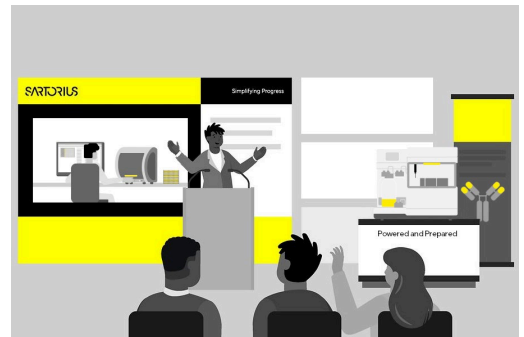


# Sartorius User Group Meeting

**Date:** Thursday, July 10<sup>th</sup>, 2025  
**Location:** University of Oxford  
Department of Biochemistry  
South Parks Rd, Oxford OX1 3QU

[Register NOW](#)



**8:30** Arrival and Registration

9:00 Welcome and Opening Remarks

9:10 Use of the iQue For Functional Vaccine Assessment | Dr Stephen Taylor, UKSHA

09:40 TBC | Dr Afreen Haider, Cancer Research Horizons

**10:10** Networking Break

10:30 A Biosensor-Based Phage Display Selection Method For Automated, High Throughput Nanobody Discovery | Phebe De Keyser, VUB Belgium

11:00 Investigation of Ligand-Induced Conformational Changes in Therapeutically Relevant Targets Using Octet BLI | Tomas Rodriguez, GSK

11:30 Probing Protein-Carbohydrate Interactions Using Glycosylated Gold Nanoparticles and BLI | Dr. Sarah-Jane Richards, University of Manchester

**12:00** Lunch Break and Networking

13:00 Advancing Life Sciences with Next-Gen Innovation | Dr. Daryl Cole, Sartorius

13:30 Application of the iQue3 for Antibody Discovery at LifeArc | Zaynab Isseljee, LifeArc

14:00 Quantifying Whole-Virus Avidity to Investigate the Receptor Binding Evolution of A(H5) Influenza Viruses | Dr. Lorin Adams, Francis Crick Institute

**14:30** Meet the Experts Sessions , Networking and Coffee Break

15:00 Application of BLI for Sequencing by Synthesis Fundamentals | Natasha Crake, Illumina

15:30 Automated Cell Killing Assay to Evaluate Cell Therapy Efficacy Via Imaging and Flow Cytometry | Dr Michela Serena, Arctoris

16:00 Design of Smart Polymers for Detection and Quantification of Viral Vectors on the BLI Octet | Dr. Emma, Daniels, Tozaro

16:30 Closing remarks