



International Society for Stem Cell Research 2018

Melbourne, Australia

June 20th-23rd

Focus Sessions

Wednesday, June 20

Room 212/213, Level 2 | 9:00 am – 12:00 pm

Focus Session: BlueRock: USING PLURIPOTENT STEM CELLS TO TREAT DISEASES OF THE NERVOUS SYSTEM

9:50 - 10:20 Developing patient-specific stem cell models to inform new epilepsy therapies.

Evangelos Kiskinis, Northwestern University, IL

Poster Sessions

Wednesday, June 20

Exhibit Hall | 6:30 - 8:30 pm

W-2012 Human Induced Pluripotent Stem Cell Derived Neurons with Chromosome 1Q21.1 Deletions Display Altered Cortical Patterning and Synaptic Regulation.

Gareth Chapman: PhD Student, Cardiff University, UK

W-2024 Accelerated Maturation and Improved Functionality of Human iPSC-Derived Neurons with the B-27TM Plus Neuronal Culture System.

Alex Hannay Thermo Fisher Scientific, MD

W-2028 High-Content Imaging of iPSC-Derived Human Neurons for Toxicity Screening.

Michael Hendrickson: BrainXell, Inc., WI



W-2130 **Multiwell microelectrode array technology for the evaluation of human iPSC-derived cardiomyocytes and neuron development and maturation.**

Mike Clements: Axion BioSystems, GA

Thursday, June 21

Exhibit Hall | 6:00 - 8:00 pm

T-1029 **Characterization of human iPSC-derived cardiomyocyte electrophysiology with the local extracellular action potential assay.**

Mike Clements: Axion BioSystems, GA

T-2027 **Rapid Generation of Mature Cortical and Spinal Astrocytes from Human iPSCs.**

Michael Hendrickson: BrainXell, Inc.

T-3019 **BrainPhys Neuronal Medium Supports the Maturation and Electrical Activity of Human Pluripotent Stem Cell (hPSC)-Derived Neurons in Long-Term Cultures.**

Vivian Lee: STEMCELL Technologies

T-3056 **Spontaneous Network Activity Development in Cultured Cortical Neurons: Comparison Between Rat Embryonic and human Pluripotent Stem Cell-Derived Systems.**

Susanna Narkilahti: University of Tampere

Friday, June 22

Exhibit Hall | 6:00 - 8:00 pm

F-1038 **Non-Integrating RNA-Based Modifications of hPSC-Derived Cardiomyocytes – Applications for Optogenetics, siRNA Knockdown and Dominant Negative Expression of Mutated Transcripts.**

Felix von Haniel: Ncardia

F-2027 **Functional Phenotypic Screening of Patient iPSC-Derived Motor Neurons – *In Vitro* HTS Disease Modeling with Micro Electrode Arrays Coupled with AI-Based Analysis Methods.**

Michael Hendrickson: BrainXell, Inc.

F-2030 Establishment of Microfluidics Platform for Compartmentalized Neuronal Culture with Axonal Propagation Velocity Measurement.

Mervi Ristola: University of Tampere

F-4006 Investigating population variability using high throughput electrophysiological phenotyping of human induced pluripotent stem cell-derived cardiomyocytes.

Melissa M. Mangala: Victor Chang cardiac research institute