

Pushing the Boundaries in Biomolecular Interaction Analysis with the Creoptix WAVEsystem and waveRAPID

From biologics to Fragment-Based Drug Discovery

Program:

- 13:45 - 14:00 Grab your coffee & network
- 14:00 - 14:15 Welcome & Introduction
- 14:15 - 14:45 Pushing the Boundaries in Biomolecular Interaction Analysis with the Creoptix WAVEsystem and waveRAPID,
Andrea Moretti, Ph.D., Field Application Scientist, Creoptix
- 14:45 - 15:00 Q&A
- 15:00 - 15:30 Interactive instrument software overview
- 15:30 - 15:45 Coffee Break
- 15:45 - 16:15 Using GCI to overcome technical challenges in sensor chip assays,
Dr. Jana Wolf, Team leader at Domainex
- 16:15 - 16:45 Measuring affinities of peptides for the SPAK and OSR1 Conserved C-Terminal (CCT) Domains
Dr. Benjamin Bax, Reader in Structural Biology Medicines Discovery Institute, Cardiff University and Karen Elvers, Research Associate, Cardiff University
- 16:45 Closing & Apéro

United Kingdom
Cambridge

Vernalis HQ
20-09-22

Where: Vernalis HQ, Granta Park, Great Abington, Cambridge CB21 6GB, United Kingdom

When: 20/09/2022