



SOPHION USER MEETING

August 21 - 22, 2024 | Hilton La Jolla Torrey Pines, San Diego, CA

Sophion North America User Meeting 2024

August 21

03:30 PM - 04:30 PM	Keynote Address - Daniel Minor Ph.D. - UCSF <i>Resisting a toxic environment - frog saxiphilins in action</i>
04:30 PM - 06:30 PM	Welcome Reception: Drinks and Appetizers

August 22

08:30 AM - 09:30 AM	Check-In/Welcome Remarks <i>News and Views from Sophion</i>
09:30 AM - 10:00 AM	Michel Jurgens and Daniel Sauter Ph.D. – Sophion Bioscience <i>Virtual Screening for modulators of GABA, , KCa, and Nav1.7 channels: Does it work?</i>
10:00 AM - 10:30 AM	Heike Wulff Ph.D. – University of California at Davis
10:30 AM – 10:45 AM	Break <i>Cryo-EM Structures of the Small-Conductance Calcium-Activated Potassium Channel</i>
10:45 AM - 11:15 AM	Miao Zhang Ph.D. – Chapman University



SOPHION USER MEETING

August 21 - 22, 2024 | Hilton La Jolla Torrey Pines, San Diego, CA

Sophion North America User Meeting 2024

August 22 cont'd

11:45 AM - 12:15 PM	<p><i>Speed is the Game: A Deadly Combination of Genomic Medicine and Automated Patch Clamp</i></p> <p>Bikash Pattnaik Ph.D. – University of Wisconsin at Madison</p>
12:15 PM – 1:15 PM	Lunch
1:15 PM – 1:45PM	<p><i>Automated Patch Camp: Tips and Tricks</i></p> <p>Pietro Scaduto Ph.D. – Sophion Bioscience</p>
1:45 PM - 2:15 PM	<p><i>Kv 7 .x channels using Qube</i></p> <p>Jennifer Wesley – Eurofins Discovery Services</p>
2:15 PM - 2:45 PM	<p><i>"Exploring the Structural Mechanics of PIEZO Ion Channels with Fluorescence Nanoscopy"</i></p> <p>Eric Mulhall Ph.D. – Scripps Research Institute</p>
2:45 PM – 3:00 PM	Break
3:00 PM - 3:30 PM	<p><i>Measuring If in atrial-like cardiomyocytes utilizing automated patch clamping</i></p> <p>Soomin Song Ph.D. – NYU IonLab</p>
3:30 PM - 4:00 PM	<p><i>Streamlining Disease Modeling and Drug Discovery using iPSC derived Induced Excitatory Neurons</i></p> <p>Scott Schachtele Ph. D. – Fujifilm CDI</p>
4:00 PM - 4:30 PM	Closing Remarks