



Neuroscience 2025

San Diego, CA. November 15-19

Booth #835

Thank you for visiting our booth and we hope you enjoy your time at Neuroscience 2025. We encourage you to visit the posters below and speak to current Maestro MEA users about how our tools helped their research. For information about Axion news and giveaways during this conference, or a closer look at our many neural applications, please visit: axionbiosystems.com/axion-biosystems-sfn-2025.



Contents

Saturday, November 15 th	1
Sunday, November 16 th	2
Monday, November 17 th	2
Tuesday, November 18 th	3
Wednesday, November 19 th	4

Are we missing your poster? Let us know!

Please send your poster with [Axion products](#) to contact@axionbio.com

Saturday, November 15th

1 PM – 5 PM

- » **Astrocytes: Disease Mechanisms I:** Contribution of astrocytic Sparcl1 to cortical synaptic dysfunction in C9orf72-FTD/ALS. R. A. Culibrk *et al.* *Barrow Neurolog. Inst.* [PSTR009.09 / C17](#)
- » **Developmental Disorders:** GxE in a dish: common genetic variants altering neurodevelopment in response to valproic acid, an environmental risk factor for autism. A. B. Marquez Gonzalez *et al.* *UNC at Chapel Hill.* [LBP006.25 / LBP055](#)

[Return to top >>](#)



- » **Neurogenesis and Gliogenesis:** Robust production of parvalbumin interneurons and fast-spiking neurons from human medial ganglionic eminence organoids. *M. P. Walker et al. Univ. of Michigan.* [LBP001.08 / LBP008](#)
- » **Nicotinic Acetylcholine Receptors:** Sar guided discovery of positive allosteric modulators of $\alpha 9/\alpha 10$ nicotinic acetylcholine receptors for the treatment of hidden hearing loss. *P. Sapkota et al. Idaho State Univ.* [PSTR005.07 / B15](#)
- » **Oscillatory, Aperiodic, and Waveform Analysis of Electrophysiological Recordings:** Developmental trajectory of rhythmic and non-rhythmic spiking in hiPSC-derived cortical tissue. *A. Hutton et al. Univ. of California San Diego.* [PSTR007.13 / B48](#)

Sunday, November 16th

8 AM – 12 PM

- » **Motorneuron-Muscle Interface and Muscle Physiology/Biochemistry:** Coupling Compartmentalized Microfluidic Platforms with MEA for Advancing Neuromuscular Junction Modeling. *N. Bagdadi. NETRI, France.* [PSTR078.07 / Q3](#)
- » **Motorneuron-Muscle Interface and Muscle Physiology/Biochemistry:** Novel microfluidic chambers for longitudinal profiling of brain cortical organoids using graphene-based optical stimulation and microelectrode arrays. *T. Zhou et al. University of California San Diego.* [LBP020.13 / LBP049](#)

1 PM – 5 PM

- » **Preclinical Pain Models and Measurements:** Digital Signature Library: using neurons as universal bio-digital sensors. *A. Toma. NETRI, France.* [PSTR117.09 / G11](#)
- » **Preclinical Pain Models and Measurements:** Neuron-as-a-sensor : a detailed method to map CIPN. *B. Maisonneuve. NETRI, France.* [PSTR117.08 / G10](#)

Monday, November 17th

8 AM – 12 PM

- » **Electrophysiology: Electrode Arrays I:** Optimization of electrical stimulation in microelectrode arrays to reduce variability in functional assays of neuronal network activity. *S. A. Chvatal et al. Axion BioSystems.* [PSTR196.12 / WW1](#)
- » **Mechanisms and Propagation of Alpha-Synuclein:** Neuroprotective mechanisms of endosulfine-alpha in cellular models of Parkinson's Disease. *G. Drafor et al. Purdue Univ.* [PSTR158.02 / D29](#)
- » **Synaptic Plasticity and Structural Remodeling: Mechanisms Shaping Brain Function and Dysfunction:** Selective Blockade of Microtubule Invasion into Dendritic Spines Illuminates Cytoskeletal Roles in Synaptic Plasticity. *H. L. Miller et al. Univ. of Wisconsin.* [PSTR152.11 / B29](#)

[Return to top >>](#)



1 PM – 5 PM

- » **Molecular Mechanisms Underlying Developmental Disorders II:** Methylmercury's impact on redox homeostasis and parvalbumin-expressing interneuron maturation. *M. Watanabe. Univ. of Hawaii.* [PSTR201.11 / A51](#)

Tuesday, November 18th

8 AM – 12 PM

- » **Spinal Pain Mechanisms and Circuitry:** Molecular and Functional Characterization of hiPSC-Derived Dorsal Horn Neurons for Central Pain Target Discovery. *V. Truong et al. Anatomic Inc.* [PSTR277.15 / L2](#)
- » **Using Pluripotent Stem Cells to Model Human Diseases:** Embodiment of the insoluble: Investigating nanoplastic neurotoxicity in models of the developing human brain. *A. Hudock et al. Univ. of California San Diego.* [PSTR255.30 / A30](#)

1 PM – 5 PM

- » **Biotechniques: Gene, Protein, or Cell Based Approaches:** Co-culture of frozen neural progenitors, neuron precursors, and astrocytes derived from the highly characterized reference scti003-a line. *A. Moosa et al. STEMCELL Technologies Inc.* [PSTR367.06 / YY11](#)
- » **Parkinson's Disease: Cellular Mechanisms:** Evaluating TMEM175 mutation phenotypes in human iPSC-derived dopaminergic neuron models for Parkinson's disease. *S. Schachtele et al. Fufifilm Cell. Dynamics.* [PSTR325.23 / D54](#)
- » **Spinal Cord Injury: Plasticity and Recovery II:** The Effects of Optogenetic Stimulation on Human Induced Pluripotent Stem Cell-Derived V2a Spinal Interneurons. *M. Dominguez et al. Gladstone Inst.* [PSTR335.14 / N1](#)
- » **Treatment and Drug Discovery in Mood Disorders:** Discovery of aminopyrazole-based multitarget antidepressants: Virtual screening, molecular docking, and preliminary SAR and toxicity studies. *I. Y. Attah et al. Univ. of Cape Coast.* [PSTR352.14 / JJ16](#)

[Return to top >>](#)



Wednesday, November 19th

8 AM – 12 PM

- » **Addictive Drugs: Neural Pathways, Networks, and Transcription:** Role of KCNQ2/3 channels in striatal response to dopamine and cocaine conditioned place preference. *E. T. Jorgensen et al. Univ. of Alabama at Birmingham.* [PSTR408.09 / JJ17](#)
- » **Disorders of Neuromuscular Junction and Muscle:** A Human Pluripotent Stem Cell-Derived Neuromuscular Junction Model Reveals Botulinum Neurotoxin A-Induced Synaptic Dysfunction. *C. A. Aguayo et al. Wake Forest Inst. For Regenerative Med.* [PSTR385.03 / H1](#)
- » **Pluripotent Stem Cell Neuronal Differentiation, Specification, Organoids, and Assembloids:** High throughput assessment of barrier function using human ipsc-derived brain microvascular endothelial cells. *R. Vaidyanathan et al. Fujifilm Cell. Dynamics.* [PSTR369.11 / A11](#)
- » **Pluripotent Stem Cell Neuronal Differentiation, Specification, Organoids, and Assembloids:** Functional analysis for the screening of iPSC-derived neural organoid response to neuroactive compounds. *B. Streeter et al. Axion BioSystems.* [PSTR369.16 / A16](#)
- » **Pluripotent Stem Cell Neuronal Differentiation, Specification, Organoids, and Assembloids:** Characterization of Human iPSC-Derived Forebrain Neurons for Organoid Modeling of Memory. *A. L. Woehr et al. Wake Forest Inst. Regen. Med.* [PSTR369.26 / A26](#)

1 PM – 5 PM

- » **Brain Injury: Therapeutic Strategies:** Sustained BDNF delivery mitigates secondary injury and enhances recovery in a rodent model of traumatic brain injury. *N. Hong et al. Dankook Univ.* [PSTR444.09 / L4](#)
- » **Molecular Mechanisms Underlying Developmental Disorders III:** Neuronal hyperactivity becomes mTORC1 independent due to epigenetic changes in Tuberous Sclerosis Complex. *W. Afshar Saber et al. Boston Children's Hosp.* [PSTR425.03 / A22](#)

[Return to top >>](#)