



You Never Walk Alone

Join the Nanion Week at Biophysics

Come See Us Booth 509

Friday, February 14th

2025 Satellite Meeting Drug Discovery for Ion Channels XXV

8:00 AM - 5:15 PM, Room 501ABC

Los Angeles Convention Center

Saturday, February 15th

Channels, Receptors & Transporters Subgroup Symposia

8:30 AM - 12:30 PM, Petree Hall C

Los Angeles Convention Center

Nanion is a proud sponsor.

Nanoscale Approaches to Biology Subgroup Symposia

8:30 AM - 12:30 PM, Room 502A;

Los Angeles Convention Center

Nanion is a proud sponsor.

Sunday, February 16th

Poster #B337 1:45 PM - 3:45 PM

Light-Induced Action Potentials & Stem Cell-Derived Sensory Neurons: an Automated Patch Clamp Investigation.

Poster #B343 1:45 PM - 3:45 PM

Block of Voltage- and Ligand-Gated Ion Channels by Toxins Using Automated Patch Clamp - a Detection Tool?

Nanion User Reception 6:30 pm

Registration: **Sign up at booth 509** to save your spot at our User Reception. Registration is required!

Monday, February 17th

Nanion Workshop: 11:30 AM - 1:00 PM, Room 404AB
Novel Insights Into Ion Channel & Transporter Research:
From Silent Translocation to Deep Learning Tools

Poster #B402 1:45 PM - 3:45 PM
Functional Characterization of Lysosomal Channels TMEM175 &
TPC2 Using Solid Supported Membrane-Based electrophysiology.

Poster #B329 1:45 PM - 3:45 PM
Investigating the Effect of Temperature & Trafficking Modulators on
hERG Current Using High Throughput Automated Patch Clamp.

Tuesday, February 18th

Poster #B103 1:45 PM - 3:45 PM
Quantitative & Mechanistic Analysis of Protein Membrane Pore
Formation on a High-Throughput Electrophysiology Platform.

Come See Us: Membrane Physiology Symposium

Napa California, April 30 - May 1 2025

Scan Here To Reserve You Spot
Space is Limited, So Act Fast!

