

Simplifying Progress



Sartorius Stedim Biotech
Capital Markets Day 2026

March 17, 2026

SARTORIUS



Note

This presentation contains statements concerning the future performance of Sartorius Stedim Biotech. 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.

Transcripts or recordings of this event are not permitted.



What we are building on



Anchored in purpose – Ambitious to shape the future

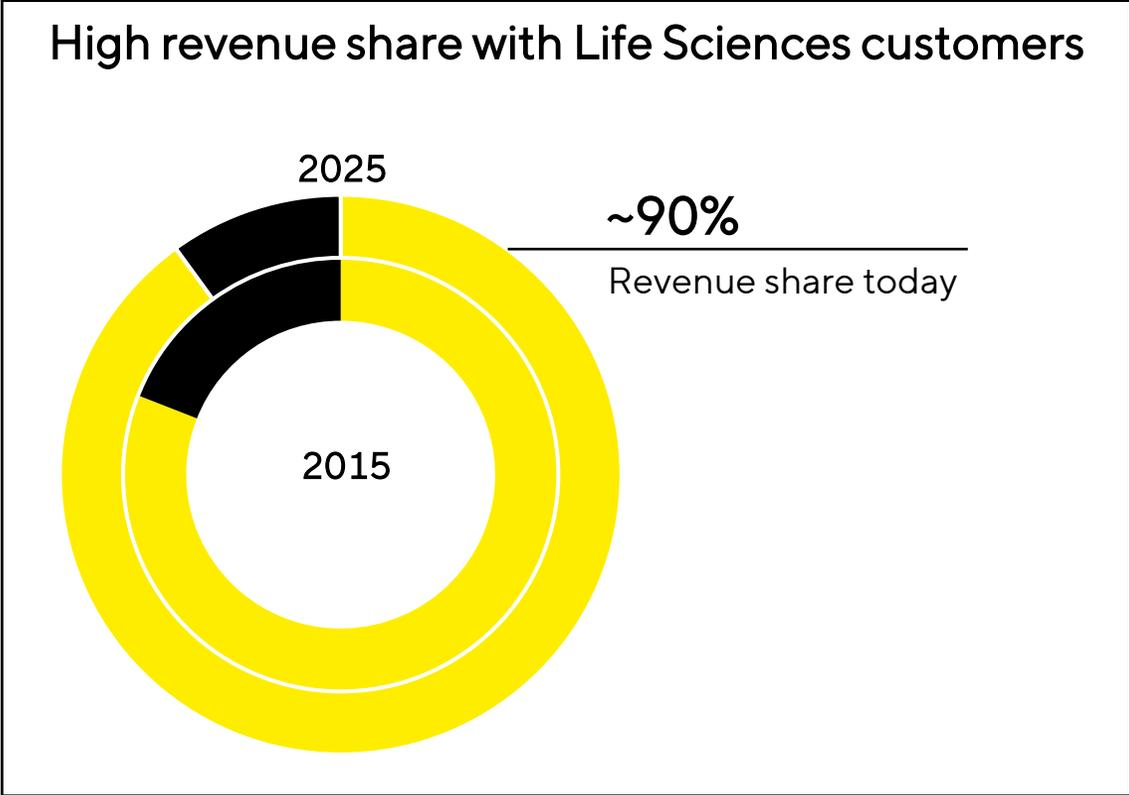
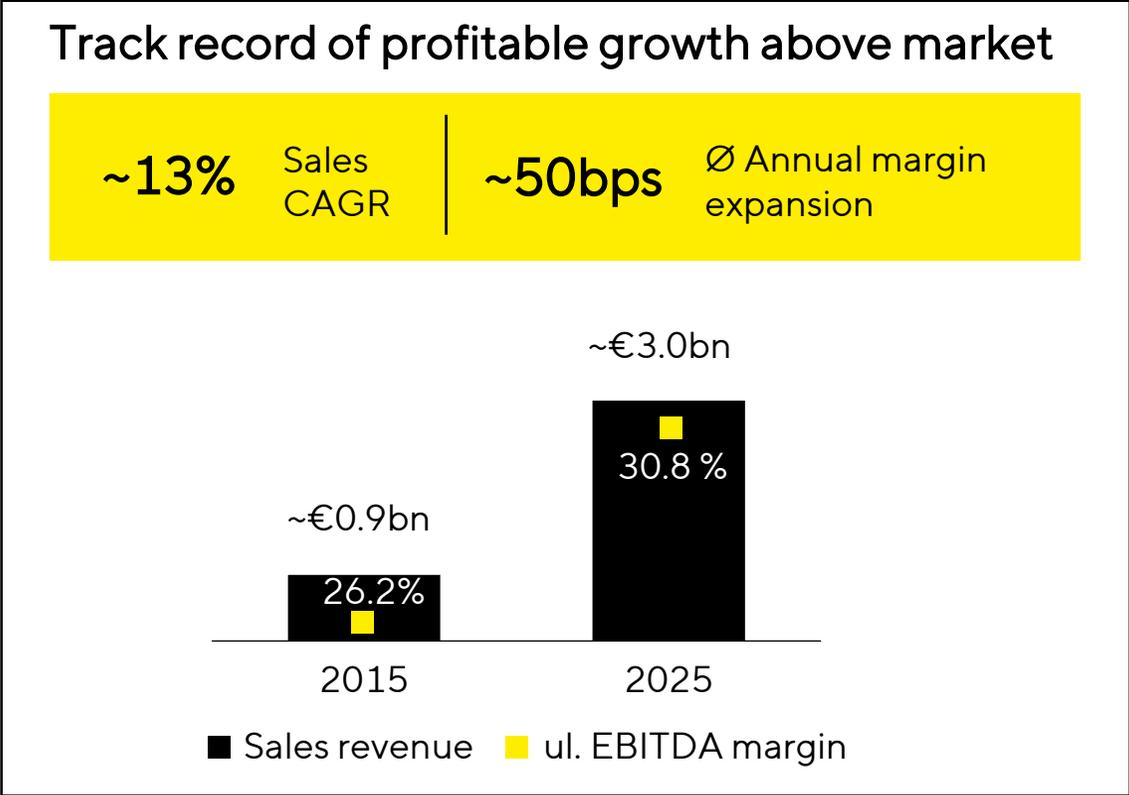
Our Purpose	Better health for more people.	
Our Mission	We empower biopharma customers to accelerate drug development and enhance manufacturing efficiency for better therapies and more affordable medicine.	
Our Promise	Simplifying Progress	



A leading biologics pure play with unique strengths in single-use

 Strong focus on biopharma	Highly differentiated end-to-end portfolio; leading market positions
 Unique know-how	Deep scientific, application and regulatory expertise
 Trusted supplier	Superior and consistent quality and global technical support
 Global footprint	Strong presence in all regions, resilient supply chains
 Long-term partnerships	With scientific research institutions, universities, and customers

Compelling business profile in a ten years perspective



Continuous portfolio expansion through innovation

Strengthened portfolio in three key areas mainly through M&A

 **Chromatography**

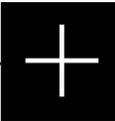
- Closed gaps in downstream offering
- True end-to-end portfolio

 **Solutions for advanced therapies**

- Built an early position in emerging market
- Focus on mission-critical, high-value applications

Strong innovation engine

 **Own product development**
in areas of core competencies



 **Cooperations**
with customers, research institutes,
start-ups, and industry-leading experts

Addressing the complex and regulated biopharma value chain

1 Cell line & process development

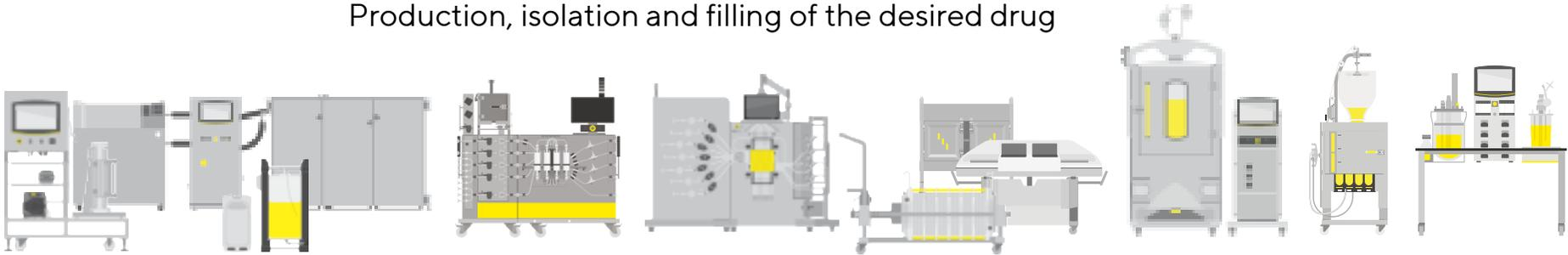
Creation of a stable cell line;
definition of a robust
manufacturing process



Sartorius offers an
integrated, fully scalable
end-to-end portfolio

2 Upstream & downstream production

Production, isolation and filling of the desired drug



A scalable single-use platform for efficient biopharma manufacturing

 **Our ambition**
More efficient bioprocessing

 Scalable end-to-end solutions for all biopharmaceutical modalities;
Market leader in single-use technologies

Design safe and flexible processes

Reduce investments

Increase productivity

Enhance product yield

Media & reagents



Bioreactors



Chromatography



Filtration



Fluid management



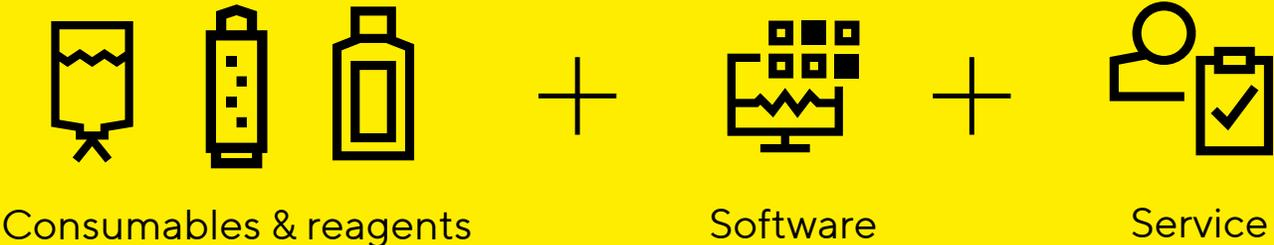
Digital solutions



Note: Highlighted products are a sample from our end-to-end portfolio

Attractive business model with high share of recurring revenue

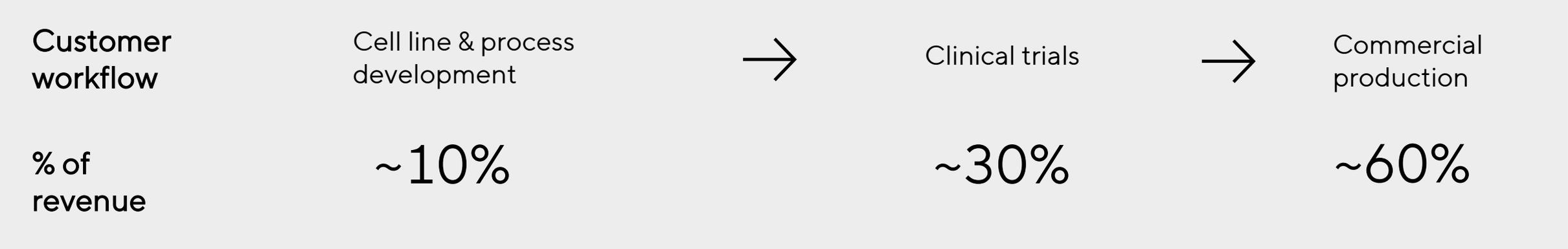
Steady recurring revenues supported by validated manufacturing processes and equipment installations



~85%
Recurring revenue

FY 2025 figures

Regulatory validation as a foundation for the resilient Sartorius business model



Note: Sales split by development phase based on Group sales with Life Science customers only (~90% of total Group sales)

Strong market fundamentals

Sustained strong market fundamentals



Growing and aging population¹

>1.6bn

Estimated number of people aged 65+ by 2050 (2x vs 2025)



Growing pharma market²

~5% p.a.

growth of global pharma market



Shift in medicine to biologics²

~57%

share of biologics in pharma market 2030e (+~22pp vs 2015)



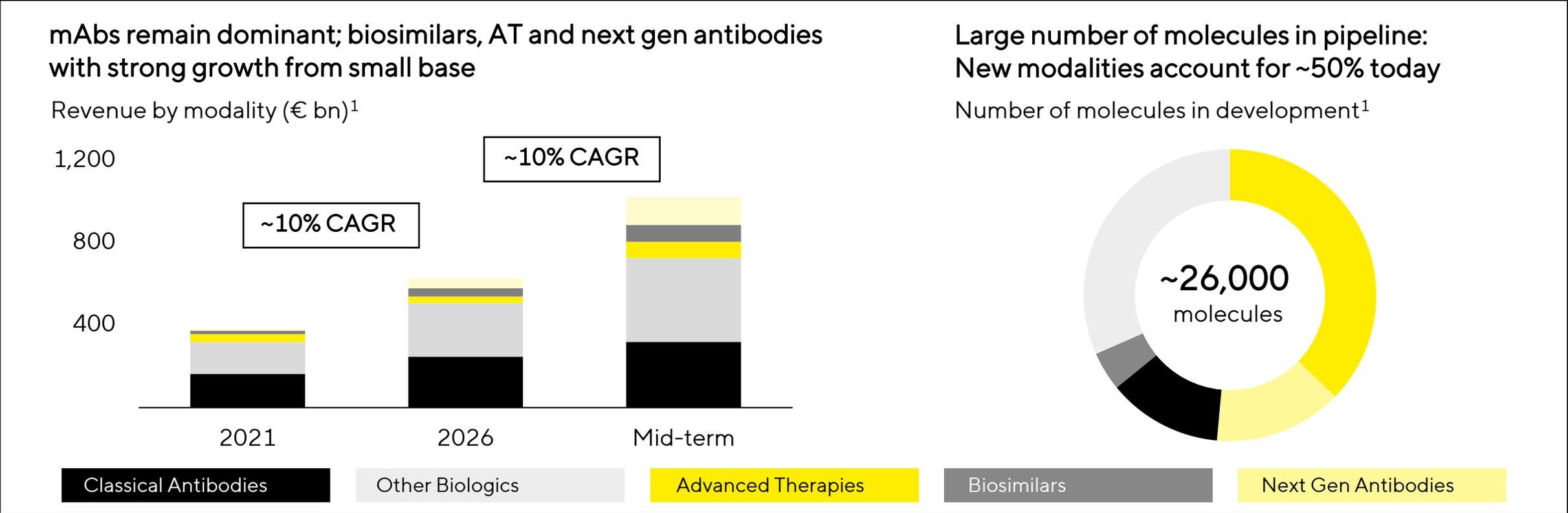
Increasing number of drug approvals^{2,3}

2x

number of approvals⁴ in 2020-2025 (~890) vs 2015-2020 (440)

Sources: 1 WHO 2 Evaluate Pharma, April 2025, GloblaData , January 2026 3 Sartorius market intelligence 4 Global first-time approvals of biologics

Broad growth of biologics; rising relevance of new modalities



¹ Global Data, IQVIA, Market Reports & Sartorius Market Intelligence January 2026

R&D and production volumes as structural market growth drivers

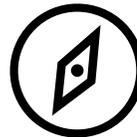
Addressable market CAGR mid-term¹

8-10%



Opportunities

- Biopharma R&D and production volumes continue to grow, driven by expanding pipelines, drug approvals, and rising patient demand
- Increasing complexity and diversity of modalities drive investment in technological innovation
- AI and technology are accelerating the pace of change



Challenges

- Pressure on healthcare systems and drug pricing
- Geopolitical dynamics and policy shifts add complexity and friction

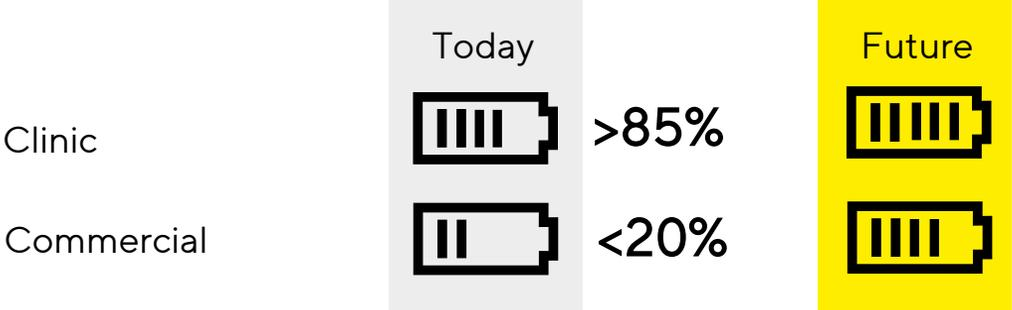
Source: Global Data, IQVIA, Market Reports & Sartorius Market Intelligence January 2026

Single-use (SU) and process intensification: Reshaping bioprocessing

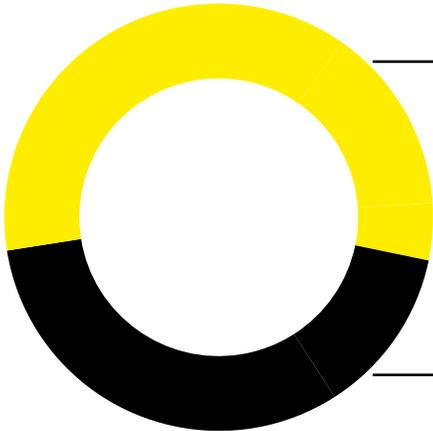
Biologics manufacturing is evolving

- More single-use, even at commercial scale
- More intensified, automated, and data-driven
- More modular, flexible, and regionalized
- Less capital-intensive, more sustainable and productive

Single-use penetration:



Modality mix supports adoption: SU is preferred platform for majority of pipeline modalities



~55%
Predominantly SU both for clinical and commercial applications

~45%
Largely hybrid manufacturing, SU prominent in clinical applications

Source: Global Data, Sartorius Market Intelligence January 2026

How we shape the future

Strengthening Sartorius Stedim Biotech's competitive edge

Financial ambition

Profitable growth above market, enhanced cash generation, reduced leverage

Strategic initiatives

Portfolio

How to add value to customers

Focus on innovation, core strengths and emerging opportunities

Customer experience

How to engage with customers

Trusted partnership, fast and reliable technical support and high service quality

Excellence & efficiency

How to operate

More cost-efficient, digital, automated and lean to provide short lead-times and high delivery reliability

Enablers

People engagement & culture

AI, data and digitalization

Empowerment

Sharpening portfolio for above-market growth

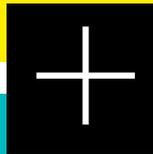


Our Goal

We stay focused on biopharma customers' workflows and strengthen our positioning through continued innovation leadership

Grow the core

Grow ambitiously and expand market leadership in our core businesses



Expanding into future businesses

Establish leading positions in emerging fields and technologies

- Focus on high-value, mission critical applications and/or high growth areas
- Capture greater share of wallet
- Focus on organic business development and partnerships
- Leverage data, AI, and automation technologies across the portfolio

Growing the core across mission-critical bioprocessing and cell analytics

 <p>Process Intensification</p>	<ul style="list-style-type: none">▪ Build on our pioneering position to lead single use manufacturing	<p>Pionic</p> 	<p>Ambr</p> 
 <p>Core Single Use Technologies</p>	<ul style="list-style-type: none">▪ Maintain leadership through continuous innovation of established product lines▪ Strengthen positioning in critical applications	<p>Sartopore Evo</p> 	<p>Celsius Pak</p> 

Advancing future businesses beyond the core: Additional growth vectors



ATS¹

Specialized solutions built on differentiated raw materials, analytics, and instrument expertise



PAT & AC/QC³

Integrated PAT and in-/at-line analytic platforms in manufacturing and quality labs



1 Advanced Therapy Solutions 2 ACM = Advanced Cell Models, NAM = New Approach Methodologies 3 PAT = Process Analytical Technology, AC/QC = Analytical Characterization & Quality Control

Enhancing customer experience across the end-to-end journey

Customer expectations across the life science industry



Short lead times, convenient product customization



Reliable delivery and high quality products



Simplified ordering and service workflows



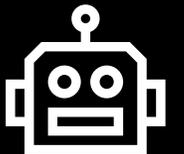
Comprehensive & accessible online solutions

Initiatives to enhance customer experience

Improving lead times, delivery performance and configurability of our product to enhance customer value.



More automated and digitalized customer journey including self-service platforms, proactive messaging and predictive relevance to our customers



Evolving into a more cost-efficient, digitalized and lean organization



Operations transformation to achieve best-in-class supply chain performance, product and service quality

<p>Factory of the future</p> <p>more automation, digital tools, data-driven quality, fewer manual steps</p>	<p>Network optimization</p> <p>resilient and efficient global production footprint</p>	<p>Simplify operations</p> <p>clearer flows, less duplication, more standardization</p>	<p>Strengthen sourcing</p> <p>reduce risk, improve availability and cost positions</p>
--	---	--	---

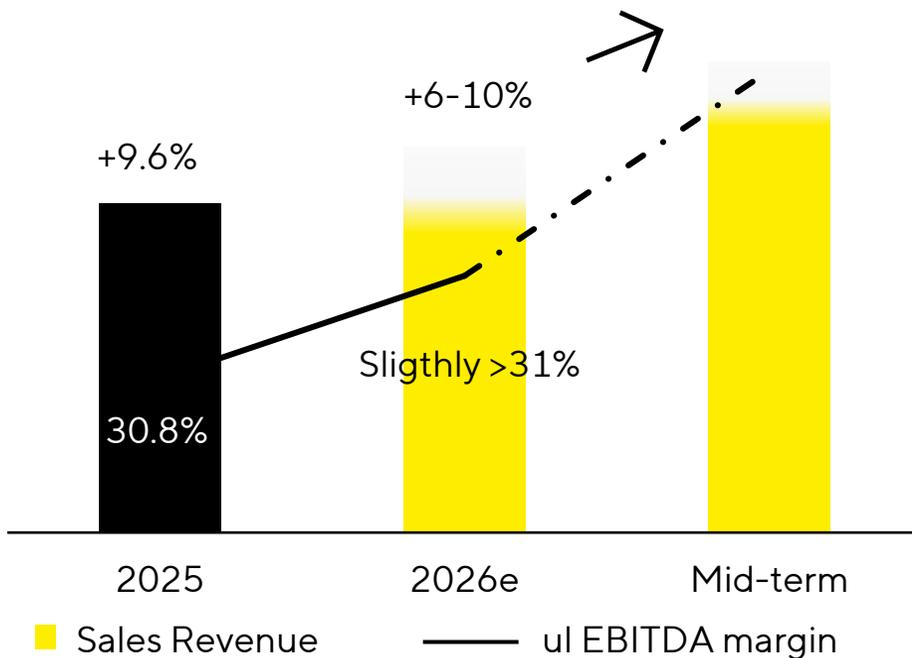


AI and digital capabilities as the foundation for next-level operational excellence

Delivering an attractive
financial profile

2026 as a transition year towards mid-term growth ambition

Sales revenue growth¹, ul. EBITDA margin



¹ in constant currencies

What we currently see



- Market is showing signs of gradual improvement
- Equipment business stabilizing, becoming less dilutive to growth
- China stable to slightly improving, forming a foundation for healthier growth post-2026
- Customer sentiment gradually improving

Sustained outperformance supported by market trends

	Addressable market CAGR mid-term ¹	SSB key drivers of outperformance	SSB mid-term ambition ²			
			Above-market growth p.a.	Organic sales revenue growth p.a. in cc	ul EBITDA margin expansion p.a.	
Sartorius Stedim Biotech	8-10%	<ul style="list-style-type: none"> Leadership in SU and process intensification Unique advanced therapies portfolio 	+100-200bps	>	9-12%	~60-85bps

SSB = Sartorius Stedim Biotech 1 Sartorius market intelligence 2 starting as of 2027

Long-term margin upside embedded in the business model

	Share of cost positions		Expected contribution to margin development	
	2019	2025	2026e	Mid-term
Sales	100%	100%		
Cost of sales	-47.7%	-50.8%	↗	↑
S&D	-16.1%	-15.2%	↗	↗
R&D	-5.5%	-4.4%	→	→
Overhead	-5.3%	-5.3%	↗	↗
Depr. / other	+3.8%	+6.4%	→	→
ul. EBITDA margin	29.3%	30.8%	slightly >31%	+60-85 bps p.a.

Cost positions excluding extraordinary items and amortization

Mid-term key assumptions

- Increasing capacity utilization; growing consumables contribution
- Cost improvements driven by increasing utilization, automation, use of AI
- Operating leverage and fixed cost degression

Bottom-line

- Double-digit ul. EPS growth (p.a.)

↑	↗	→
Positive	Slightly positive	Neutral

Higher efficiency through continuous improvement mindset and scalable growth platforms

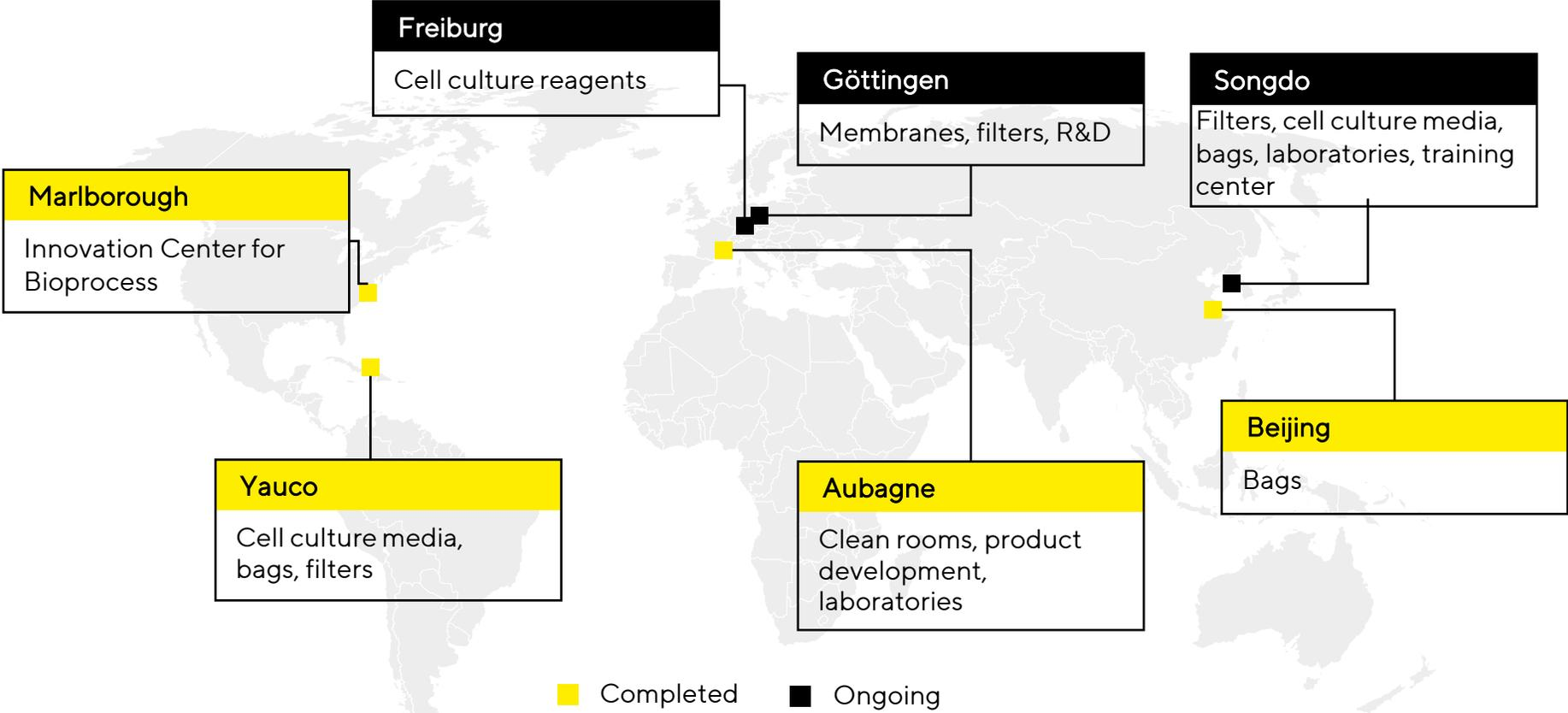
 Efficiency program 2024 as a base	Cost savings of ~€85m with positive margin effects in 2025
 Continuous efficiency initiatives	based on Lean/CI thinking to improve cost structures from year to year
 Powerful digital tools	across commercial, operational and people processes enable efficiency gains
 Under-proportional cost growth	further fueled by operating leverage of infrastructure

Substantial investments support resilient footprint and future growth

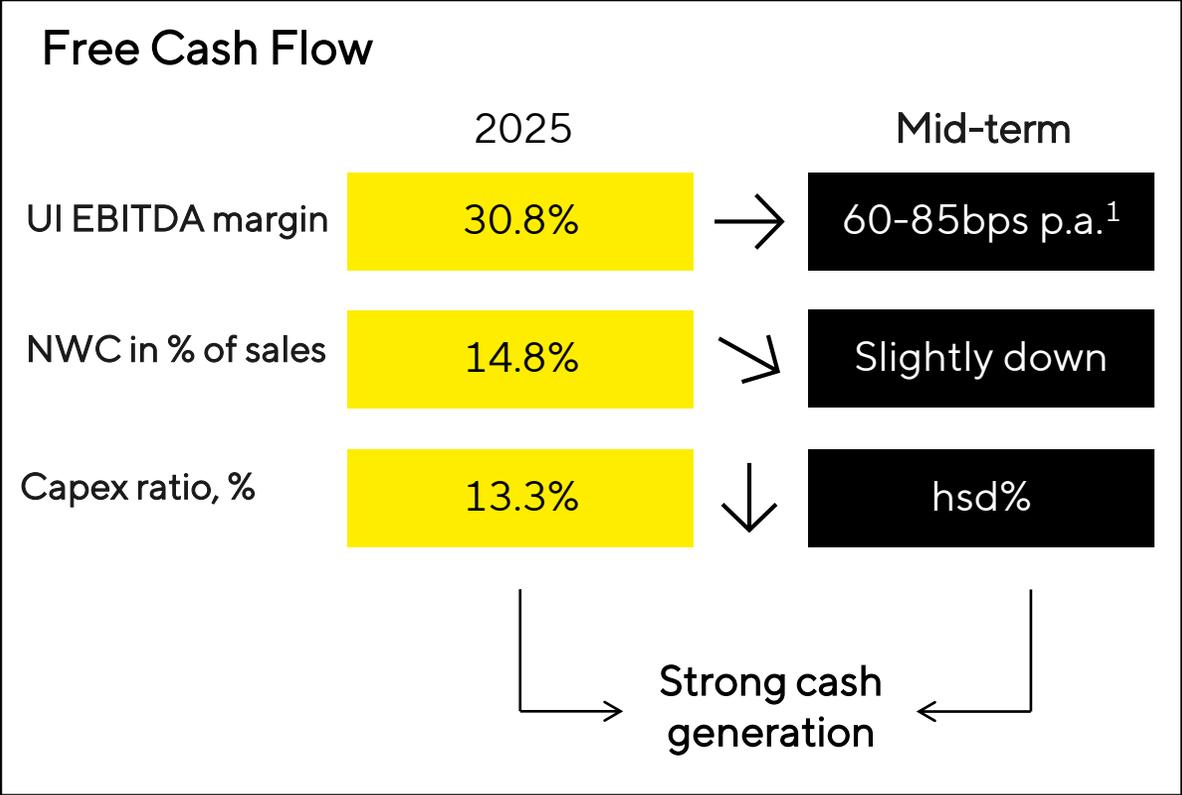
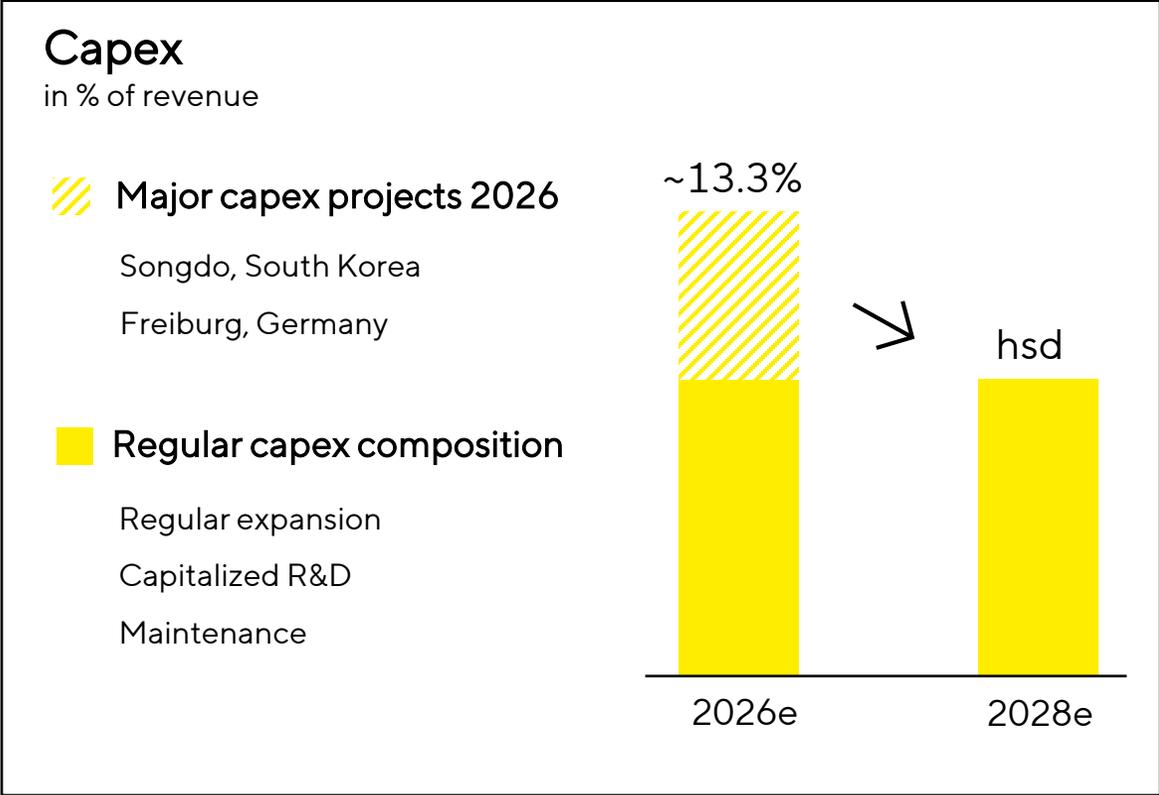
Key Focus Areas

-  Global resilience
-  Redundant capacities
-  Operational excellence
-  Customer proximity

~€2.1bn
Capex in 2020-2025

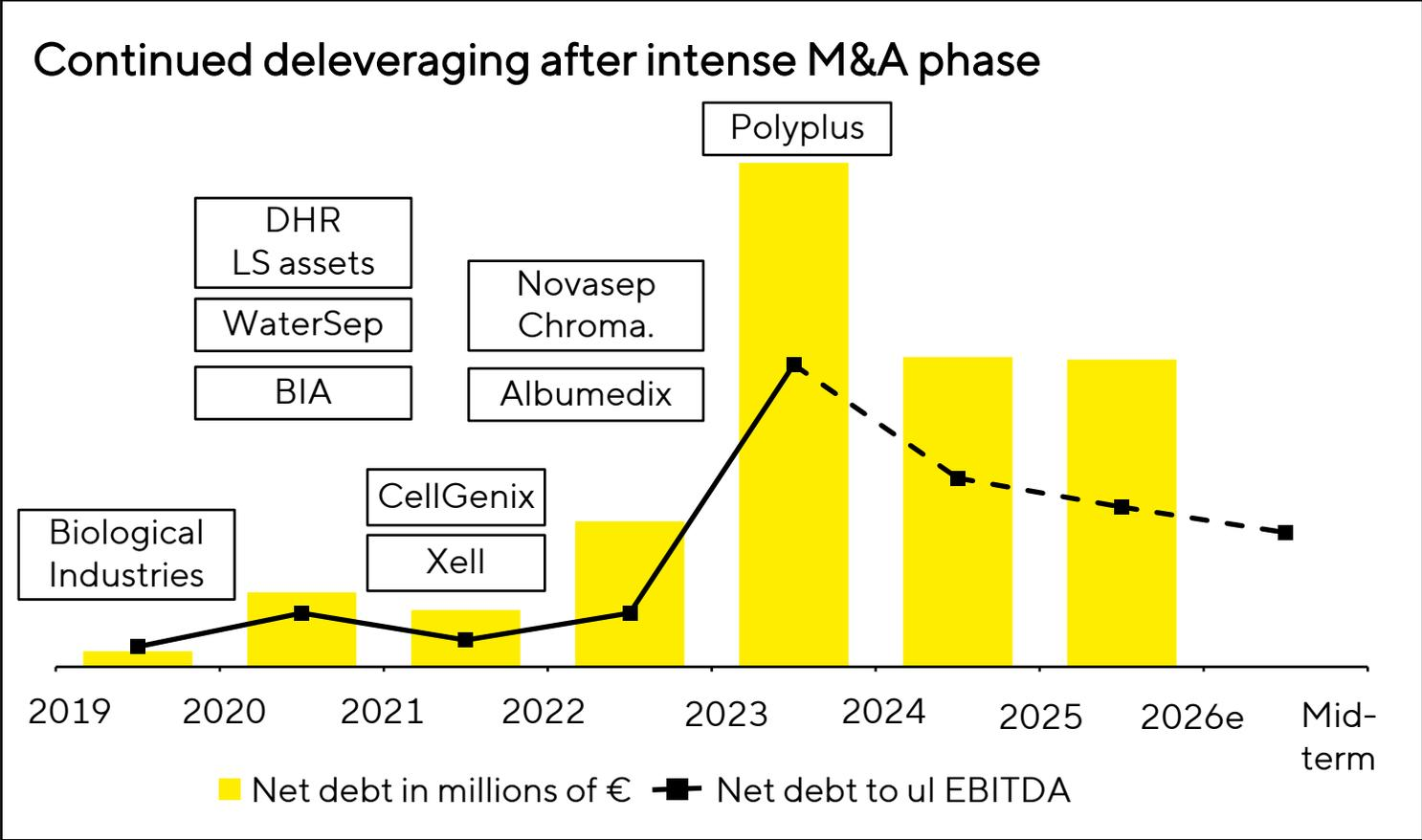


Strong cash generation and financial flexibility



¹ starting as of 2027

Strong focus on deleveraging



- ### Disciplined capital deployment
- Organic deleveraging the top priority
 - Capex discipline and focused on essential growth and productivity projects
 - M&A selective; bolt-on possible
 - Dividends broadly in line with payout ratio in previous years

Sustainable value creation remains well on track

Proven strengths that continue to matter

- Addressing attractive biopharma end markets with high entry barriers and strong structural growth drivers
- Biologics pure-play with a focus on single-use, uniquely positioned to capture above-market growth
- Strong focus on innovation leadership in relevant applications
- Sticky business model with a high share of recurring revenues and significant operating leverage potential

Strategic focus and disciplined execution to drive future growth and returns

- Roadmap to leverage leadership in the core business while expanding future growth platforms
- Targeted initiatives to achieve best-in-class customer experience and operational excellence
- Remain focused on delivering above-market growth in the mid-term based on strong market fundamentals, alongside continuous margin expansion
- Capex normalization and improving cash generation, supporting ongoing deleveraging

Contacts and Financial calendar



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



March 26, 2026

Annual General Meeting

April 23, 2026

Publication Q1 Results 2026

May 19, 2026

RBC Global Healthcare Conference, New York

May 26, 2026

dbAccess European Champions Conference, Frankfurt

June 3, 2026

BNPP Exane CEO Conference, Paris

June 18, 2026

J.P. Morgan Healthcare Forum, London

June 24, 2026

Jefferies German & Swiss Corporate Conference, Baden-Baden

July 23, 2026

Publication H1 Results 2026

October 22, 2026

Publication 9M Results 2026

Simplifying Progress

SARTORIUS