

Sophion North America User Meeting

Harvard Medical School - August 24 - 25th 2023

Day 1 – August 24th

10:00 AM – 10:30 AM	Coffee and Registration
10:30 AM – 10:35 AM	Samantha Cramer - Sophion Bioscience Welcome Remarks
10:35 AM – 11:05 AM	Michael Bohnen - Columbia University hiPSC Derived Cardiomyocyte Recordings on Qube384
11:05 AM – 11:35 AM	Michael Sonnested - Sophion Bioscience Sophion Bioscience Update
11:35 AM – 12:05 PM	Jonas Friard - Seguela Lab - McGill University Comparison of Human Sensing Ion Channel properties using QPatch 16x
12:05 PM – 1:05 PM	Lunch
1:05 PM – 1:35 PM	Brooke Ahern - US Army DEVCOM CBC Discrepancies of Nav1.4 Inhibition by Toxins is due to Structural Variability
1:35 PM – 2:05 PM	Kyle Jensen - Sophion Bioscience Cannabinoid Modulation of Voltage gated Calcium Channels
2:05 PM – 2:35 PM	Reza Ghovanloo - Waxman Lab - Yale University Cannabinol functionally-selectively inhibits threshold sodium currents and reduces excitability in neurons of the dorsal root ganglion
2:35 PM – 3:00 PM	Coffee Break
3:00 PM – 3:30 PM	Shanti Amagasu - Latigo Biotherapeutics Implementation of a robust Kv1.x QPatch II assay
3:30 PM – 4:15 PM	QPatch Compact Features Demonstration A review of the newly updated QPatch Compact
4:15 PM – 5:00 PM	Drinks Reception

Sophion North America User Meeting

Harvard Medical School - August 24 - 25th 2023

Day 2 – August 25th

10:00 AM – 10:30 AM	Welcome and Registration
10:30 AM – 11:00 AM	Patrick Walsh - Anatomic Incorporation iPSC Derived Cell Allays on Automated Patch Clamp
11:00 AM – 11:30 AM	Andrew Cook - Eurofins Discovery Services Robust High Throughput CiPA Screening at Physiological Temperature
11:30 AM – 12:00 PM	Sooyeon Jo & Patric Vaelli - Bruce Bean Lab - Harvard University Characterization of the Nav1.8 inhibitors VX-150 and VX-548 by Qube and QPC
12:00 PM – 12:30 PM	Brent Dawe - Janssen Pharmaceuticals