

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,

Edward A. FitzGerald, Ph.D., Field Application Scientist, Creoptix

14:45 - 15:00 Q&A session

15:00 - 15:30 Interactive instrument software overview

15:30 - 15:45 Coffee break

15:45 - 16:15 Grating-coupled Interferometry and biophysical methods enabling structure based drug discovery on membrane protein targets,

Michael Hennig, Ph.D., CEO of leadXpro

16:15 - 16:30 Q&A session

16:30 - Closing & Apéro

**Switzerland
Schlieren**

**DiNAQOR HQ
07-09-22**

Where: DiNAQOR AG,
Wagistrasse 25, 8952 Schlieren

When: 07/09/2022