

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

From biologics to Fragment-Based Drug Discovery

Creoptix® invites you to join us for a **1 hour seminar** at the **University of Copenhagen** to learn about Molecular Interaction Analysis using **Grating-Coupled Interferometry (GCI)**.

GCI is a surface-based, **label-free** biosensing technique.

When target molecules (e.g. proteins) are attached to the sensor surface, binding of analytes can be detected by understanding the refractive index change close to the sensing surface.

With **high sensitivity**, the ability to resolve extremely **rapid dissociating kinetics** and innate compatibility with high molecular weight ratios, the Creoptix® WAVEsystem's GCI technology has a **broad range of applications**.

Denmark
Copenhagen

University of
Copenhagen
26-09-22

Join us if...

- You're interested in label-free technology
- You're working with kinetics of intermolecular interactions
- You're curious about the newest biotech developments

Where: University of Copenhagen, Novo Nordisk Foundation Center for Protein Research, Biophysics Unit, Blegdamsvej 3B, Copenhagen, Denmark

When: 26/09/2022, 10:00 am

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

