

# Sartorius Day Basel 2020

February 18, 2020 | Gare du Nord, Basel, Switzerland



## Agenda

**09:00 – 09:45**

Open Doors | Coffee  
Bar du Nord

**09:45 – 10:00**

Welcome  
Konzertsaal

### Session 1

**ambr® Upstream**

**10:00 – 10:30**

Introduction: ambr® Upstream  
Karl Scheibenbogen, Application Specialist Fermentation, Sartorius Stedim Biotech GmbH, Goettingen, Germany  
Konzertsaal

**10:30 – 11:00**

ambr® 250 Reduced Scale Model Fit-For-Purpose Evaluation for Process Development and Characterization of Bacterial Fed-Batch Fermentation

Claudia Roder, Team Leader USP Development, Janssen Vaccines, Branch of Cilag GmbH International, Bern, Switzerland  
Konzertsaal

**11:00 – 11:30**

Exhibition | Hands on Session  
Konzertsaal | Bar du Nord

**Session 2**

**ambr® Downstream**

**11:30 – 12:00**

Introduction: ambr® Downstream

Susan Prochaska, Application Specialist Separation, Sartorius Stedim Switzerland AG, Tagelswangen, Switzerland  
Konzertsaal

**12:00 – 12:30**

User Experience With ambr® crossflow at CSL Behring: Automated pH Adjustment During Crossflow Filtration Using ambr®

Fiona Daly, Scientist, R&D | Process Development Group 4, CSL Behring AG, Bern, Switzerland  
Konzertsaal

**12:30 – 13:30**

Lunch  
Bar du Nord

Register now:

[www.sartorius.com/pr/sartorius-day-basel](http://www.sartorius.com/pr/sartorius-day-basel)



## Agenda

### Session 3

#### Option 1: Umetrics

**13:30 – 14:00**

Introduction: Umetrics

Winfried Geis, Application Specialist Umetrics, Sartorius Stedim Biotech GmbH, Goettingen, Germany

Konzertsaal

**14:00 – 14:30**

ambr<sup>®</sup> crossflow and DoE as a Tool for Optimization of EnPCs<sup>®</sup> Purification

Victoria Demina, Senior Scientist, Neuway Pharma GmbH, Bonn, Germany

Konzertsaal

**14:30 – 15:00**

MVDA in Bioprocessing: Selected Case Studies

Timo Schmidberger, Senior Data Scientist, Sartorius Stedim Biotech GmbH, Goettingen Germany

Konzertsaal

#### Option 2: Sartorius Lab Instruments & BioAnalytics

**13:30 – 14:00**

Introduction: Sartorius Lab Instruments & BioAnalytics

Ines Hassad, Sales Representative, Sartorius Stedim Switzerland AG, Tagelswangen, Switzerland

Bar du Nord

**14:00 – 14:45**

IncuCyte<sup>®</sup> Live Cell Analysis Imaging System: Technique and Application Overview for Phenotypic Cell Based Assays

Nicolas Barsuglia, Technical Sales Specialist, Switzerland | France, Essen BioScience Ltd., United Kingdom

Bar du Nord

**15:00 – 15:30**

Coffe Break | Exhibition | Hands on Session

Konzertsaal | Bar du Nord

### Session 4

#### Cellca & BioOutsource

**15:30 – 16:00**

Introduction: Cellca & BioOutsource

Julia Uhlmann, Application Specialist Cellca & BioOutsource, Sartorius Stedim Biotech GmbH, Ulm, Germany

Konzertsaal

**16:00 – 16:30**

Quality and Flexibility to Guarantee Success

Dr. Stefan Warmuth, Head CMC – Therapeutic Program Leader, Numab Therapeutics AG, Wädenswil, Switzerland

Konzertsaal

**16:30 – 16:45**

Closing

Konzertsaal

**16:45 – 19:00**

Apero

Bar du Nord

## Venue

Gare du Nord  
Schwarzwaldallee 200  
4058 Basel, Switzerland  
[www.garedunord.ch](http://www.garedunord.ch)

## Your Contact

Max Günther Müller  
Sartorius Stedim Biotech GmbH  
[Sartorius\\_Day\\_Basel\\_2020@Sartorius.com](mailto:Sartorius_Day_Basel_2020@Sartorius.com)