



# Nanion Automated Patch Clamp User Meeting 2017

**Friday, June 9<sup>th</sup> 2017**

8:15 AM to 5:30 PM

**Hyatt Regency Hotel, 575 Memorial Dr.  
Cambridge MA 02139**

---

**Friday, June 9<sup>th</sup> 2017**

8:15AM to 9:00AM

**Breakfast and Registration –Charles View Ballroom 16<sup>th</sup> Floor**

9:00AM to 9:15AM

**Welcome & Introduction: Rodolfo Haedo**, Senior Vice President  
Nanion Technologies

9:15AM to 10:00AM

**-Keynote- Dr. Chris Miller, Professor. Biochemistry, Brandeis  
University** "A Weird Ion Channel for a Very Weird Ion"

10:00AM to 10:30AM

**Dr. Alfred George, Professor and Chair,  
Department of Pharmacology Northwestern University**  
"Automated electrophysiology of iPSC-derived cardiomyocytes"

10:30AM to 11:00AM

**Dr. Tim Strassmaier Senior Scientist Nanion** "Pursuit of TRP channel  
targets with the SyncroPatch384PE"

11:00AM to 11:15AM

**Break (Coffee and snacks) Posters Foyer Area 16<sup>th</sup> Floor**

11:15AM to 11:45AM

**Andrew Allen, and Dr. David Baez** Broad Institute "The Agony  
and the Ecstasy: Characterization of Disease-Related Ion  
Channel Variants using the SyncroPatch 384®"

---

11:45AM to 12:15PM **Dr. Bjorn Knollmann, Director, Vanderbilt Center for Arrhythmia Research** “Modeling heart disease with human iPSC-derived cardiomyocytes – First Experiences with CardioExcyte96”

12:15PM to 1:15PM **Lunch (16<sup>th</sup> Floor)**

**-Afternoon Session-**

---

**Friday, June 9<sup>th</sup> 2017**

1:15PM to 2:00PM **-Keynote Presentation- Dr. Owen McManus, CTO Q-State Biosciences Inc.** “Phenotypic Assays of Neuronal function to support Drug Discovery”

2:00PM to 2:30PM **Dr. Maria Barthmes Application Scientist Nanion Munich** “Targeting Transporters: High throughput screening and characterization of membrane transporters and pumps with the SURFE2R 96SE”

2:30PM to 3:00PM **Dr. Matt Fuller Senior Scientist, Icagen** “Advancing Early Drug Discovery with the SyncroPatch 384PE at Icagen”

3:00PM to 3:30PM **Dr. Greg Kaczorowski, CEO Kanalis** “Lessons Learned in Drugging Ion Channels from Prosecuting the Novel ROMK Target

3:30PM to 4:00PM **-Poster and Coffee Break- Foyer Area 16<sup>th</sup> Floor**

4:00PM to 4:30PM **Dr. Srinivasan Venkatachalan, Assist Principal Scientist, Chromocell** “Ion channel Therapeutic Discovery and Development effort at Chromocell Corporation using Patchliner”

4:30PM to 5:00PM **Dr. Carlos Vanoye Research Assoc. Professor Northwestern University** “ High Throughput Functional Analysis of Ion Channel Gene Variants ”

---

