

Nanion User Meeting 2024 Agenda

Dates: October 22-23rd, 2024

Venue: Nanion Technologies HQ,

Tuesday, October 22nd

10:00 - 12:00 h CEST

User Workshop: SyncroPatch 384 (Registration required)

11:45 - 13:00 h CEST

Registration & welcome bites

13:00 - 13:15 h CEST

Niels Fertig (Nanion Technologies)

Welcome words

13:15 - 13:40 h CEST

Martin Gunthorpe (Autifony Therapeutics)

K_v3 channel positive allosteric modulators as a precision medicine approach for the treatment of progressive myoclonic epilepsy

13:40 - 14:05 h CEST

Joachim Wegener (University of Regensburg & Fraunhofer EMFT)

Profiling GPCR pharmacology using impedance monitoring of cell-based assays

14:05 - 14:30 h CEST

Christian Grimm (Ludwig-Maximilian University)*

A small GTPase is a direct effector of the intracellular Ca²⁺ channel TPC2 regulating melanoma progression through modulation of the Wnt signaling pathway

14:30 - 14:35 h CEST

Poster flash talks (x2, 2 min each)

14:35 - 16:00 h CEST

Instrument demos & Networking

16:00 - 16:25 h CEST

Alexandre Santinho (Oria Bioscience)

Inside out: Pioneering the future of organellar target screening

16:25 - 16:50 h CEST

Adriana Bizior (SB Drug Discovery)

Utilising SURFER technology to accelerate transporter drug discovery

16:50 - 17:15 h CEST

Surabhi Rajendra Kokane (Stockholm University)

PI-(3,5)P2 mediated oligomerization of the endosomal Na⁺/H⁺ exchanger NHE9

17:15 - 17:40 h CEST

Ulrich Hammes (Technical University Munich)

Substrate recognition and transport mechanism of the pin-formed auxin exporters

17:40 - 17:45 h CEST

Poster flash talks (x2, 2 min each)

17:45 - 18:45 h CEST

Posters & networking

19:00 - 23:00 h CEST

Dinner hosted by Nanion Technologies

Fräulein Wagner, Am Bavariapark 16, 80339 Munich

*This talk will be only available to the attendees on-site. It will not be broadcasted to the virtual attendees

Wednesday, October 23rd

08:45 - 09:15 h CEST

Coffee & tea

09:15 - 09:40 h CEST

Stefan Kubick (B4 PharmaTech)

Cell-free systems for the production and functional characterization of membrane proteins

09:40 - 10:05 h CEST

Gregor Anderluh (National Institute of Chemistry)

Pore forming toxins: from molecular mechanisms to nanosensing

10:05 - 10:30 h CEST

Tobias Ensslen (Hahn-Schickard Institute)

Real-Time Peptide Differentiation: Insights from Nanopore Technology

10:30 - 11:00 h CEST

Coffee break

11:00 - 11:25 h CEST

Jamie Bhagwan (Axol Bioscience)

Chamber-specific pharmacological responses of axoCells™ hiPSC-derived Atrial and Ventricular Cardiomyocytes on the FLEXcyte96 platform

11:25 - 11:50 h CEST

Jieun An (NEXEL)

Utilizing Cardiosight®-S with the CardioExcyte96 for Cardiotoxicity Screening

11:50 - 12:15 h CEST

Jamie Vandenberg (Victor Chang Cardiac Research Institute)

High throughput phenotyping of cardiac ion channel variants

12:15 - 13:30 h CEST

Lunch

13:30 - 13:55 h CEST

Janina Sörmann (University of Copenhagen)*

Leveraging Machine Learning tools to design hASIC1a modulators

13:55 - 14:20 h CEST

Rajnish Ranjan (EPFL)

CHANNELOME: Towards a paradigm shift in drug screening

14:20 - 14:45 h CEST

Karen Elvers & Iwan Williams (Cardiff University)

Enabling Drug Discovery with the SyncroPatch 384

14:45 - 15:10 h CEST

Alexandr Ilyaskin (Friedrich-Alexander-Universität)

Structure-based analysis of proteolytic and ligand-mediated ENaC activation

15:10 - 15:20 h CEST

Ali Obergrussberger (Nanion Technologies)

Closing words