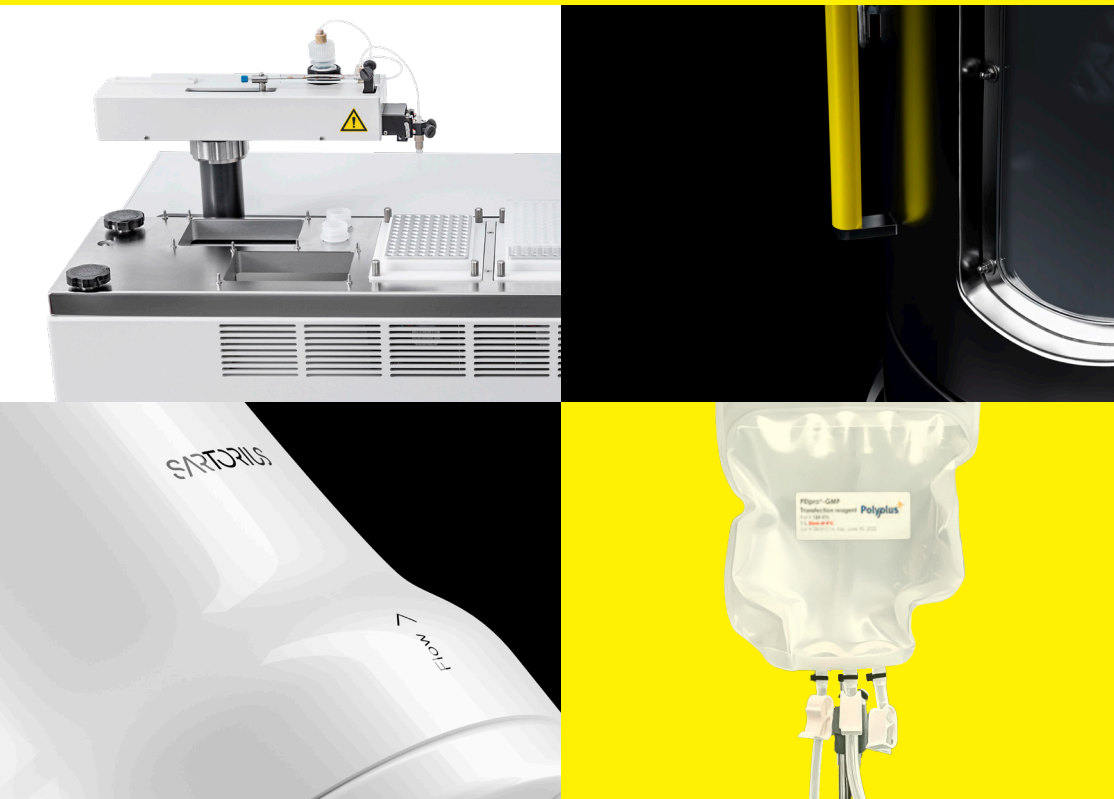


Sartorius and Sartorius Stedim Biotech Capital Markets Day

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. Extensive Q&A sessions will complement the presentations.



Agenda | May 16, 2024

08:00 – 10:00 Registration for on-site participants

08:30 – 09:30 Site tour

09:55 – 10:00 Opening remarks
Dr. Joachim Kreuzburg, CEO

Webcasted

10:00 – 10:40 Group vision & strategy
Dr. Joachim Kreuzburg, CEO

Webcasted

Q&A session

10:40 – 11:20 Financial perspective
Dr. Florian Funck, CFO

Webcasted

Q&A session

11:20 – 11:40 Coffee break

11:40 – 12:40 Deep dive Lab Products & Services Division
Dr. Alexandra Gatzemeyer, Head of LPS Division
& LPS leadership team

Q&A session

12:40 – 13:30 Lunch break

13:30 – 14:30 Deep dive Bioprocess Solutions Division | Sartorius Stedim Biotech
Dr. René Fáber, Head of BPS Division & CEO SSB
& SSB leadership team

Q&A session

14:30 – 15:00 Digital tools in biopharma research & production
Dr. Oscar-Werner Reif, CTO

Q&A session

15:00 Closing remarks
Dr. Joachim Kreuzburg, CEO