

Nanion User Meeting 2025 Agenda

Dates: October 28-29th, 2025

Venue: Ganghoferstraße 66b, 80339 München

Tuesday, October 28th

09:45 - 12:00 h CEST

User instrument Workshops (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

Christian Grimm (Ludwig-Maximilians University)

High-throughput organellar electrophysiology to allow early Drug discovery from native Lysosomes

13:40 - 14:05 h CEST

Ann-Katrin Piper (University of Wollongong)

Potential Role of the CLCA1-ANO1 Axis in Supporting Survival of Gastric Cancer Circulating Tumour Cells

14:05 - 14:15 h CEST

Poster flash talks (x3, 3 min each)

14:15 - 14:30 h CEST

Short break

14:30 - 15:15 h CEST

Instrument demos - Rotation 1

15:15 - 16:00 h CEST

Instrument demos - Rotation 2

16:00 - 16:15 h CEST

Short break

16:15 - 16:40 h CEST

Jan Behrends (University of Freiburg)

Simultaneous high-resolution fluorescence and voltage clamp measurements on free-standing membranes on a chip

16:40 - 17:05 h CEST

Anasua Mukhopahyay (University of Fribourg)

Single-Molecule Detection of Protein Biomarkers in Neurodegenerative Disease

17:05 - 17:30 h CEST

Marianna Misioni (EPFL University)

Fluorescence-Based Approaches to Study Biological Nanopore Mechanisms

17:30 - 17:40 h CEST

Poster flash talks (x3, 3 min each)

17:40 - 18:45 h CEST

Posters & networking

19:00 - 23:00 h CEST

Dinner hosted by Nanion Technologies

Wednesday, October 29th

08:45 - 09:15 h CEST

Coffee & tea

09:15 - 09:40 h CEST

David Colameo (University of Zurich)

Exploring Heterogeneous Cell Populations Using Automated Patch-Clamp and Correlative Imaging

09:40 - 10:05 h CEST

Michel de Waard (University of Nantes)

Photosensitive natural peptides for the optical control of ion channels

10:05 - 10:30 h CEST

Stephan Pless (University of Copenhagen)

Leveraging deep learning tools to design miniprotein modulators of ion channels

10:30 - 11:00 h CEST

Coffee break

11:00 - 11:25 h CEST

Francesco Tadini-Buoninsegni (University of Florence)

Investigating charge transport by P-type ATPases with the SURFER technology

11:25 - 11:50 h CEST

Ines Benhammouche (Aarhus University)

Structural and mechanistic insights into PIN8 and its inhibition by Morphactins

11:50 - 12:15 h CEST

Felix Baerenz (Sanofi)

High-throughput SLC inhibition assays: Combining SSM electrophysiology & ADE-MS

12:15 - 13:30 h CEST

Lunch

13:30 - 13:55 h CEST

Terence Hébert (McGill University)

GPCR-based drug discovery using iPSC-derived cardiomyocytes and cardiac fibroblasts

13:55 - 14:20 h CEST

Bettina Lickiss (innoVITRO)

In Vitro Modeling of Atrial Fibrillation Using Human iPSC-Derived Cardiomyocytes for Therapeutic Screening

14:20 - 14:35 h CEST

Julia Erl (Regensburg University)

Monitoring the Reversibility of GPCR Signaling by Combining Photochromic Ligands with Label-free Impedance Analysis

14:35 - 14:50 h CEST

Tobias Ensslen (Hahn-Schickard)

Real-Time Monitoring of Pore-Forming Toxin Activity in Live Cells Using High-Throughput Impedance Assays

14:50 - 15:00 h CEST

Elena Dragicevic (Nanion Technologies)

Closing words