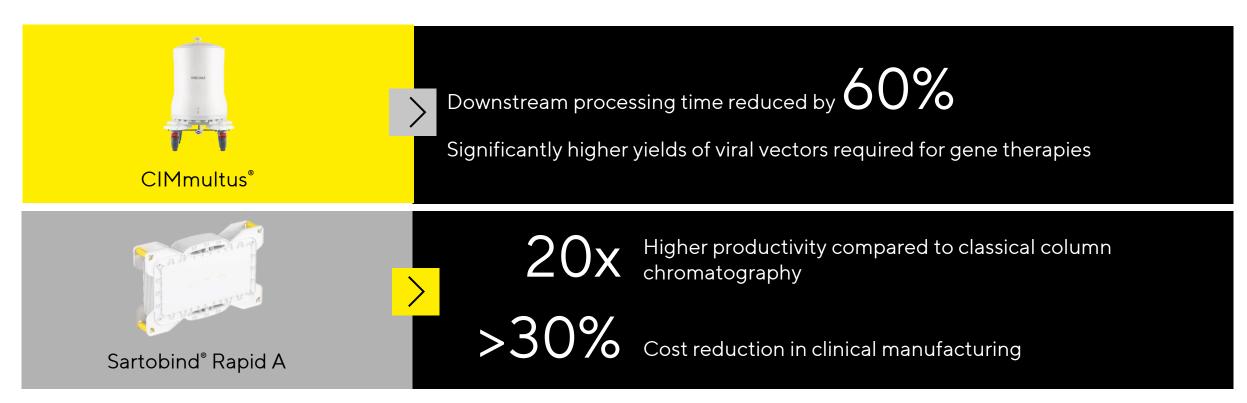




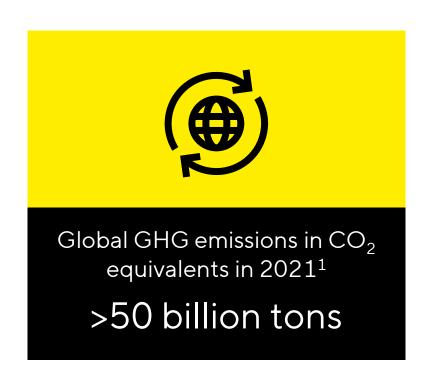
Saves more than **5** weeks in the selection of cell clones

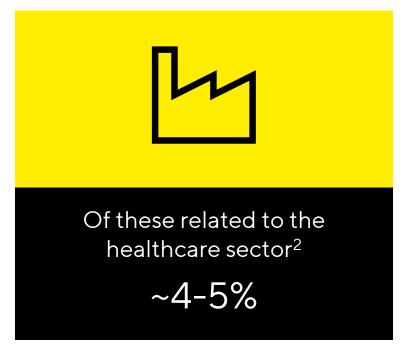
... to increase yield and efficiency in biopharma production ...

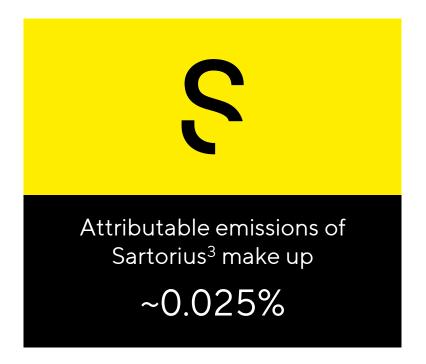
Example: new, innovative bioprocess technologies



A strong lever of impact on green aspects: Our customers' operations



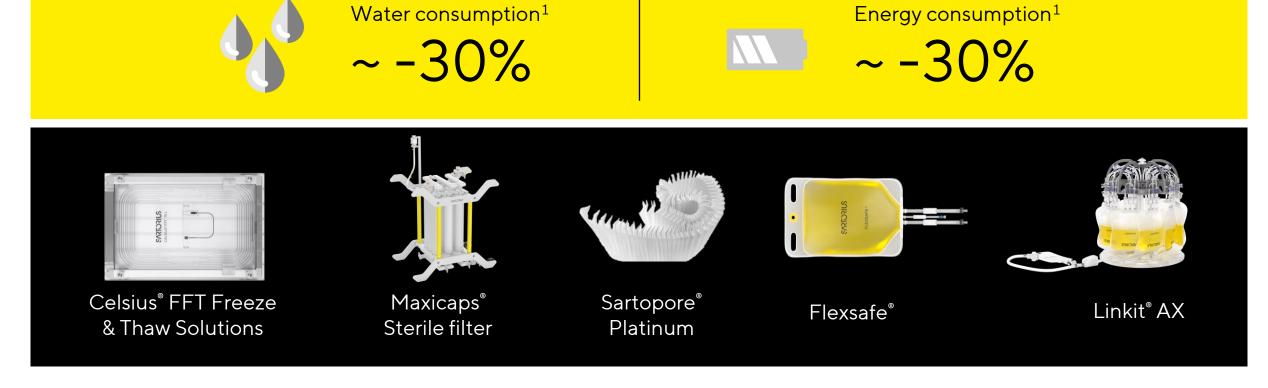




1 UN Environment Programme "Emissions Gap Report 2022", preliminary estimate 2 According to World Economic Forum June 2021 3 Own calculations



SU technologies help to reduce water and energy consumption at the customers' sites



1 Based on data from Flanagan et al (2011). LCA of a mAb process at 2,000 L scale.

Process intensification leads to higher efficiency and less eco impact

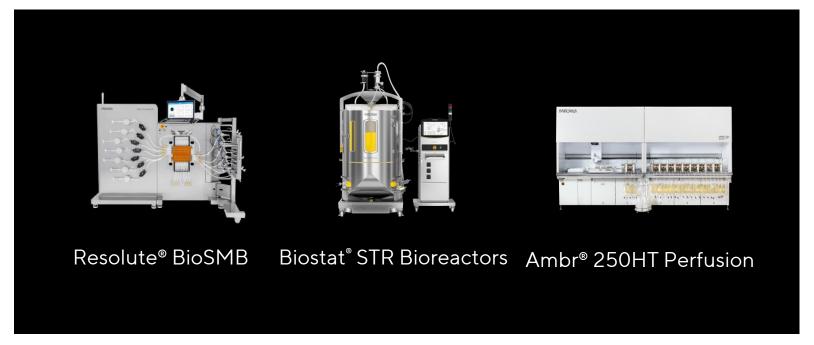


Process Mass Intensification¹

-50%

PMI = Total raw materials, consumables and water (kg)

Product (kg)



1 Data from in-house case study: 300 kg/year and 1,500 kg/year throughputs. PMI measured for a traditional stainless steel fed batch process (SS 15kL), and a process intensification scenario (SU dynamic perfusion process with downstream multi-column chromatography. Data was generated using BioSolve modeling software.



Data analytics solutions support customers in further CO₂ reduction

Success Story

Implementation of Umetrics® Suite data analytics tools at Michigan University helps increase process efficiency and reduces carbon footprint significantly

Phase 1 | prevented

780

tons CO₂e

Phase 2 | aims to prevent

10,000

tons CO₂e



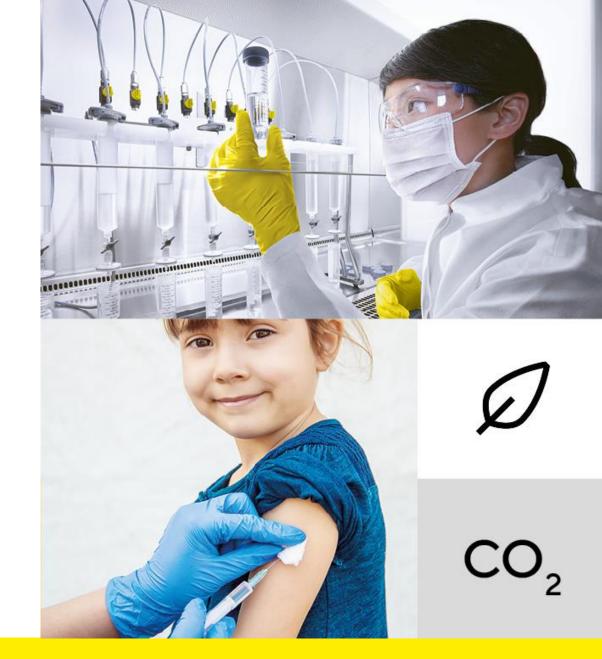
Agenda

Overview | Mission

Green sustainability

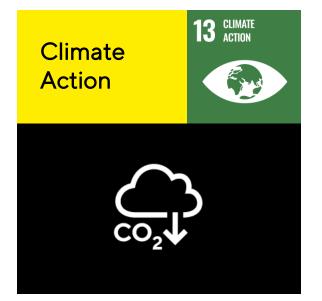
Social responsibility

Governance



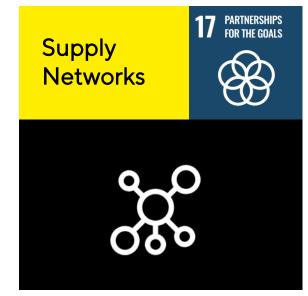


Sartorius' environmental fields of action at its own operations









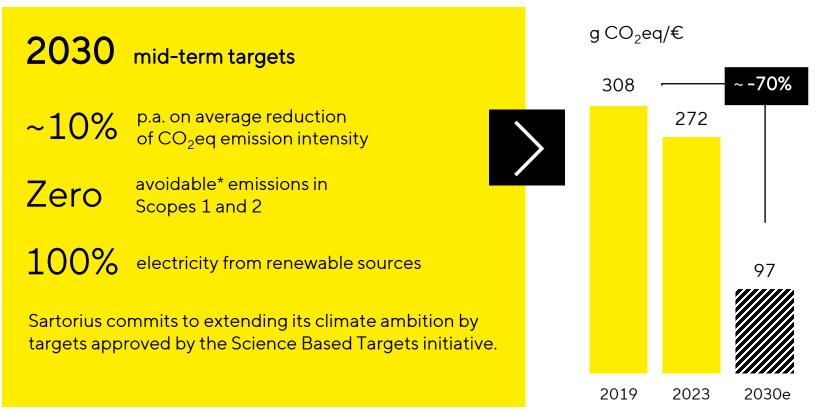
Only a small portion of emissions under own control





Ambitious mid- and long-term climate targets





2045 long-term target Net-zero emissions Decarbonization of the entire supply chain in collaboration with suppliers and customers

^{*}Process emissions from membrane manufacturing are currently considered unavoidable based on available technology.



Additional sustainability target: Climate neutral by 2045

Long-term target complements existing climate strategy

- Decarbonization in collaboration with suppliers and customers
- 100% electricity from renewable sources by 2030
- Electricity from hydropower at German sites in Göttingen, Guxhagen and Ulm since 2021





Sartorius commits to extending its medium-term climate related ambition by targets approved by the Science Based Targets initiative.



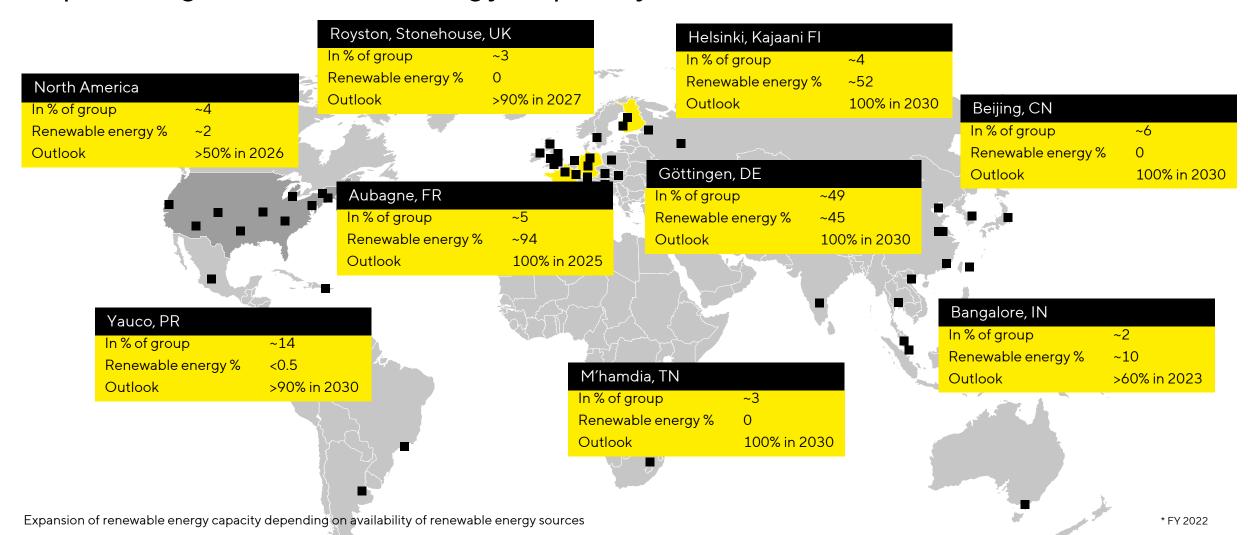
CO₂ reduction levers across the value chain

Upstream **Production** Downstream (Scope 3) Scope 1&2 (Scope 3) **P** Other (capital goods, Purchased Upstream incl. Down-stream Product use & **Business** Manufacturing employee commuting, goods & services int. transportation travel & warehousing transportation end of life buildings) Eco-design products & Eco-design Reduce travel Improve energy Eco-design Eco-design products Buildings & packaging packaging efficiency & packaging construction packaging Sustainable Efficiency in travel Produce clean Process intensification Capital goods Transport Transport Frequency Frequency manufacturing energy Distance Distance Commute Fright mode Fright mode Supplier engagement Procure sustainably Fuels Fuels renewable energy Production process steps





Expanding renewable energy capacity at Sartorius' sites worldwide





Flexsafe®- first film designed for recycling



A milestone on the way to circularity

- From 20ml to 2,000l, Flexsafe® is the first film in the industry designed for recycling in a standard polyethylene waste stream
- Used in bioprocessing for non-hazardous applications, it can technically be recycled into raw materials with quality equivalent to pristine grades of polymers.

Partners









Materials and circularity: Transitioning towards a circular economy



Our ambitions

We aim to optimize the selection and use of materials to improve our products' environmental footprint.

- Optimize use of resources
- Avoid packaging waste
- Recycle production waste
- Life cycle analysis

2030 Circularity targets

>75%

Sales revenue with products designed according to circularity principles

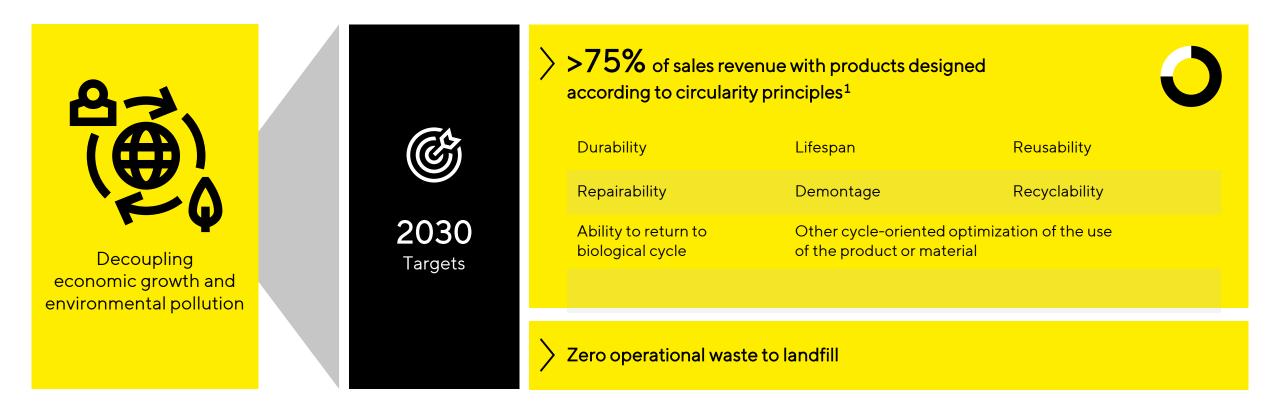
Zero

operational waste to landfill



1 including product and transport packaging

Circular economy: New, medium-term goals defined



1 including product and transport packaging





Current projects regarding the single-use technology life cycle

| Feedstock Beginning of life | Production | Product usage | Product end of life |
|--|-----------------------------------|------------------|---------------------|
| Resins Fossil Naphtha Crackers → Resins | Product Packaging | $\Rightarrow \\$ | |
| Alternative Feedstock 1 | Scrap Reduction Recycling Program | PMI 4 | Recycling Options 5 |
| Less Material 2 | | | |









¹ Figure taken from: Vachette, Elisabeth & Fenge, Christel & Cappia, Jean Marc & Delaunay, Lucie & Greller, Gerhard & Barbaroux, Magali. (2014). Robust and convenient single-use processing: The superior strength and flexibility of flexsafe bags. 12.



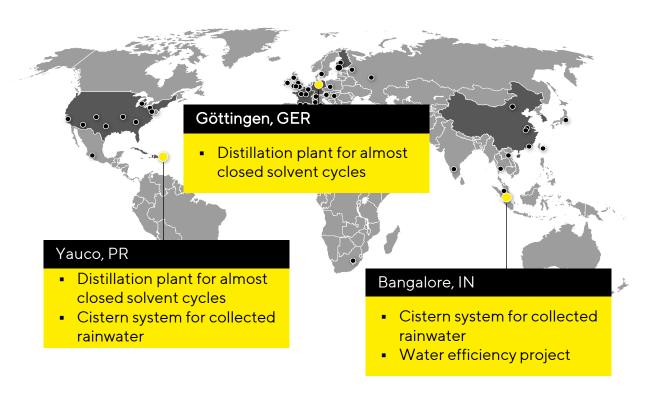
Water and wastewater: Improve water efficiency in own operations



The ambitions

- Reduce use of drinking water, increase use of rainwater
- Decrease wastewater, increase water recycling
- Optimization of production cycles with solvents





Solvent and water reuse at the Yauco plant



4 distillation columns
Processing capacity: 16,500 kg/day of spent solvent (oldware)

3.9 million kg of oldware processed for reuse

 $11.9 \ \text{million litres}$ of water processed for reuse



90% recovery of 2-pyrrolidone (2P) from oldware



98% recovery
of Isopropanol (IPA) from oldware

Base April 2022



Supply chain due diligence act: Risk assessment





1 Coverage in terms of spend as of June 2023



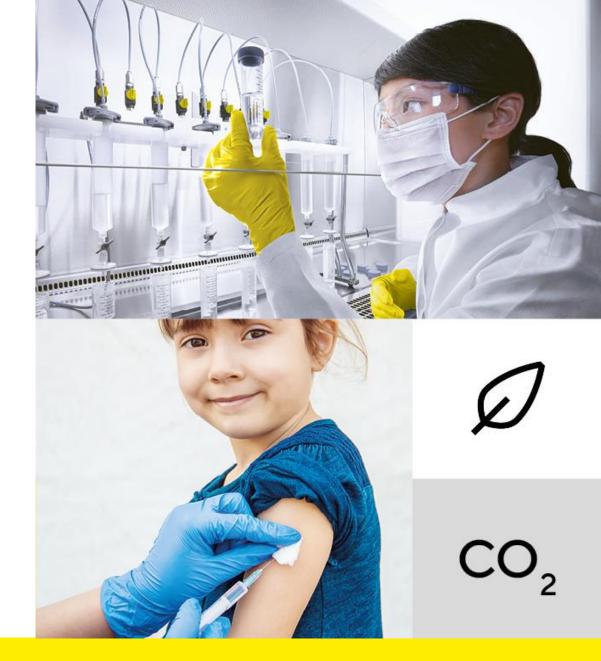
Agenda

Overview | Mission

Green sustainability

Social responsibility

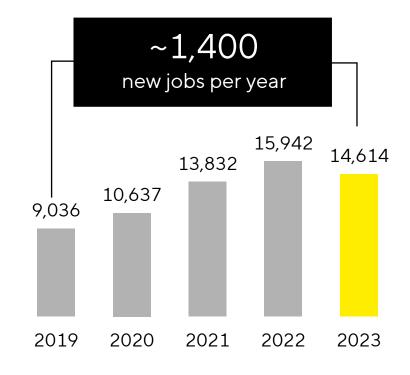
Governance



Strong increase in personnel over 5 years: 62% more employees

After strong team growth, current focus on efficiency





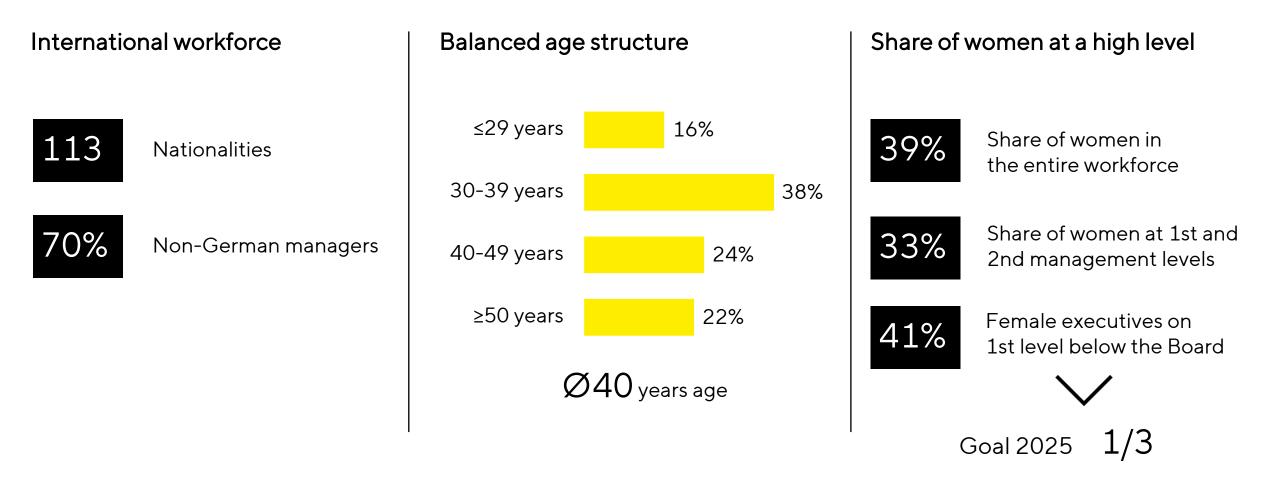
Many opportunities to grow with the company



of management positions filled internally



An increasingly diverse team that brings in a variety of perspectives





High employee satisfaction reflected internally and externally

87%

say they contribute to the company's mission

77%

are proud to work for Sartorius

82%

agree that Sartorius respects and strengthens diversity

3.6 ★★★★☆

Jglassdoor

Intern. competitors \emptyset 3.7

4.0 ★★★★☆

kununu

DAX40 Ø 3.8



As of Q1 2024

Promoting science, education and innovation – also beyond our walls

Cell Culture Facility at Penn State | Supporting groundbreaking research with \$1.5 million

Life Science Factory |
Promoting startups and attract
entrepreneurs in life sciences

Creating a stage for science communicators since 2021







Engaged in humanitarian aid and disaster relief globally

action medeor provides medical care to humanitarian organizations and health centers all over the world.

Our latest commitments include

- Support for earthquake victims of in Türkiye and Syria
- Humanitarian aid for the people of Ukraine
- Healthcare support in Guatemala
- Pharmaceutical training in West Africa



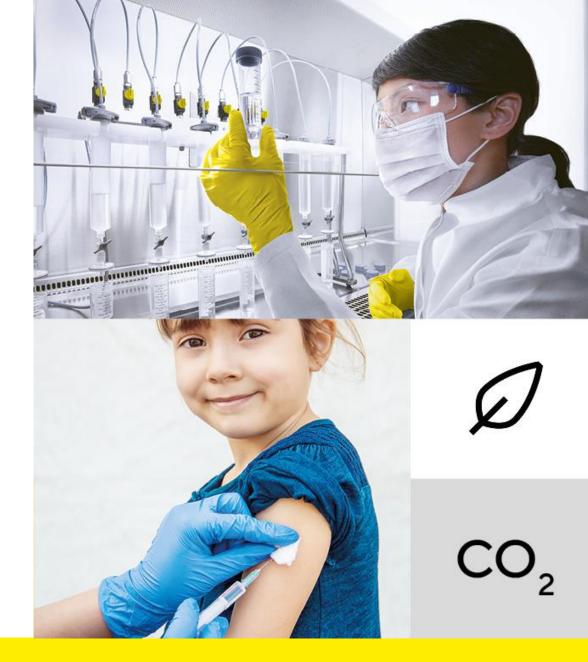
Agenda

Overview | Mission

Green sustainability

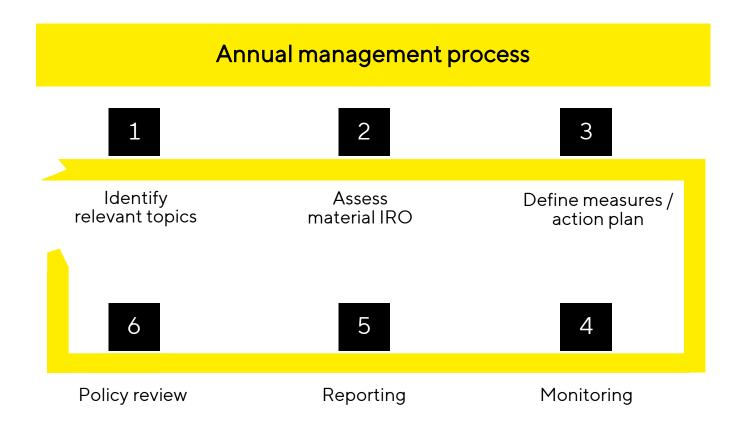
Social responsibility

Governance



Sartorius actively manages its sustainability impacts, risks and opportunities at its own operations

Existing and upcoming EU regulation Corporate Sustainability Due Diligence Directive (CSDDD) Taxonomy Regulation Corporate Sustainability Reporting Directive (CSRD)





Multiple data points to be consistently tracked consistently



Challenges

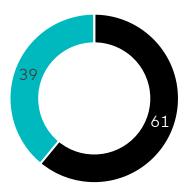
- > 80 disclosure requirements
- > 1.000 data points
- 17 GHG emission categories
- Data from multiple sources
- Concepts for data calculation and estimates
- Data governance and enrichment projects
- Continuous data improvement

Reporting partially based on external data which can be subject to changes and specifications; therefore, continuous and retrospective restatements of certain data might be necessary.



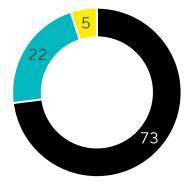
Alignment of 2023 turnover & expenditures with EU taxonomy





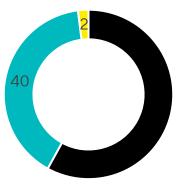
- Not Taxonomy-eligible
- Taxonomy-eligible but not aligned
- Taxonomy-aligned

Total CapEx in %



- Not Taxonomy-eligible
- Taxonomy-eligible but not aligned
- Taxonomy-aligned

Total OpEx in %



- Not Taxonomy-eligible
- Taxonomy-eligible but not aligned
- Taxonomy-aligned

All figures rounded

1 Sartorius does not have any taxonomy-eligible economic activities for which it would currently have been possible to assess compliance with the technical screening criteria as legislation only addresses particularly relevant economic activities for the achievement of the environmental objectives in the field of climate change mitigation and climate change adaptation, i.e. only a limited range of industries is covered.



Sustainability ratings and partnerships

Listed in

Memberships and cooperation in internationally renowned initiatives and platforms

DAX 50 ESG

Sustainability ratings







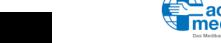






































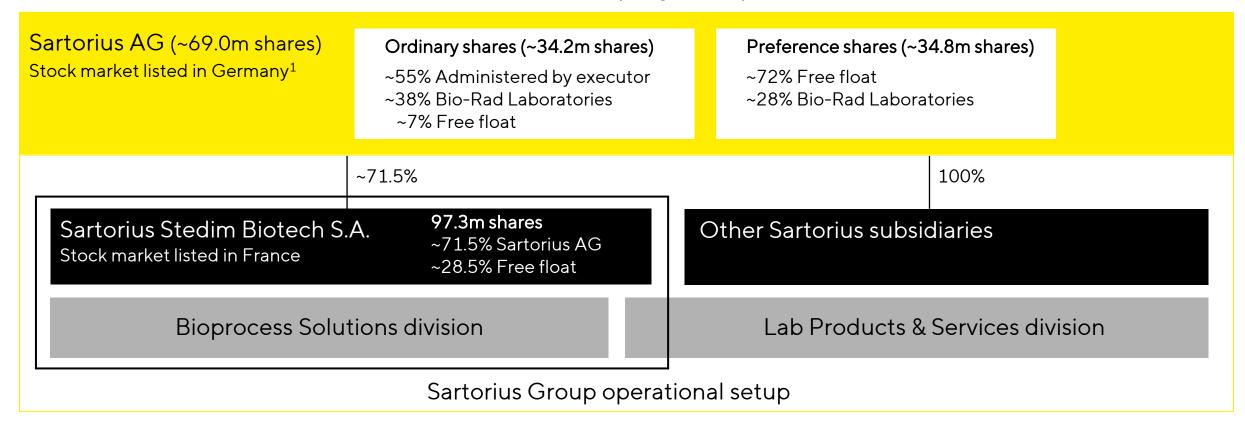






Sartorius legal structure

Sartorius Group legal setup



¹ The ownership relates to outstanding shares and thus excludes treasury shares. Information on shareholdings and shares in free float pursuant to the disclosure requirements of Sections 33 et seq. of the German Securities Trading Act (WpHG) and the shareholders' own disclosures. The legal disclosure requirements refer only to voting shares and not to non-voting preference shares.



Executive Board of Sartorius AG



Joachim Kreuzburg, CEO

Joined Sartorius in 1999 Member of the Board since 2002

Group Strategy, Corp Research, HR, Legal & Compliance, Communications, Sustainability



René Fáber

Joined Sartorius in 2002 Member of the Board since 2019 Bioprocess Solutions Division



Alexandra Gatzemeyer

Joined Sartorius in 2005 Member of the Board since May 1, 2023

Lab Products & Services Division



Florian Funck, CFO

Joined Sartorius in 2024 Member of the Board since April 1, 2024

Finance, Information Technology, Data Management, Corporate Sourcing

Two-tier Board system of Sartorius AG – Supervisory Board and its committees

Twelve members, half of the members are shareholders` representatives and half are employees' representatives

Executive Task Committee

Dr. Lothar Kappich (Chairman)

Annette Becker

Prof. Dr. Klaus Rüdiger Trützschler

Manfred Zaffke

Audit Committee

Prof. Dr. Klaus Rüdiger Trützschler (Chairman)

Dietmar Müller

Dr. Lothar Kappich

Manfred Zaffke

Nomination Committee

Dr. Daniela Favoccia

Dr. Lothar Kappich

Prof. Dr. Klaus Rüdiger Trützschler

Conciliation Committee

Dr. Lothar Kappich (Chairman)

Annette Becker

Prof. Dr. Klaus Rüdiger Trützschler

Manfred Zaffke



One-tier Board of Sartorius Stedim Biotech S.A. subgroup







Dr. René Fáber CEO



Romaine Fernandes Employee representative



Pascale Boissel



Susan Dexter



Anne-Marie Graffin Lead Independent

Director



Lothar Kappich



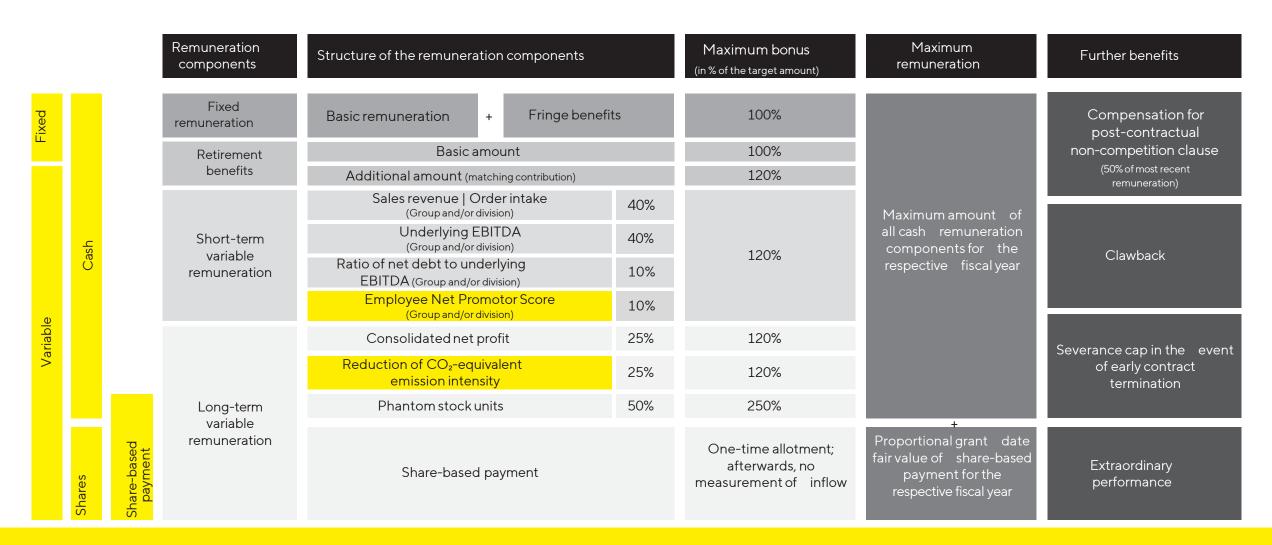
Henri Riey

Share of independent members

| Audit Committee | • • • • | 75% |
|---|---------|-----|
| Remuneration & Nomination Committee | • • • | 75% |
| Independence | • • | 43% |



ESG aspects included in Sartorius AG executive remuneration





Corporate responsibility & further financial information

To Sartorius, sustainability means operating responsibly over the long term - with respect to business partners, employees and society as well as to natural resources. Learn more about our commitment:











Please click below for further financial information:







Corporate Governance Report 2023

Earnings Release 9M 2023

Half Year Report H1 2023

Contacts and Financial calendar



Petra Kirchhoff

Head of Corporate Communications & IR

+49.551.308.1686

petra.kirchhoff@sartorius.com



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



Mirko Koch

Manager Investor Relations
+49.551.308.2402
mirko.koch@sartorius.com

Sartorius Corporate Administration GmbH Otto-Brenner-Straße 20 37079 Göttingen, Germany

- April 28, 2024
 Publication Q1 2024 Results
- May 16, 2024
 Capital Markets Day | Goettingen
- May 30, 2024
 Kepler Cheuvreux ESG Conference | Virtual
- Jul 19, 2024
 Publication Half-Year Report 2024
- Oct 17, 2024
 Publication 9M 2024 Results



Disclaimer

This presentation 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 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.

Throughout this presentation, differences may be apparent as a result of rounding during addition.

