

# SARTORIUS

## Simplifying Progress

### Investor Conference Call Q1 2026 Results

Michael Grosse, Florian Funck, René Fáber  
Sartorius | Sartorius Stedim Biotech | April 23, 2026



# Agenda

Sartorius Group  
Q1 2026 results | FY 2026 outlook

Sartorius Stedim Biotech Group  
Q1 2026 results | FY 2026 outlook

Questions & Answers



# Q1 marks positive start to 2026

## Q1 2026 performance

### Group

- Substantial sales revenue growth; robust growth contribution of consumables in both divisions
- Resilient ul EBITDA margin despite tariff headwinds and investments in future growth initiatives
- Strong free cash flow development; continued deleveraging

### Bioprocess Solutions

- Revenue up ~8%<sup>1</sup>, reflecting strength of recurring business
- Soft start of equipment business as expected; positive outlook for remainder of the year

### Lab Products & Services

- Revenue up ~5%<sup>1</sup>; continued positive development since H2 2025
- Strong consumables and positive bioanalytics business contributing to growth

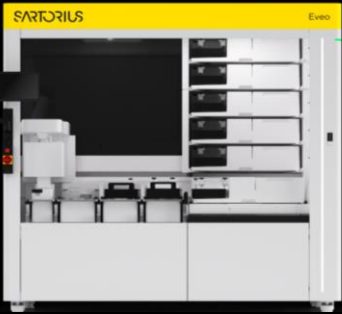
## 2026 guidance confirmed

- Group sales revenue growth<sup>1</sup> of ~5% to ~9%
- Ul. EBITDA slightly above 30%

<sup>1</sup> Constant currencies

# High customer interest underscores relevance of Sartorius' innovations across biologics development and manufacturing

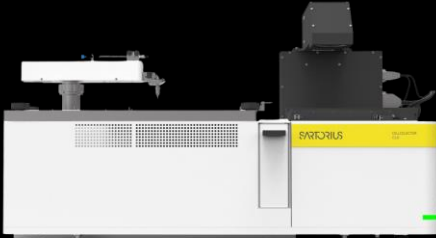
## More efficient cell therapy manufacturing



Evoe Cell Therapy Platform

Automated, multi-parallel production for up to 4x higher yield

## More efficient cell line development



CellCelector CLD

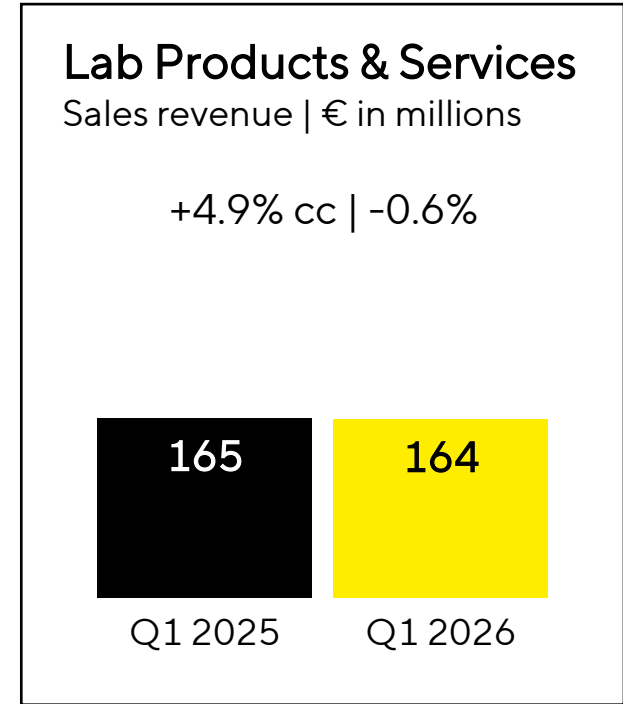
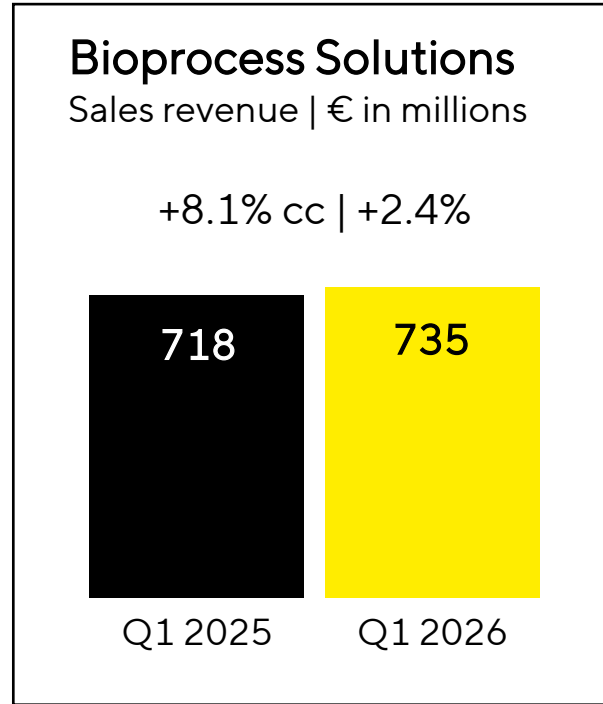
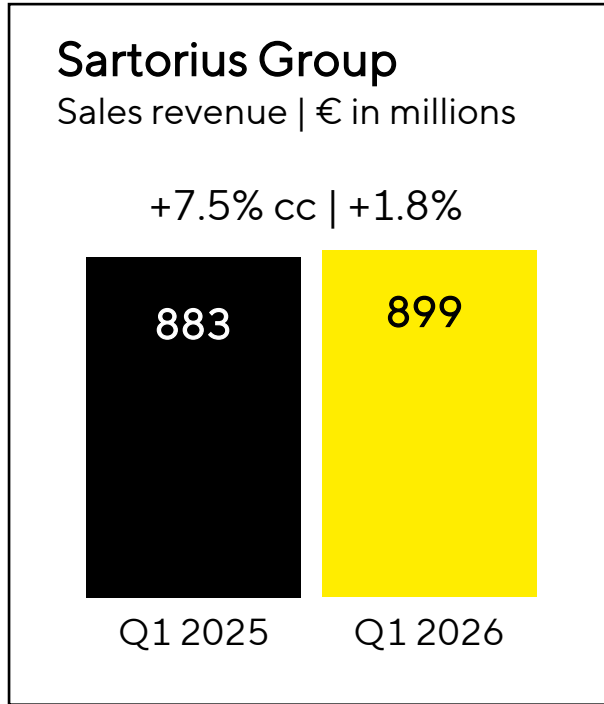
Automated clone selection and advanced imaging



Genetically engineered CHO host cell line

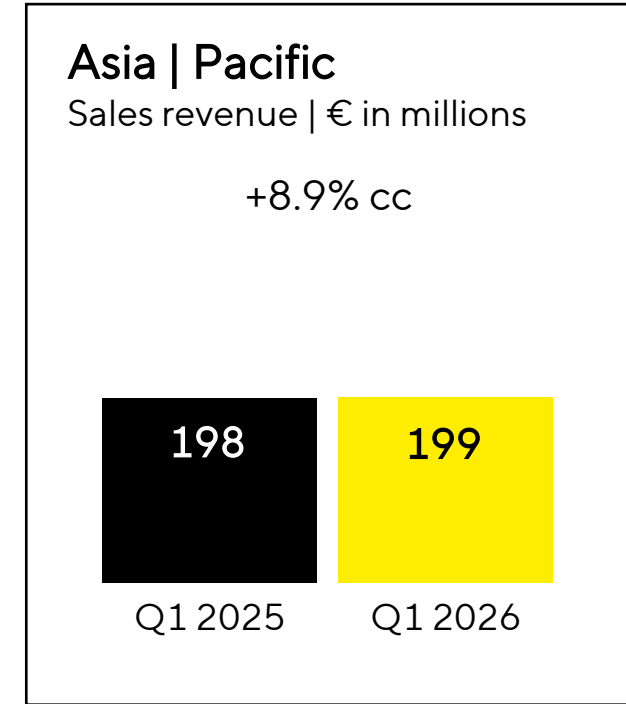
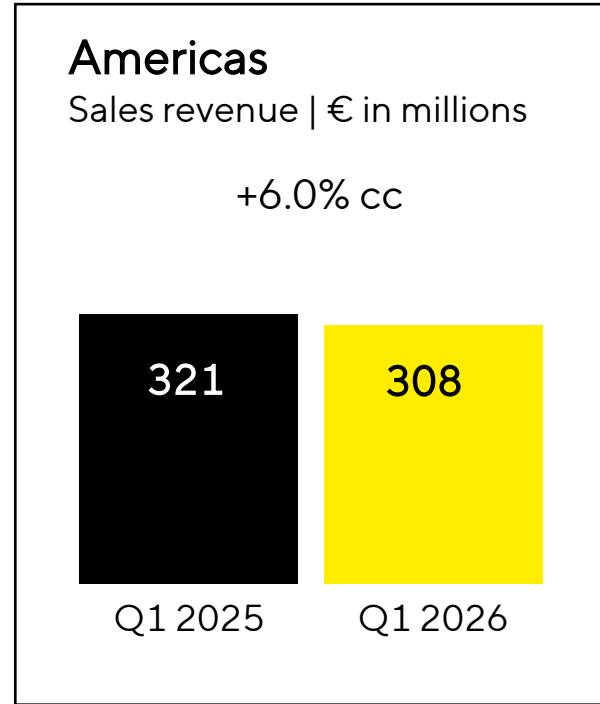
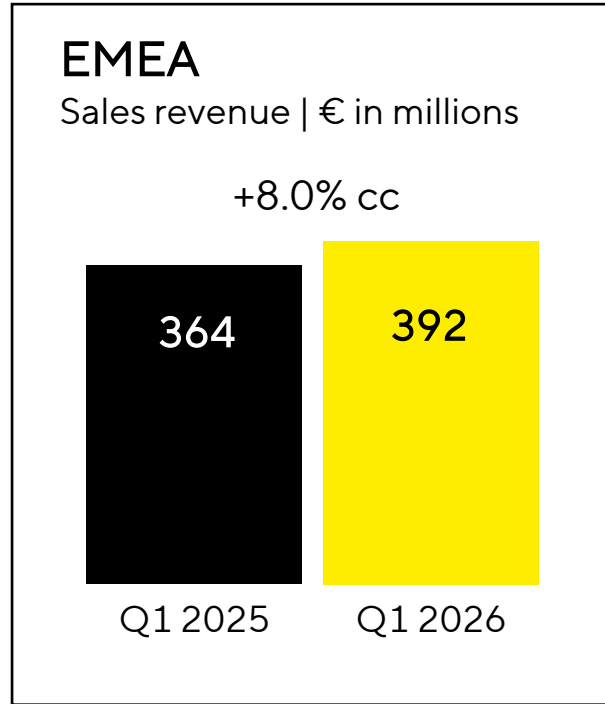
Faster clone development and up to 3x higher productivity

# Substantial sales growth driven by continued consumables strength



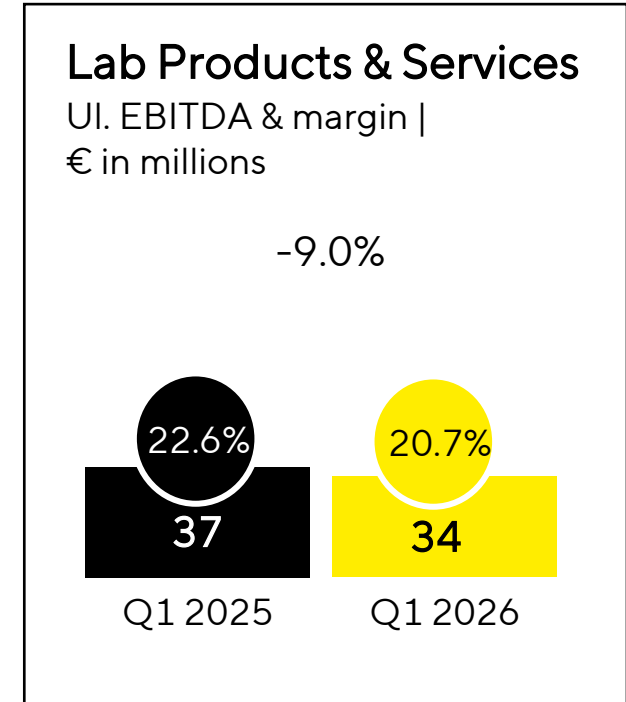
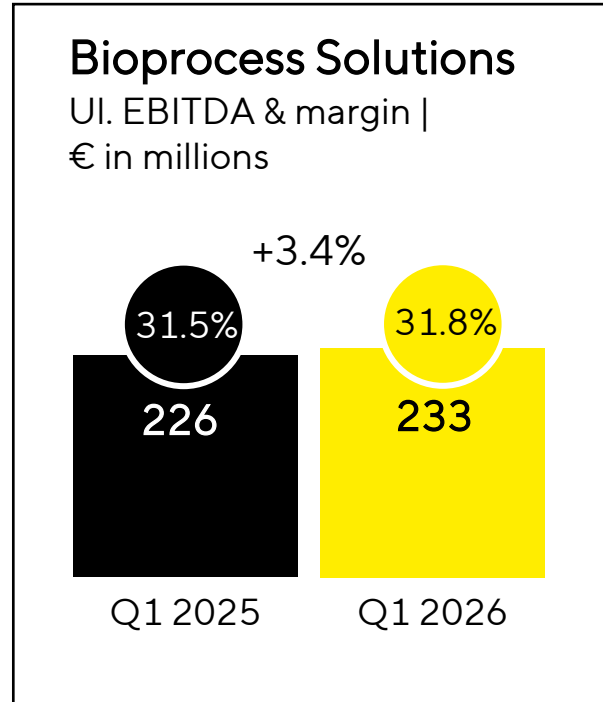
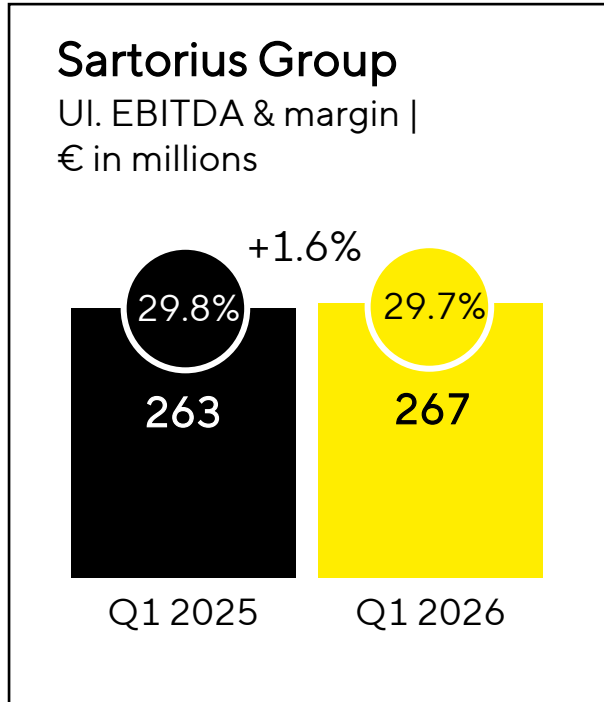
- Continued strong recurring business
- BPS equipment business showed expected soft start to the year; positive momentum of bioanalytics business in LPS
- MATTEK contributes ~2.8pp to LPS growth

# All regions contribute positively to business development



- BPS: Strong growth in APAC and EMEA; Americas with solid growth against a high PY comps
- LPS: Substantial expansion in the Americas against low PY comps; solid expansion in APAC while development in EMEA modest

# Continued solid profitability



- Positive volume effects and operating leverage partly offset by product mix effects and tariff impacts; LPS additionally impacted by investments in future growth initiatives

# Sound cash flow development

Sartorius Group € in millions	Q1 2025	Q1 2026	in %
Underlying net profit <sup>1,2</sup>	85	<b>83</b>	-1.9
Underlying EPS (ord.)	1.22	<b>1.20</b>	-2.0
Underlying EPS (pref.)	1.23	<b>1.21</b>	-1.9
Reported net profit <sup>2</sup>	48	<b>56</b>	+16.0
Operating cash flow	139	<b>189</b>	+35.8
Free cash flow	61	<b>113</b>	+84.6
Capex ratio (in %)	8.6	<b>8.6</b>	+0.0pp

- Underlying net profit reflects higher depreciation for capacity expansion
- Reported net profit increased strongly, driven by FX effects and lower extraordinaries
- Strong increase in operating cash flow, supported by higher EBITDA and lower tax payments compensating growth-related rise in working capital.
- Free cash flow up significantly, driven by higher operating cash flow and stable capex at prior-year level, in line with plan to support future growth

1 Underlying net profit = net profit adjusted for extraordinary items, amortization and based on a normalized financial result and a normalized tax rate 2 After non-controlling interest

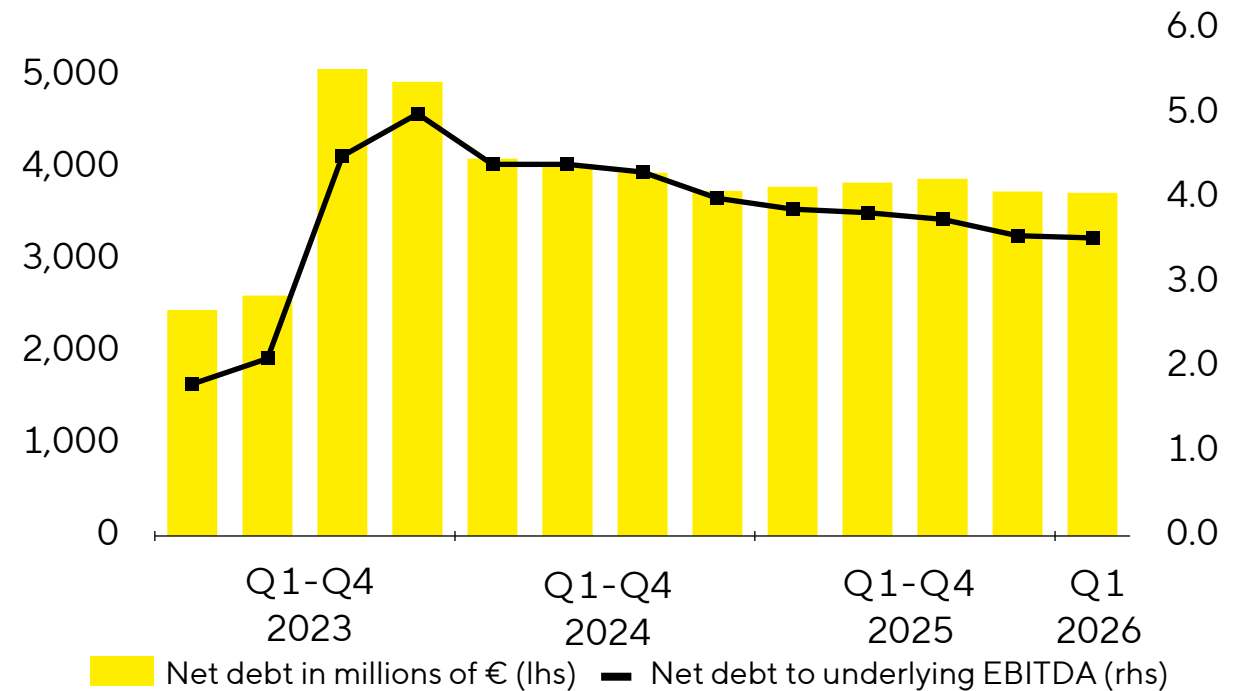
# Deleveraging ongoing; commitment to investment-grade rating unchanged

## Key financial indicators

Sartorius Group € in millions unless otherwise specified	Dec. 31, 2025	Mar. 31, 2026
Equity ratio in %	39.8	<b>39.4</b>
Net debt	3,741	<b>3,727</b>
Net debt   underlying EBITDA	3.55	<b>3.53<sup>1</sup></b>

<sup>1</sup> Includes underlying pro forma EBITDA of acquisitions completed during the last 12 months

## Net debt and net debt to underlying EBITDA



# 2026 guidance confirmed

Guidance 2026	Sartorius Group	Bioprocess Solutions	Lab Products & Services
Sales revenue growth <sup>1</sup>	~5% to ~9%	~6% to ~10%	~2% to ~6%
Underlying EBITDA margin	Slightly above 30%	Slightly above 32%	Slightly below 21%

- Group sales growth includes ~1pp contribution from MATTEK acquisition and US tariff-related surcharges; LPS revenue growth includes ~1.5pp contribution from MATTEK
- Capex ratio at ~PY level
- Net debt/underlying EBITDA to decrease to slightly above 3x

#### For financial modelling on Group level

Depreciation	260mn to 270mn €
Normalized financial result	-170mn to -180mn €
FX effect on reported vs. cc growth	~ -2pp

Due to the continued high dynamics and volatility across the life science industry, the forecast remains subject to greater uncertainty, which is reflected in the current guidance range. Potential changes in U.S. tariffs are likewise not included.

<sup>1</sup> In constant currencies

# Agenda

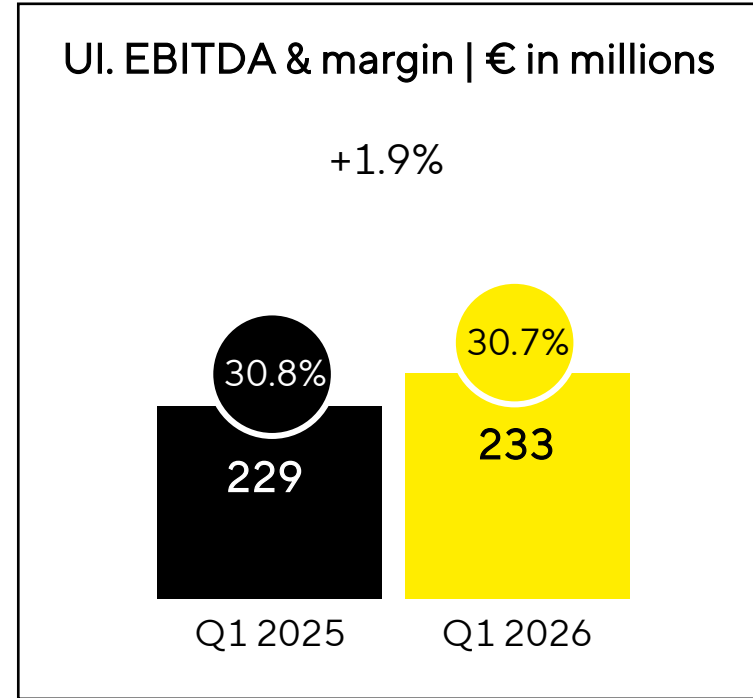
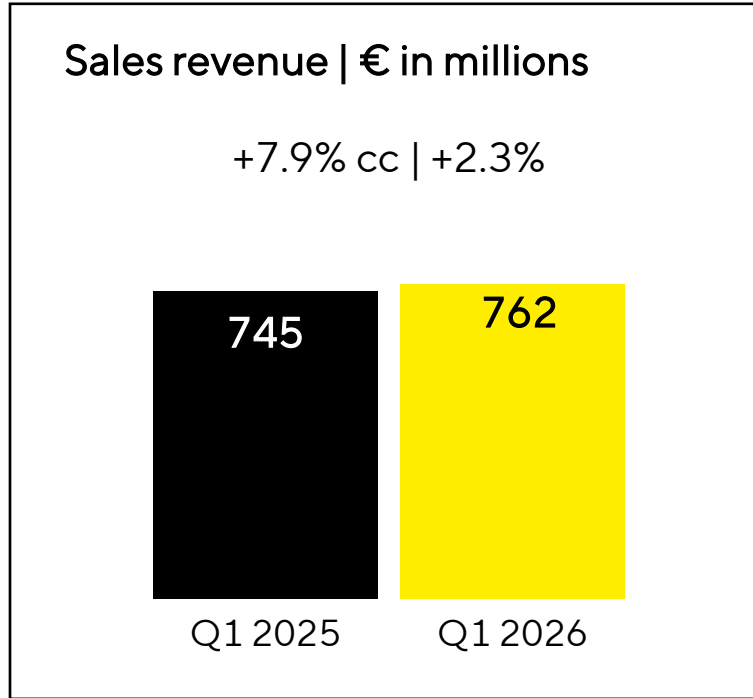
Sartorius Group  
Q1 2026 results | FY 2026 outlook

Sartorius Stedim Biotech Group  
Q1 2026 results | FY 2026 outlook

Questions & Answers

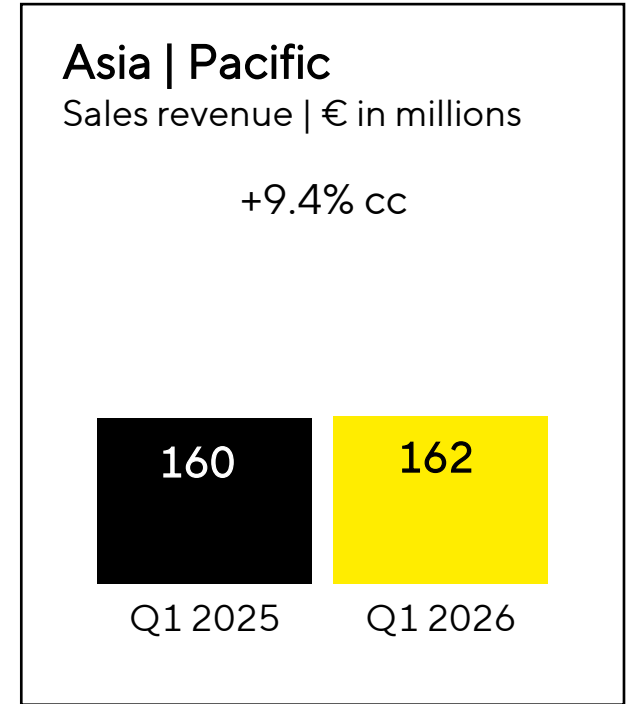
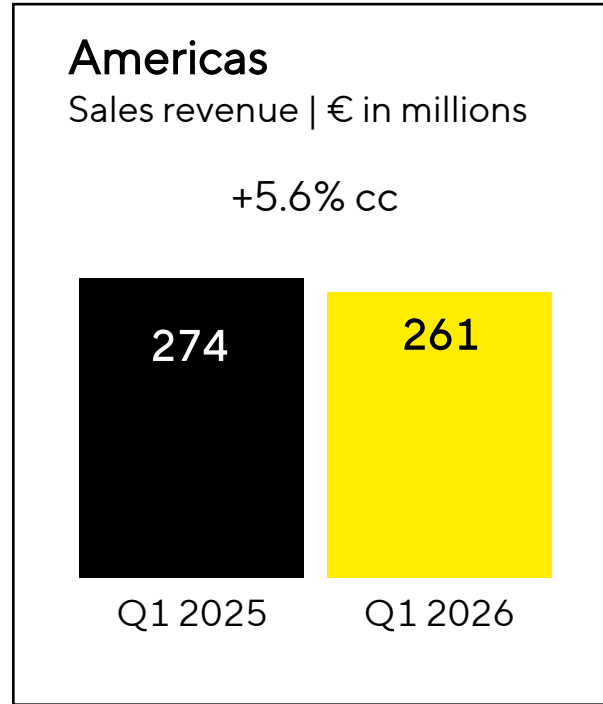
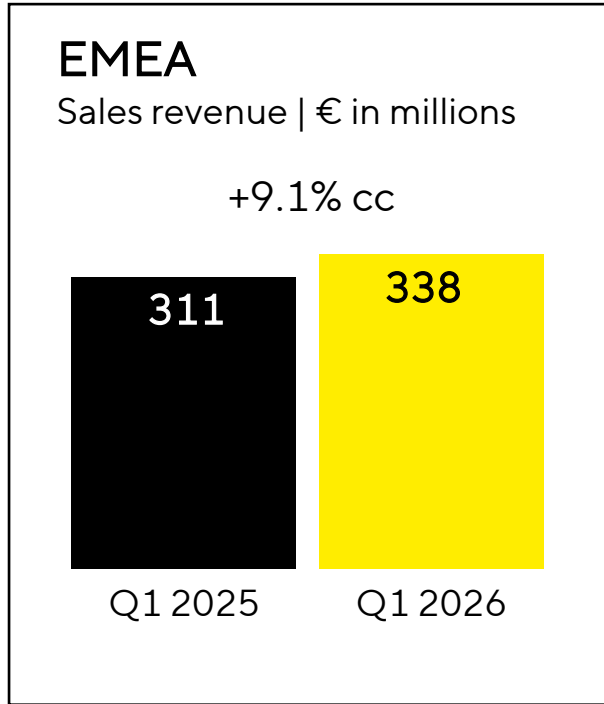


# Substantial sales growth and resilient margin



- Continued strong recurring business; equipment business showed expected soft start to the year
- Positive volume effects and operating leverage offset by product mix effects and tariff impact

# All regions contribute positively to business development



- Strong growth in APAC and EMEA; Americas with solid growth against a high PY comps

# Sound cash flow development

Sartorius Stedim Biotech € in millions	Q1 2025	Q1 2026	in %
Underlying net profit <sup>1,2</sup>	113	<b>114</b>	+0.5
Underlying EPS	1.16	<b>1.17</b>	+0.5
Reported net profit <sup>2</sup>	86	<b>88</b>	+3.0
Operating cash flow	120	<b>193</b>	+61.4
Free cash flow	55	<b>124</b>	+126.9
Capex ratio (in %)	8.8	<b>9.1</b>	+0.3pp

- Underlying net profit reflects higher depreciation for capacity expansion
- Reported net profit up, driven by FX effects and lower extraordinaries
- Strong operating cash flow, supported by higher EBITDA and lower tax payments compensating growth-related increase in working capital.
- Free cash flow rose significantly, driven by higher operating cash flow and capex slightly above prior-year level, in line with plan to support future growth

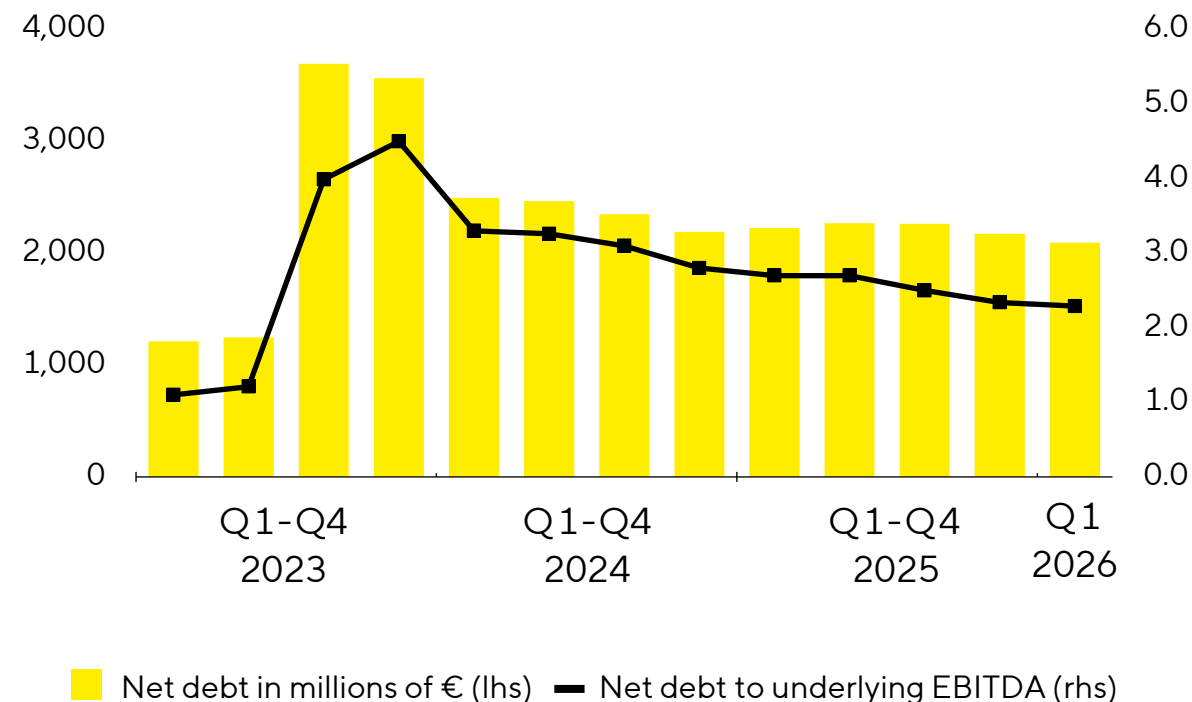
1 Underlying net profit = net profit adjusted for extraordinary items, amortization and based on a normalized financial result and a normalized tax rate 2 After non-controlling interest

# Strong balance sheet; deleveraging ongoing

## Key financial indicators

Sartorius Stedim Biotech € in millions unless otherwise specified	Dec. 31, 2025	Mar. 31, 2026
Equity ratio in %	51.7	50.6
Net debt	2,173	2,093
Net debt   underlying EBITDA	2.38	2.28

## Net debt and net debt to underlying EBITDA



# 2026 guidance confirmed

Guidance 2026	Sales revenue growth <sup>1</sup>	Underlying EBITDA margin
Sartorius Stedim Biotech Group	~6% to ~10%	Slightly above 31%

- Sales growth includes ~1pp contribution from US tariff-related surcharges
- Capex ratio at ~PY level
- Net debt/ul. EBITDA to decrease to slightly above 2x

For financial modelling on SSB level	
Depreciation	200mn to 210mn €
Normalized financial result	-105mn to -115mn €
FX effect on reported vs. cc growth	~ -2pp

Due to the continued high dynamics and volatility across the life science industry, the forecast remains subject to greater uncertainty, which is reflected in the current guidance range. Potential changes in U.S. tariffs are likewise not included.

<sup>1</sup> In constant currencies

# Questions & Answers



# Contacts and Financial calendar

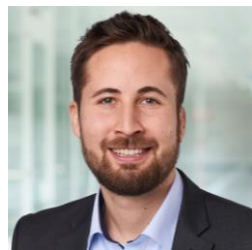


**Petra Müller**

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



**May 19, 2026**

RBC Global Healthcare Conference, New York

**May 21, 2026**

Berenberg Roadshow, Luxembourg

**May 26, 2026**

dbAccess European Champions Conference, Frankfurt

**May 28, 2026**

Jefferies Roadshow, Stockholm

**June 3, 2026**

BNPP Exane CEO Conference, Paris

**June 18, 2026**

J.P. Morgan Healthcare Forum, London

**June 24, 2026**

Jefferies German & Swiss Corp. Conference, Baden-Baden

**July 23, 2026**

Publication H1 Results 2026

**October 22, 2026**

Publication 9M Results 2026

# Disclaimer

---

This presentation contains statements concerning the future performance of the Sartorius and the Sartorius Stedim Biotech Groups. 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 substantially diverging 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.

---

# Simplifying Progress

**SARTORIUS**