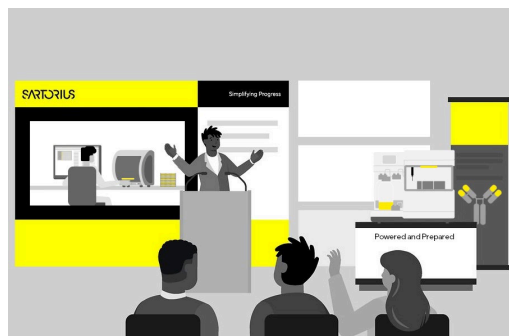


# Sartorius User Group Meeting

Wednesday, October 2<sup>nd</sup>, 2024, in Malmö, Sweden

Register NOW



|       |   |
|-------|---|
| 8:30  | Arrival and Registration  |
| 9:00  | Welcome and Opening Remarks   |
| 9:10  | Antibody Kinetics on Challenging Targets   Gilbert van den Tillaart, Genmab B.V.  |
| 09:40 | Surface Plasmon Resonance (SPR) in Fragment-Based Drug Discovery (FBDD) – Targeting Oxidative Stress Proteins   Anders Bach, University of Copenhagen |
| 10:10 | Stabilizing Ni-NTA for Regeneration in a Quantitation Assay   Gemma Lindemann, Novonesis  |
| 10:20 | Networking Break  |
| 10:50 | iQue Based Screening for Improved Drug Selection for AML   Sandra Gordon, University of Copenhagen  |
| 11:20 | Octet as Facilitator for CMC Development   Lena Sørensen Hall, AGC Biologics AS   |
| 11:50 | Measurement of Antibody Avidity to Pertussis Toxin Using a BLI Method   Hanna Khan, University of Gothenburg  |
| 12:00 | CellCelector Flex: High-Throughput Isolation of Single Cells, Colonies and Organoids   Jurij Kintz, Sartorius   |
| 12:10 | Lunch Break and Networking  |
| 13:10 | Application of High-Throughput Flow Cytometry in Early Drug Discovery at AstraZeneca   Mei Ding, Astra Zeneca   |
| 13:40 | Expansion of the use of Octet in CMC and Research at Symphogen   Lotte Fynbo Hansen, Symphogen AS   |
| 14:10 | Clone Selection by Off-Rate Ranking   Pi Camilla Poulsen, Agilent Technologies Denmark APS  |
| 14:20 | Meet the Experts Sessions , Networking and Coffee Break   |
| 15:30 | End-to-End mAb Discovery: A Phage Display Campaign Supported by Sartorius Equipment   Cynthia de Bont, IPA (ImmunoPrecise Antibodies)                 |
| 16:00 | (Title TBA)   Caroline Bang Løppke, University of Copenhagen  |
| 16:20 | Cell Line Development Outsourcing: The Road from Virtual to Actual   Christopher Lupton, Sartorius  |
| 16:30 | Closing remarks   |