

nanji[on

**Novel Insights Into Ion
Channel and Transporter
Research:**

From Silent Translocation to Deep
Learning Tools

Monday, February 17
11:30 AM – 1:00 PM
Room 404AB

Novel Insights Into Ion Channel and Transporter Research: From Silent Translocation to Deep Learning Tools



11:30 AM – 11:40 AM
Dr. Tim Strassmaier
Director of Scientific Operations
Nanon Technologies, USA

Novel Insights into Ion Channel & Transporter Research: From Silent Translocation to Deep Learning Tools

For over 2 decades, Nanion Technologies has been providing diverse solutions for electrophysiologists worldwide. We aim to implement innovative technologies in the fields of ion channel automated patch clamp (APC) electrophysiology, monitoring of cell viability and contractility, as well as electrogenic transporters, with various throughput capabilities. Our symposium will start with an introduction by Dr. Strassmaier who will describe the latest advances in Nanion's assays and product portfolio, followed by our speakers, whose work focuses on ion channel and transporter research.



11:40 AM – 12:05 PM
Dr. Janina Sörmann
Postdoctoral Reseracher
University of Copenhagen

Leveraging Deep Learning Tools to Design hASIC1a Modulators

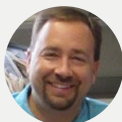
Stroke is a leading cause of death and adult disability. Acid-sensing ion channels (ASICs) in the central nervous system contribute to tissue damage after ischemic stroke but are underexplored therapeutic targets due to the lack of potent, selective, and stable inhibitors. During this presentation, Dr. Sörmann will describe de novo designed miniproteins capable of inhibiting hASIC1a by integrating deep learning tools in the drug discovery process and combined with functional analysis on the SyncroPatch 384. These computationally designed modulators offer high specificity, efficacy, and cost-effective synthesis, broadening accessibility for stakeholders.



12:05 PM – 12:30 PM
Dr. Rocco Zerlotti
Application Scientist
Nanon Technologies, Germany

Benefits of SSME for the Investigation of the Lysosomal Ion Channels TMEM175, TPC2, and TRPML1

Solid-supported membrane-based electrophysiology (SSME) is a technique for studying transport and electrogenic activities in transporters, pumps, and ion channels, including those located in inner membranes such as mitochondria and lysosomes. Using a fluidic system for rapid solution exchange, transport or ion conductance are driven by substrate concentration gradients at zero membrane potential. Dr. Zerlotti will present a recent study focused on investigation of lysosomal channels TMEM175, TPC2 and TRPML1. Measurements of K^+ and H^+ permeability in TMEM175, revealed two K^+ permeability coefficients and a PH/PK ratio consistent with literature. For TPC2, saturable Na^+ conductance with an EC_{50} of ~ 40 mM was observed, enhanced by the potentiator TPC2-A1P. SSME offers valuable insights into lysosomal channel function, complementing traditional patch-clamp techniques.

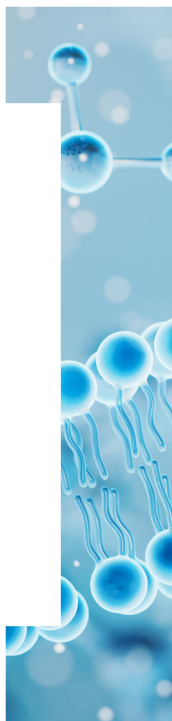


12:30 PM – 12:55 PM
Dr. Dan Burden
Professor, Analytical Chemistry
Wheaton College

Integrated Optical and Electrical Techniques for Investigating Unsupported Lipid Bilayer Dynamics at the Single-Molecule Level

This presentation examines the microelectrode cavity array (MECA) platform for studying biomolecule dynamics in lipid bilayers, emphasizing simultaneous optical and electrical measurements. The platform detects single

accelerate your research



**Find your nearest Nanion
representant on our website:
www.nanion.de**

nan|i[on