



Sartorius and Sartorius Stedim Biotech Capital Markets Day

Our Capital Markets Day will offer you insights in the latest business developments and future plans to continue Sartorius' and Sartorius Stedim Biotech's journey of profitable growth.



Agenda | March 17, 2026

07:30 – 09:00 Registration for on-site participants

08:00 – 08:50 Site tours

Strategy Session incl. Q&A

09:00 – 11:20 Strategy update & Financial perspectives

Executive Board

Webcasted

11:20 – 11:35 Coffee break

Spotlight Sessions incl. Q&A

11:35 – 12:05 Leveraging AI to drive product innovation and operational excellence

CTO & Operations Management

12:05 – 12:45 Lunch break

12:50 – 13:25 Pioneering the future of Single-Use bioprocessing

Bioprocessing Management

13:25 – 14:00 Empowering novel modalities through innovation

Advanced Therapies Solutions Management

14:00 – 14:30 Advanced cell models for faster and smarter drug discovery

Advanced Cell Models Management

14:30 Closing remarks

CEO