

ADME-Tox Strategies to Accelerate Drug Discovery

Wednesday, September 7, 2022

9 am – 4 pm

Symbion, Fruebjergvej 3,
2100 København Ø, Denmark

Agenda

- ▶ **09:00 – 10:20** Registration, Networking and Refreshments
 - ▶ **10:20 – 10:30** Welcome and Introduction
 - ▶ **10:30 – 11:15** Anders Sonesson PhD, Ferring Pharmaceuticals: *ADME Strategies for Peptide Drugs*
 - ▶ **11:15 – 12:00** Paul Walker PhD, Cyprotex Discovery Ltd: *The Use of Toxicogenomics (TGx) in Early Pre-clinical Safety Assessment*
 - ▶ **12:00 – 13:00** Buffet Lunch
 - ▶ **13:00 – 13:45** Janne Koch PhD, LEO Pharma A/S: *LEO Pharma Early Toxicology Strategy*
 - ▶ **13:45 – 14:30** Phil Butler PhD, Cyprotex Discovery Ltd: *From Small Molecules to New Modalities – The Changing Landscape of ADME Screening*
 - ▶ **14:30 – 16:00** Refreshments and Networking
-

Places are limited so please RSVP to reserve your place by contacting enquiries@cyprotex.com