

AGENDA

DAY 1			
TIMING	DURATION (min)	TITLE	SPEAKER
9:00-9:50	50	Registration and coffee	
9:50-10:00	10	Welcome and opening remarks	Péter Krajcsi, PhD, Chief Scientific Officer, SOLVO Biotechnology, Hungary
10:00-10:40	40	Keynote: Ki determination in vitro and its impact on modelling and prediction of DDI	Yuichi Sugiyama, PhD, Head of Sugiyama Laboratory, RIKEN, Japan
Session 1: Drug disposition and drug-drug interaction, Chair: Péter Krajcsi, PhD, Chief Scientific Officer, SOLVO Biotechnology			
10:40-11:10	30	Drug disposition classification systems - a comparative review of BDDCS, ECCS and ECCCS	Birk Poller, PhD, Lab Head, PK Sciences Department, Novartis, Switzerland
11:10-11:40	30	Assessment of unbound target site-concentration in brain and lungs: methodological considerations and practical implications	Irena Loryan, M.D., Ph.D., Researcher, Translational PKPD, Dept. Pharmaceutical Biosciences, Uppsala University
11:40-11:55	15	Poster presentation: Answering to the new FDA DDI guidance: Increasing the impact of IC50 studies through customization and modeling	Zhizhou Fang, PhD, Laboratory Head Drug Transporter, Senior Scientist II, DMP, Abbvie Ludwigshafen, Germany
11:55-13:00	65	Lunch Break	
13:00-13:40	40	Keynote presentation: Endogenous transporter substrates as clinical biomarkers	Dietmar Weitz, PhD, Principal Scientist, Drug Metabolism and Pharmacokinetics FF, Sanofi-Aventis Deutschland GmbH R&D, Germany
Session 2: Biomarkers, Chair: Krisztina Herédi-Szabó, PhD, Principal Scientist, ADME-Tox Services, SOLVO Biotechnology			
13:40-14:10	30	Mass spectrometry-based quantification of transporter proteins and metabolizing enzymes: an update on advantages and challenges	Stefan Oswald, PhD, Assistant Professor at the Department of Clinical Pharmacology, University Medicine Greifswald, Germany
14:10-14:40	30	A transporter, a nuclease and a metabolite: the trio orchestrating connective tissue calcification	András Váradi, PhD, Institute of Enzymology, RCNS, Hungarian Academy of Sciences, Budapest, Hungary
14:40-14:55	15	Poster presentation: Determination of reference values of MDR-ABC transporter activities in CD3+ T lymphocytes by using SOLVO MDQ Kit™	Katalin Filkor, PhD, Lead Scientist, MDQuest, Szeged, Hungary
14:55-15:25	30	Coffee Break	
15:25-16:05	40	Keynote presentation: Epacadostat: Transport and Metabolism of A First in Class IDO1 Inhibitor for Immuno-Oncology	Yan Zhang, Ph.D, Associate Director, Incyte Corporation, Drug Metabolism Pharmacokinetics & Clinical Pharmacology, US
Session 3: Disease and therapy, Chair: Roelof de Wilde, Director of Sales, SOLVO Biotechnology			
16:05-16:35	30	Transporters in antimalarial drug action and resistance	Sanjeev Krishna, MA, MBChB, FRCP, DPhil, ScD, FMedSci, Professor of Molecular Parasitology and Medicine, St. George's, University of London, UK
16:35-16:50	15	Poster presentation: Optimization of an in vitro assay for identification of NTCP inhibitors	Pieter Van Brantegem, Drug Delivery and Disposition, KU Leuven Department of Pharmaceutical and Sciences, Leuven, Belgium Pharmacological
16:50-17:05	15	Small Currents – Big Potential. Electrophysiology for transporters using the "SURFE2R" technology	Maria Barthmes, PhD, Application Scientist & Product Manager, SURFE2R Family, Nanion Technologies GmbH, Germany
17:05-19:00	115	POSTER SESSION	
19:00-22:00	180	DINNER	
DAY 2			
TIMING	DURATION (min)	TITLE	SPEAKER
8:20-9:00	40	Keynote presentation: Studying disposition of uremic waste products: can we develop novel therapeutic strategies to treat chronic kidney disease?	Roos Masereeuw, PhD, Professor and Head of the Experimental pharmacology group at Utrecht Institute for Pharmaceutical Sciences, The Netherlands
Session 4: Organ impairment/toxicity, Chair: Zsuzsanna Gaborik, PhD, Head of Laboratory, SOLVO Biotechnology			
9:00-9:30	30	Hepatotoxicity testing - primary hepatocyte and hepatic cell line models	Lysiane Richert, PhD, Professor in Cell Biology and Toxicology, Scientific Director, KaLy-Cell, France
9:30-10:00	30	Hepatic bile salt transport	Péter Krajcsi, PhD, Chief Scientific Officer, SOLVO Biotechnology, Hungary
10:00-10:30	30	Coffee Break	
10:30-11:10	40	Keynote presentation: Effect of cigarette smoke extract on the function and expression of membrane transporters in alveolar epithelial cells	Mikihisa Takano, PhD, Professor, Graduate School of Biomedical & Health Sciences, Hiroshima University, Japan
Session 5: Xenobiotics and endobiotics, Chair: Emese Kis, PhD, Director of R&D and Production, SOLVO Biotechnology			
11:10-11:40	30	Intestinal transport of fat-soluble vitamins	Emmanuelle Reboul, PhD, Researcher at the French National Institute of Agronomical Research (INRA), Research axis of the Micronutrition Team "Bioavailability of fat-soluble micronutrients", Center for Cardiovascular and Nutrition Research, Aix-Marseille University, France
11:40-12:10	30	Thyroid hormone transport	Theo Visser, PhD, Professor at Erasmus University Medical Center, Rotterdam, The Netherlands
12:10-12:25	15	Poster presentation: The role of human MRP2 in cellular efflux of microcystins, a type of cyanobacterial toxins	Gurjut Kaur, Human and Environmental Toxicology, University of Konstanz, Konstanz, Germany
12:25-12:35	15	Closing remarks	Péter Krajcsi, PhD, Chief Scientific Officer, SOLVO Biotechnology, Hungary
12:35-13:30	65	Lunch	