

Data Analytics User Meeting Conference Agenda

Wednesday, September 21

Start	End	Topic	Speaker
8:00 a.m.	8:25 a.m.	Registration and Coffee	
8:30 a.m.	8:55 a.m.	Welcome	Data Analytics, Sartorius
9:00 a.m.	9:20 a.m.	Creating a New Paradigm in Process Monitoring within a Drug Substance Manufacturing	Anthony Maguire, Consultant Statistician, Eli Lilly
9:25 a.m.	10:05 a.m.	Customer Keynote The Advanced Analytics Endgame: Deploying Data Science Programs to Maximize Value Streams in Biotech Manufacturing	Will Penland, Head of Data Analytics, Global Manufacturing Science and Technology, Boehringer Ingelheim
10:10 a.m.	10:40 a.m.	Coffee	
10:45 a.m.	11:25 a.m.	Customer Keynote Improving Equipment Reliability Through Condition-Based Maintenance	Chris Garvin, Director Data Sciences, Amgen
11:30 a.m.	11:50 a.m.	Enabling End to End Technology Transfer and Process Improvements Through Digitalization Efforts During a Pandemic	Gwyneth Lando, Data Scientist in Global Advanced Analytics, Janssen
11:55 a.m.	12:15 p.m.	Real-Time MVDA Monitoring of Biopharmaceutical Manufacturing Process With SIMCA®-online	Baptiste Duhem, Digital Transformation Lead, Takeda
12:20 p.m.	12:40 p.m.	Process Monitoring and Troubleshooting at Sanofi (Vaccines) Toronto	Ramila Peiris, Head of Data Science, Manufacturing Technology and Michael Celarek, Data Scientist, Process Specialist, Sanofi
12:45 p.m.	1:30 p.m.	Lunch	
1:35 p.m.	3:30 p.m.	Umetrics® Products – The Next Generation	Jonas Elfving, Product Manager and Jon Gabrielsson, Product Manager, Sartorius
3:35 p.m.	5:30 p.m.	Meet the Experts	Open Exhibition
6:00 p.m.	Open End	Excursion and Dinner	

Thursday, September 22

Start	End	Topic	Speaker
8:30 a.m.	9:10 a.m.	Downstream Chromatography and BIA Separation	Chris McCready, Principle Research Scientist, Sartorius
9:15 a.m.	9:35 a.m.	How to Get Your User to Use What You Make	Kelly Richards, Data Scientist, Lonza
9:40 a.m.	10:00 a.m.	Combining Multiple Bioprocesses to Create Generalized SIMCA®-online Batch Monitoring Models	Alex Delafontaine, Process Sciences Associate, Regeneron
10:00 a.m.	10:25 a.m.	Coffee	
10:30 a.m.	10:50 a.m.	Combining Software From Umetrics® Suite With Sartorius Products (BioPAT® MFCS, Chromatography and Ambr®) To Maximize Value for Our Customers	Andrew Hines, Account Manager, Sartorius and Johan Hultman, Manager OEM and Embedded Solutions, Sartorius
10:55 a.m.	11:15 a.m.	Using Multivariate Analysis to Decrease Expense and Carbon Emissions in a Power House Application	Nathan Verhanovitz, Performance Engineer Power and Water Infrastructure Planning and Facilities, Michigan State University
11:20 a.m.	11:40 a.m.	SIMCA®-online – The Journey	Tim Michaelson, Principal Data Scientist, International Paper
11:45 a.m.	12:05 p.m.	Chemometrics Novices Must Be Able to Use Multivariate Models	Mary Beth Seasholtz, Technology Principal in Analytics & Chemometrics, Dow
12:10 p.m.	1:10 p.m.	Lunch	
1:15 p.m.	2:05 p.m.	Partner Sessions <ul style="list-style-type: none"> ▪ In-Process Media Monitoring Using the Bioaccord™ LC-MS and MVDA From Multi-Parallel Ambr® 15 Microbioreactor System ▪ Empowering Process Optimization: Rapid At-Line Amino Acid Monitoring Coupled With Integrated Multivariate Analytics ▪ Emerson and Sartorius Collaborate to Deliver Spectral PAT Capabilities ▪ Cross-Scale Raman Spectroscopy in Bioprocess 	<p>Mark Wrona, Consulting Scientist, Waters Corporation</p> <p>Kerin Gregory, REBEL Product Manager, 908 Devices</p> <p>Michalle Adkins, Director, Life Sciences Consulting, Emerson</p> <p>Nicholas MacLean, Applications Scientist, Endress+Hauser</p>
2:10 p.m.	2:30 p.m.	Albemarle Intelligence – Scaling AI Around Your Global Manufacturing Plants	Jonathan Alexander, Manufacturing AI & Advanced Analytics Manager, Albemarle
2:35 p.m.	3:00 p.m.	Coffee	
3:05 p.m.	3:45 p.m.	Customer Keynote What Life Sciences Companies Can Learn from Big Tech	Dr. Dave Latshaw II, Co-Founder, Biophy
3:50 p.m.	4:15 p.m.	Closing Remark	Data Analytics, Sartorius



For more information, visit

www.sartorius.com/data-analytics