

SARTORIUS

Sartorius Food Biotech Summit

October 22 – 24, 2025 | Göttingen, Germany



Wednesday, October 2

Start	Topic
12:30 p.m.	Registration & Snacks
1:00 p.m.	Welcome and Remark
1:15 p.m.	Food Biotech Biomanufacturing: Challenges & End-to-End Solutions
Session 1: Media & Analytics	
1:45 p.m.	CCM & Analytical Tools - Proof of Concept for (cost-effective) Medium
2:15 p.m.	Industry Experts Presentations
3:30 p.m.	Coffee Break & Networking
Session 2: Process & Technologies	
4:00 p.m.	Application Tour (Group 1) Ambr® Demonstration (Group 2) Virtual Reality Experience (Group 3)
4:45 p.m.	Application Tour (Group 2) Ambr® Demonstration (Group 3) Virtual Reality Experience (Group 1)
5:30 p.m.	Clean and Consistent: The Future of Media Preparation
6:00 p.m.	Conclusion of Day 1
6:30 p.m.	Finger Food and Networking
8:30 p.m.	Travel back to Hotel

Sessions

Session 1 Media & Analytics	Session 2 Process & Technologies	Session 3 Innovation & Services	Practical Demonstrations
--------------------------------	-------------------------------------	------------------------------------	--------------------------

Thursday, October 23

Start	Topic
8:30 a.m.	Welcome & Coffee
	Session 2: Process & Technologies
9:00 a.m.	From Cells to Solutions: Exploring Bioreactor Technologies
9:45 a.m.	Industry Experts Presentations
10:15 a.m.	Coffee Break & Networking
11:15 a.m.	Application Tour (Group 3) Ambr® Demonstration (Group 1) Virtual Reality Experience (Group 2)
12:30 p.m.	Lunch
1:30 p.m.	Downstream Processing Solutions
2:30 p.m.	Industry Experts Presentations
3:30 p.m.	Coffee & Snacks
4:00 p.m.	Round Table Discussion
5:00 p.m.	Conclusion & Remarks
5:30 p.m.	Travel back to Hotel
7:30 p.m.	Dinner: AMAVI Güterbahnhofstraße 10 D-37073 Göttingen

Friday, October 24

Start	Topic
8:30 a.m.	Welcome & Coffee
	Session 3: Innovation & Services
9:00 a.m.	Digital Solutions: Bytes to Bites: Software & DOE in Meat Alternatives
9:30 a.m.	Confidence Validation & Testing
10:00 a.m.	Coffee Break & Networking
10:30 a.m.	Industry Experts Presentations
12:00 p.m.	Wrap-up, Closure
12:30 p.m.	Lunch
2:00 p.m.	Travel

Sessions

Session 1 Media & Analytics	Session 2 Process & Technologies	Session 3 Innovation & Services	Practical Demonstrations
--------------------------------	-------------------------------------	------------------------------------	--------------------------