

SARTORIUS

Sartorius Data Analytics Summit

Turning Data into Value-Driving Decisions

October 20-22, 2026
Aubagne, France



Key Topics for the Summit

DOE and MVDA:

Translating structured data analysis into usable process knowledge.

AI-enabled decision-making:

Turning analytics outputs into confident process decisions.

PAT and process monitoring:

Connecting real-time data with process insight and control.

Digital twins:

Applying predictive models to support scale-up, optimization, and troubleshooting.

Machine learning:

Revealing relationships in process data to guide faster development.

Event landing page:

[Sartorius Data Analytics Summit](#)

What You Don't Want to Miss

Hands-on Experience

Participate in our pre-conference workshops and gather practical strategies to turn data into measurable impact.

Expert Guidance

Engage in product demos and guided tours of our Application Centre to see innovations in action.

Stay Competitive

Learn how to scale analytics across your organization.

Networking with Peers

Connect with industry leaders and fellow professionals to exchange ideas and experiences.

Industry Expertise

Learn directly from real customer success stories and expert-led sessions.

Unforgettable Atmosphere

Visit our exhibition, get inspired and enjoy delicious food, including our two evening events.

Event landing page:

[Sartorius Data Analytics Summit](#)

Preliminary Agenda | Pre-Conference Workshops

Tuesday, October 20

| Start | End | Topic |
|-------|-------|---|
| 8:30 | 9:30 | Coffee Registration |
| 9:30 | 11:00 | Workshops: <ul style="list-style-type: none">▪ BUILD Create models and establish workflows while exploring machine learning and Design of Experiments.▪ INNOVATE Use data and digital twins for rapid prototyping, simulations and AI exploration. |
| 11:00 | 11:30 | Coffee Break |
| 11:30 | 13:00 | Workshops: <ul style="list-style-type: none">▪ DEPLOY Move your models into GxP environments for automated real-time process control.▪ ELEVATE Advance your analytics strategy with tools and practices that boost performance across your bioprocess. |
| 13:00 | 14:00 | Lunch Break |
| 14:00 | 15:30 | Workshops: <ul style="list-style-type: none">▪ BUILD Create models and establish workflows while exploring machine learning and Design of Experiments.▪ INNOVATE Use data and digital twins for rapid prototyping, simulations and AI exploration. |
| 15:30 | 16:00 | Coffee Break |
| 16:00 | 17:30 | Workshops: <ul style="list-style-type: none">▪ DEPLOY Move your models into GxP environments for automated real-time process control.▪ ELEVATE Advance your analytics strategy with tools and practices that boost performance across your bioprocess. |
| 18:30 | 22:00 | Welcome Reception |

Preliminary Agenda | Day 1

Wednesday, October 21

| Start | End | Topic | Speaker |
|-------|-------|--|--|
| 8:30 | 9:00 | Coffee Registration | |
| 9:00 | 9:15 | Welcome | Barry Heavey, Head of Data Analytics at Sartorius |
| 9:15 | 9:45 | Opening Keynote | Michael Grosse, CEO of Sartorius |
| 9:45 | 10:15 | Speaker Presentation: Upstream Modelling & Downstream Application | TBA |
| 10:15 | 10:45 | Speaker Presentation: Data to Insight: AI, Machine Learning and PAT to Accelerate Bioprocess Development | Stephen Goldrick, Lecturer in Bioprocess Digitalization at University College London |
| 10:45 | 11:15 | Coffee Break | |
| 11:15 | 11:45 | Speaker Presentation: Hybrid Modelling in Cell Line Development and Process Optimization | Ali Safari, Sr. Data Scientist at Sartorius |
| 11:45 | 12:15 | Speaker Presentation: Use Cases of Data Analytics in Development | |
| 12:15 | 13:00 | Speaker Presentations: Top 10 Reasons Why You're Stuck in Pilot Purgatory, preventing ROAI & Economies of Scale from your AI & Transformation Initiatives | Jonathan Alexander, Global Manufacturing AI & Advanced Analytics Manager at Albemarle |
| 13:00 | 14:00 | Lunch Break Exhibition | |
| 14:00 | 14:40 | Speaker Presentation: From Powder to Tablet: A Modeling Framework for Continuous Manufacturing | Maria Cerrato Oliveros, Senior Manager at Johnson & Johnson Innovative Medicines |
| 14:40 | 15:00 | Regulatory Compliance Session | |
| 15:00 | 16:00 | Product Updates: How can the Umetrics® Ecosystem support Actionable, Confident Decision Making | |
| 16:00 | 16:30 | Coffee Break | |
| 16:30 | 17:00 | Speaker Presentation: Data Analytics in Manufacturing | |
| 17:00 | 18:00 | Meet the Experts Exhibition | |
| 18:30 | 22:00 | Networking Dinner | |

Preliminary Agenda | Day 2

Thursday, October 22

| Start | End | Topic | Speaker |
|-------|-------|---|---|
| 8:30 | 9:00 | Coffee Registration | |
| 9:00 | 9:15 | Welcome Day 2 | |
| 9:15 | 10:00 | Speaker Presentation | Michael Mølmann, CEO of MugAln Group |
| 10:00 | 10:30 | Speaker Presentation: User Adoption of SIMCA®-online | TBA |
| 10:30 | 11:00 | Coffee Break | |
| | | Data Analytics Roadmap | |
| 11:00 | 11:45 | <ul style="list-style-type: none">▪ How the Umetrics® Ecosystem helps transform fragmented, high-dimensional data into structured, actionable insights—enabling faster, more confident decision-making.▪ Learn how Sartorius supports the deployment and scaling of analytics and digital twins into real-world operations, driving measurable impact across development, manufacturing, and the enterprise. | |
| 11:45 | 12:30 | Closing Keynote | TBA |
| 12:30 | 12:45 | Thank You | Barry Heavey, Head of Data Analytics at Sartorius |
| 12:45 | 13:45 | Lunch | |
| 14:00 | 14:30 | Application Centre Tour (Optional) Departure | |

SARTORIUS