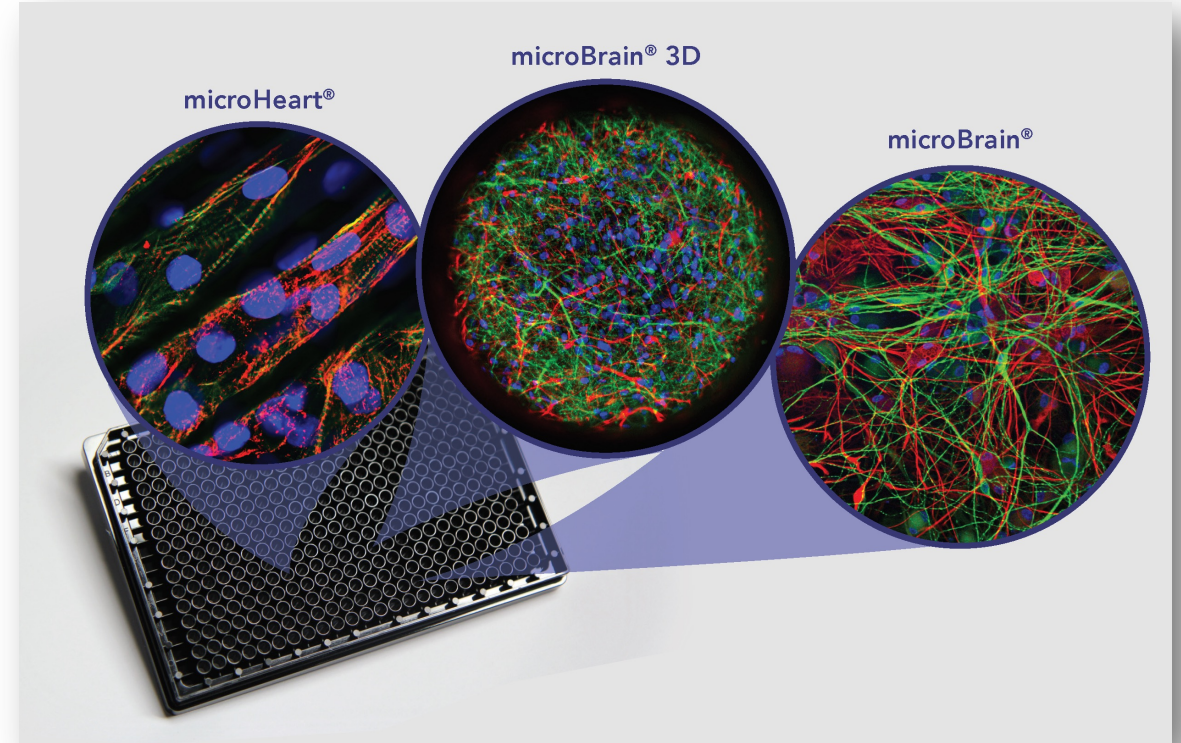


Functional and mechanistic neurotoxicity profiling using human iPSC-derived neural 3D cultures

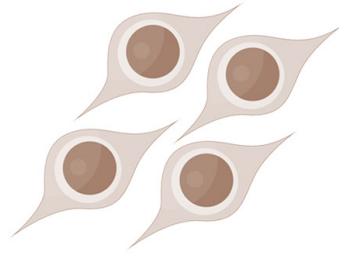
Cassiano Carromeu, PhD - StemoniX

StemoniX[®]

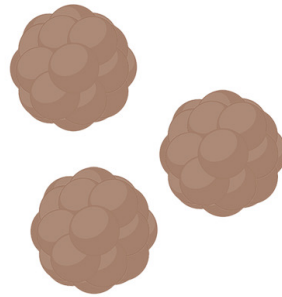
- Expertise in human iPSC-derived neural and cardiac platforms
- Focus on adapting platforms to HTS and producing assay-ready solutions to accelerate drug discovery
- Neural platform: microBrain[®]



Models available for Neuroscience research



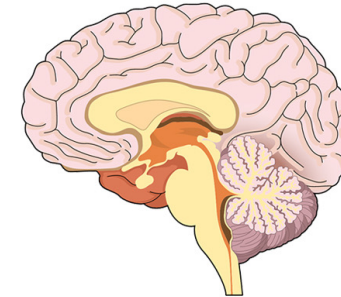
Cellular models



Organoids



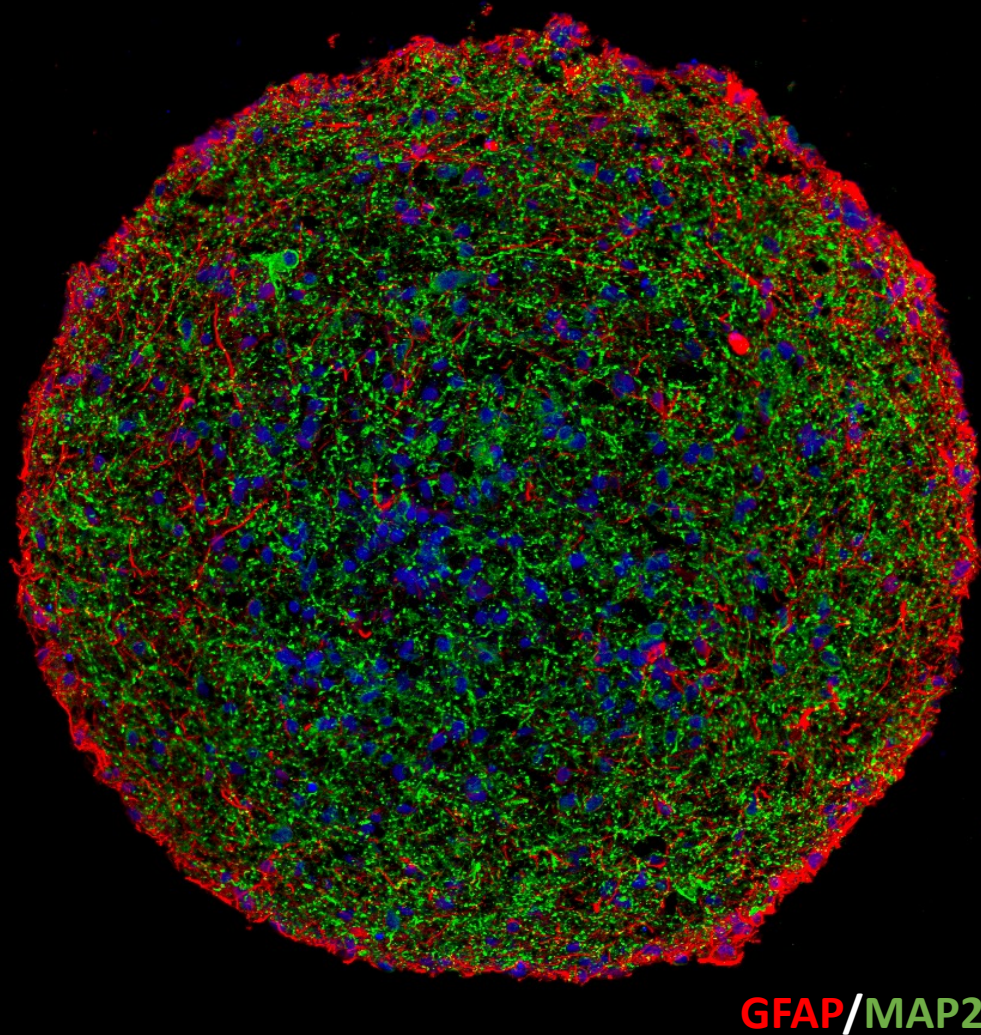
Animal models



Brain



microBrain[®] 3D

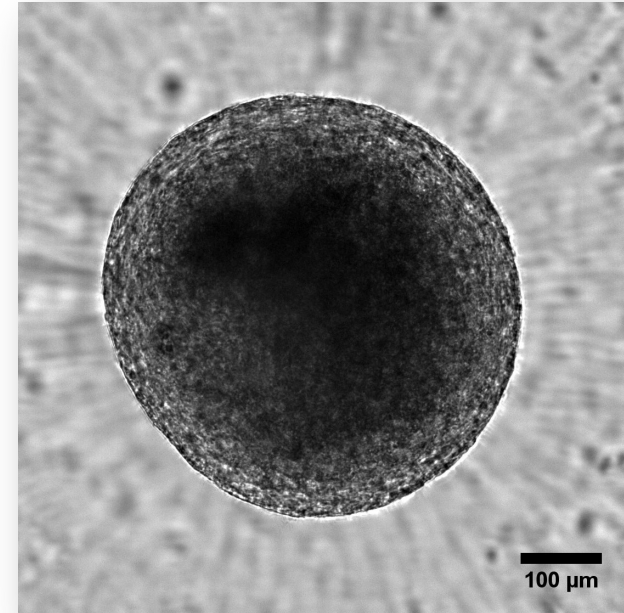
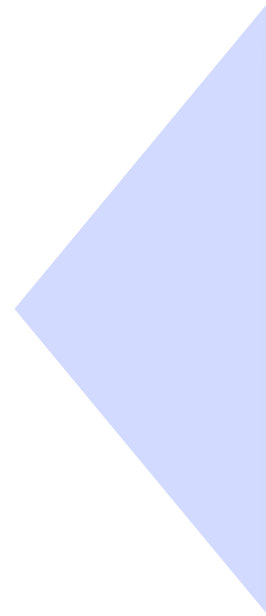
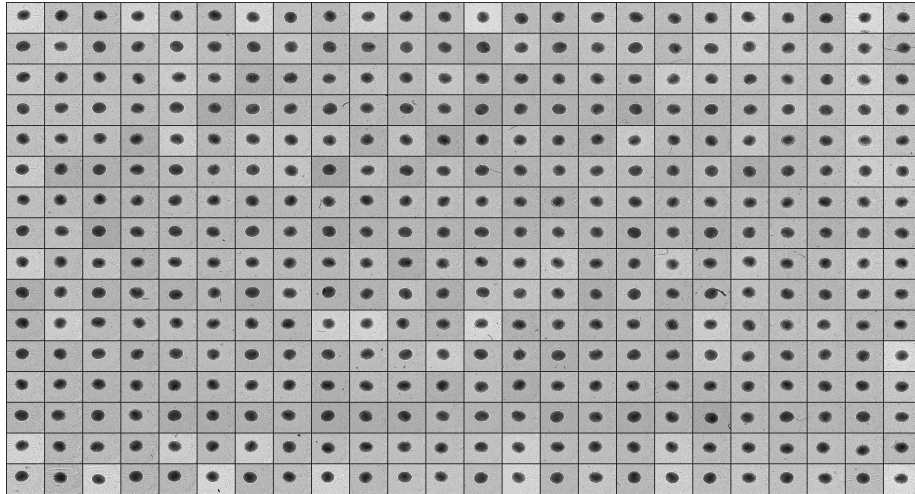


Main Features

- Single donor human iPSC line
- Balanced co-culture of neurons and astrocytes
- Display key neuronal and astrocytes markers
- Spontaneous synchronized activity
- Amenable to High Content Screening and High Content Imaging (384-well format)

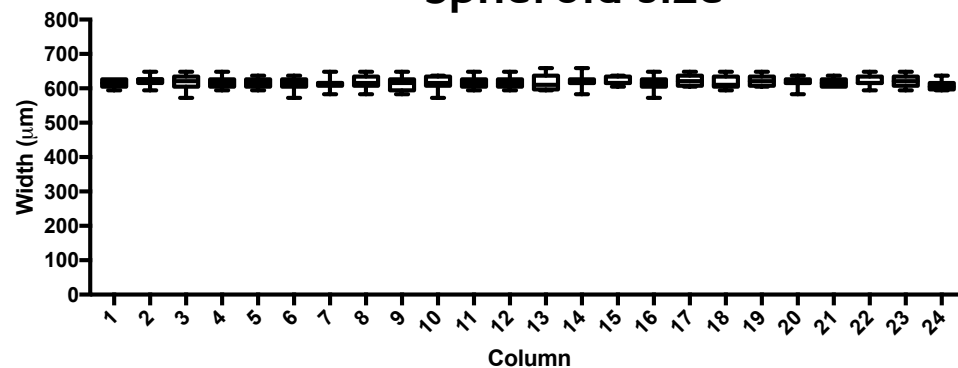
microBrain 3D spheroids are homogenous

384w microBrain 3D plate



microBrain 3D spheroid

Spheroid size

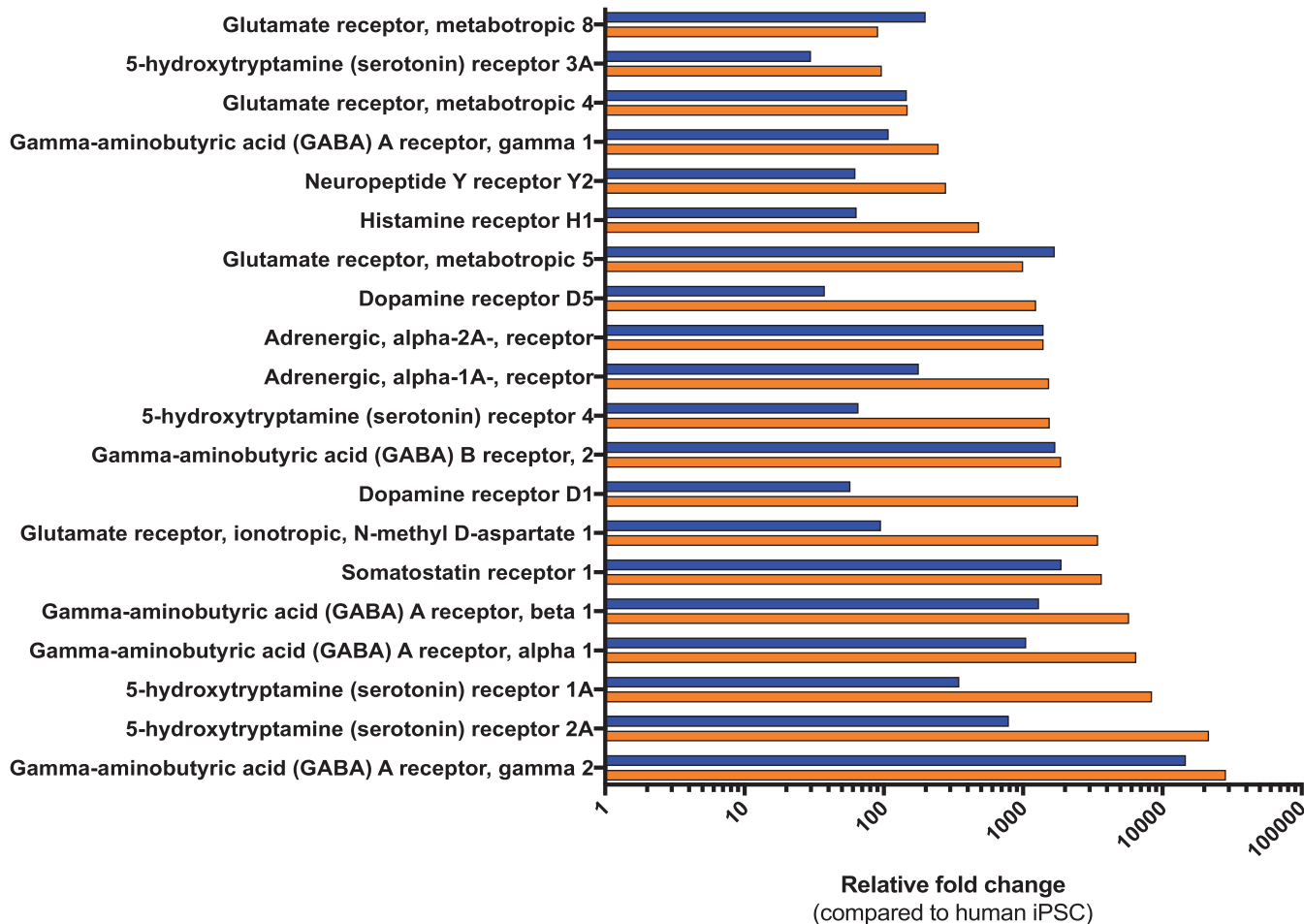


Features:

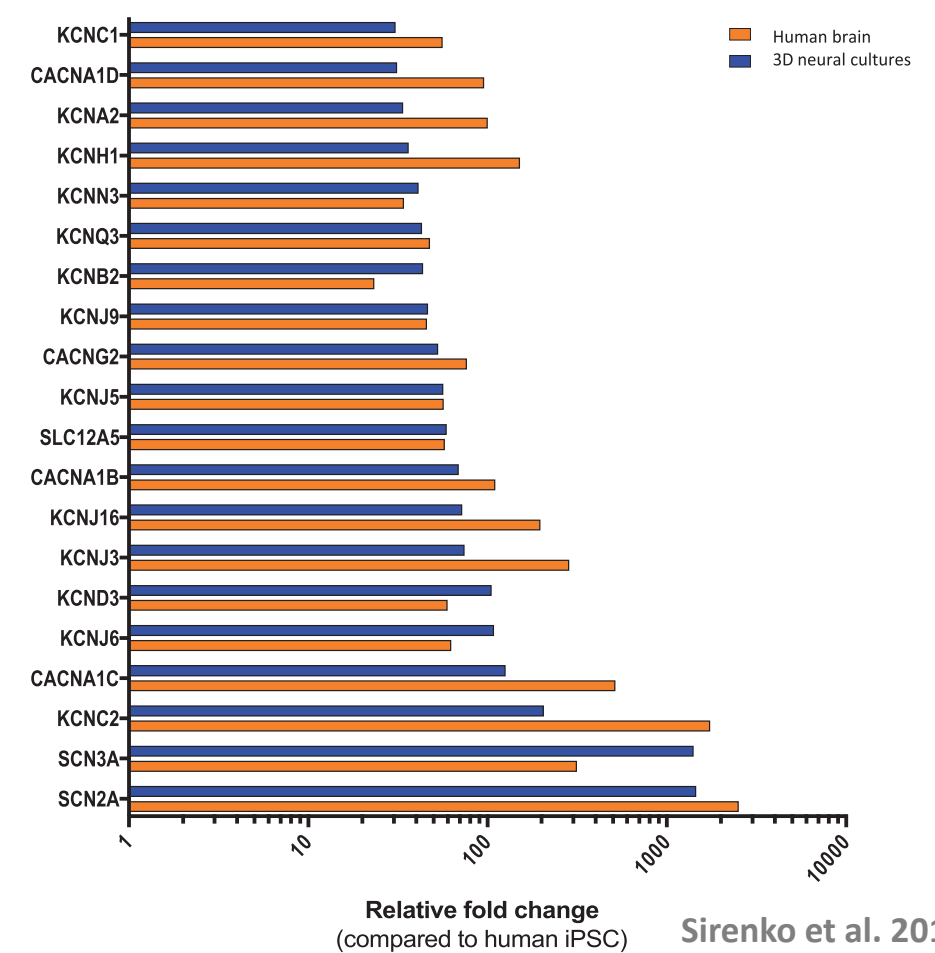
- 1 spheroid per well
- Consistent size across the plate
- Able to culture for weeks

Gene expression profile of microBrain 3D

Neurotransmitter profiling



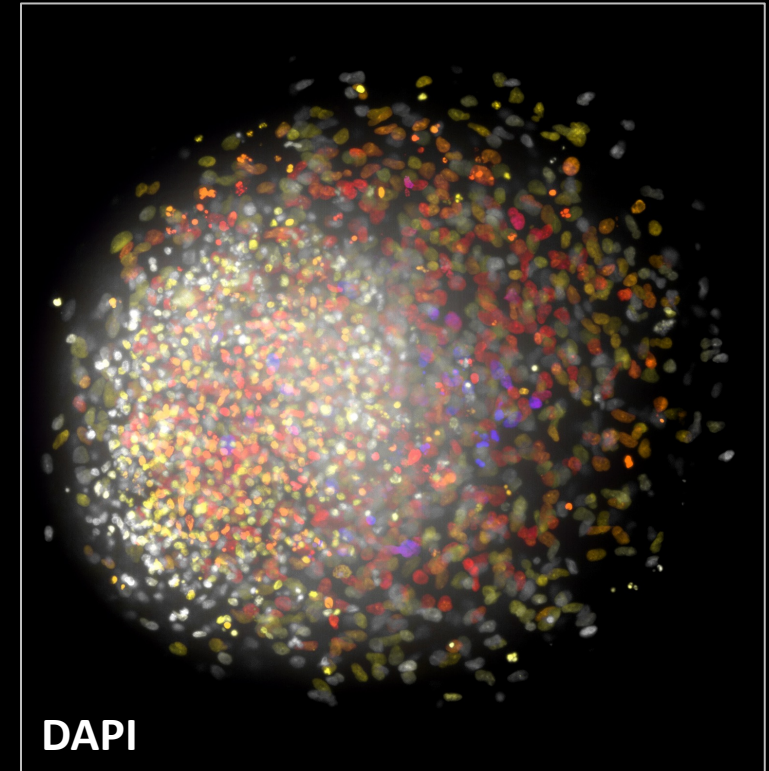
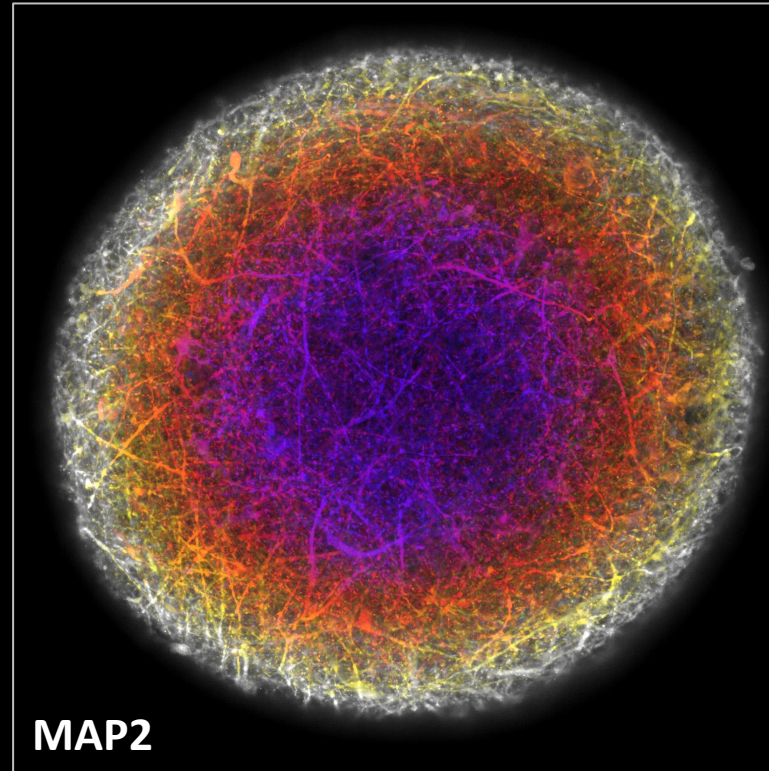
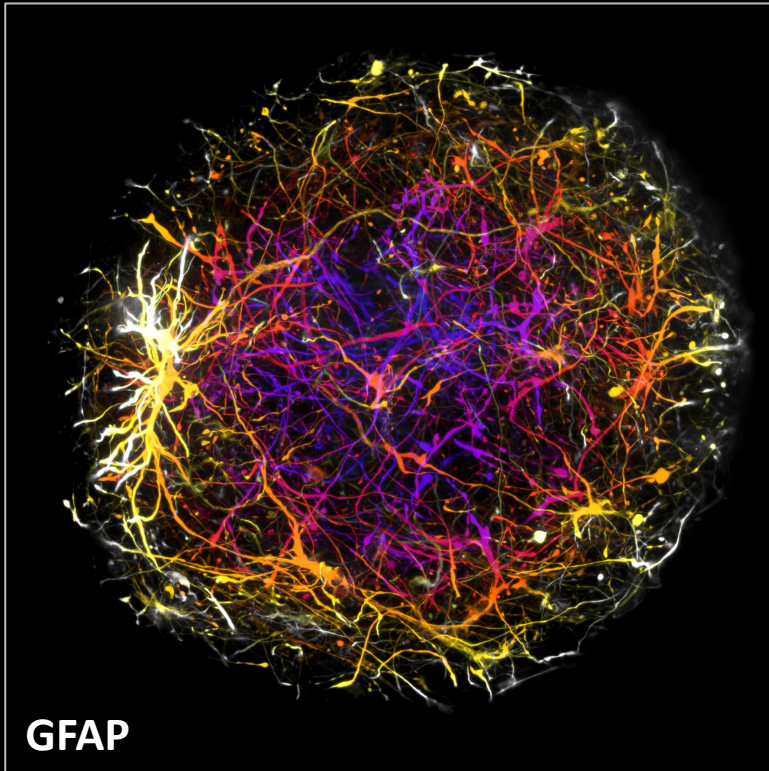
Ion channel profiling



Sirenko et al. 2018

microBrain 3D neural composition

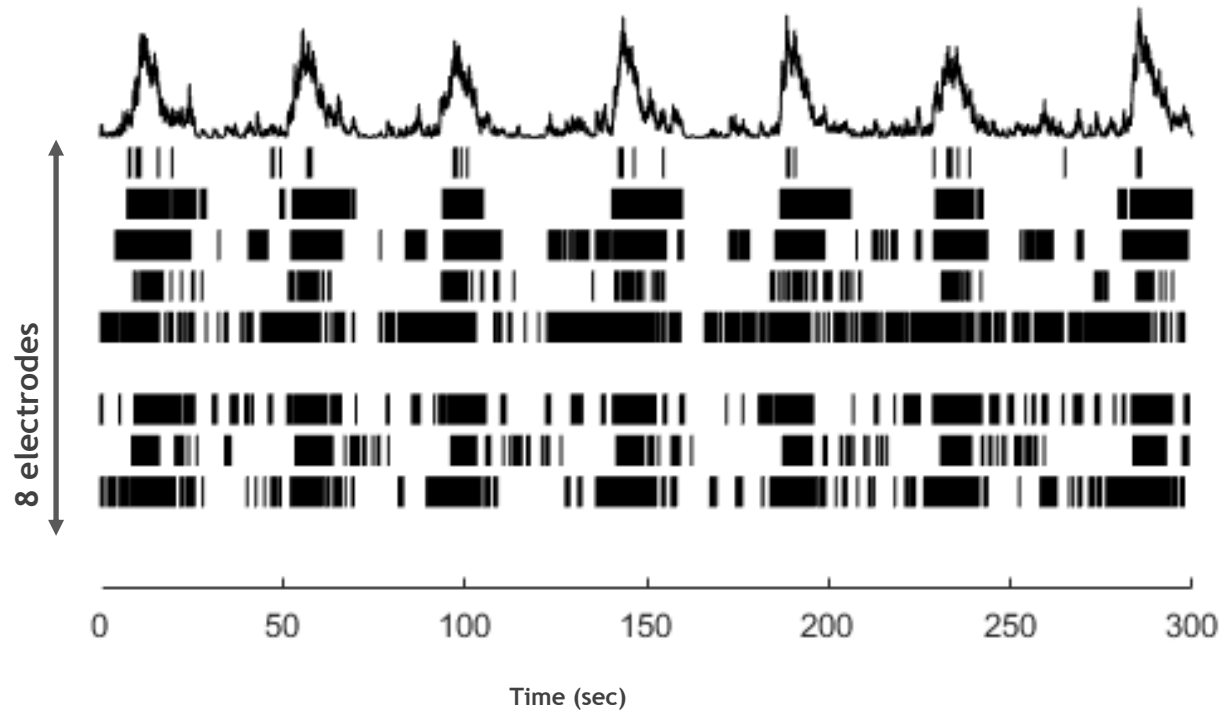
- Immunostaining with neuronal and glial markers on microBrain spheroids
- Confocal images using ImageXpress[®] Micro Confocal system (IXM-C)
- Optical clearing protocol applied after staining



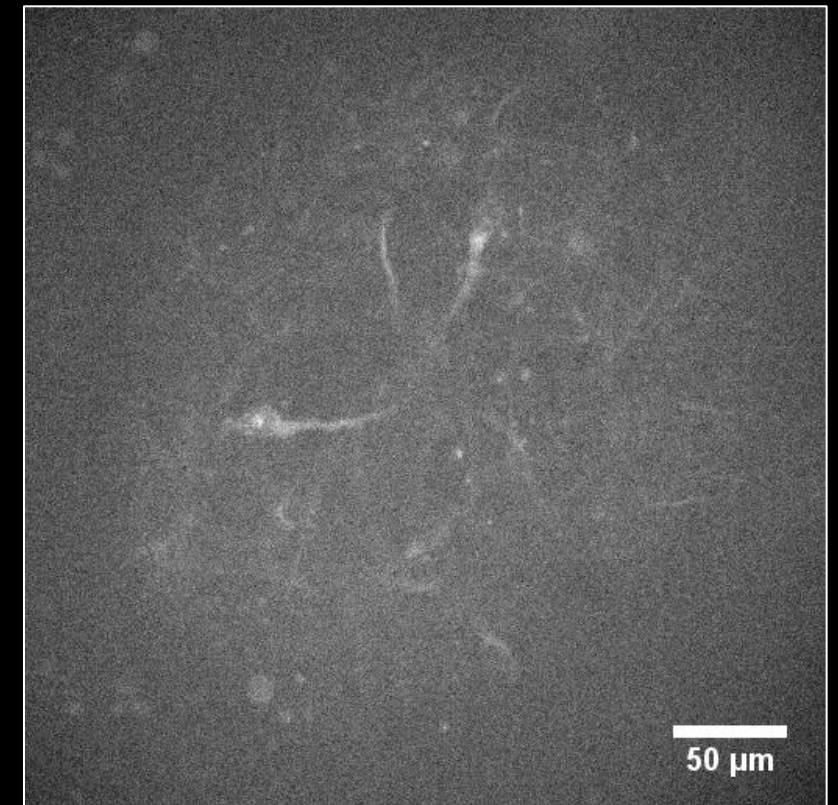
microBrain 3D demonstrates spontaneous synchronized neural activity

- Spheroid electrical activity is confirmed via MEA recordings
- Neuronal origin of activity is confirmed via synapsin-targeted Ca^{2+} measurements with high-speed confocal microscopy (IXM-C)

Plated microBrain 3D on MEA (Maestro)



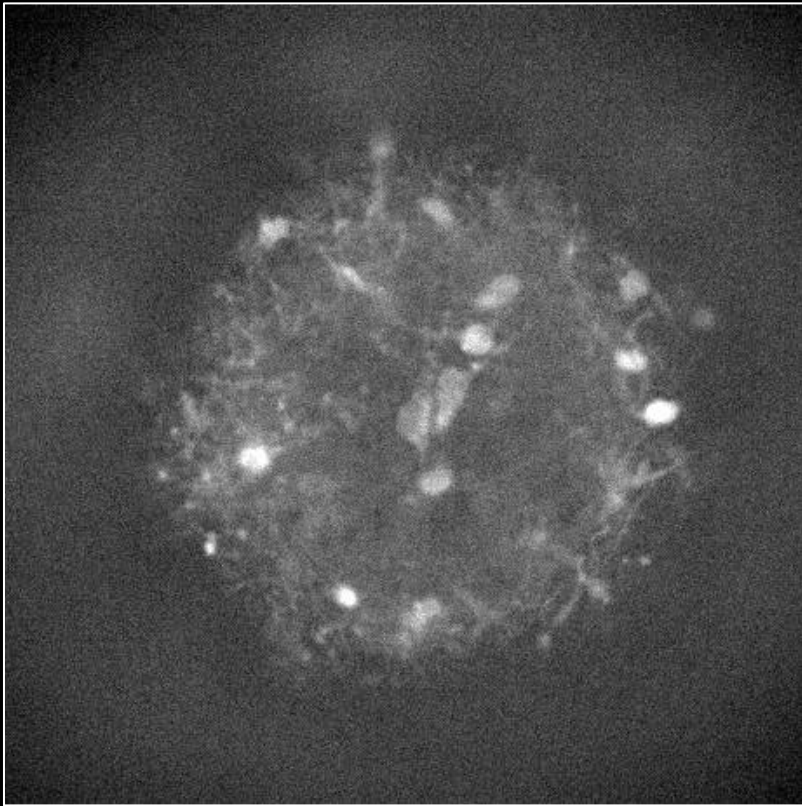
Synapsin driven Ca^{2+} indicator



microBrain 3D activity can be monitored in high throughput format

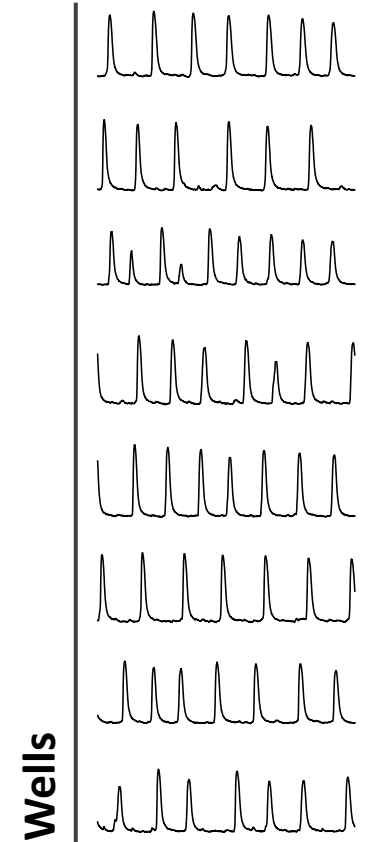
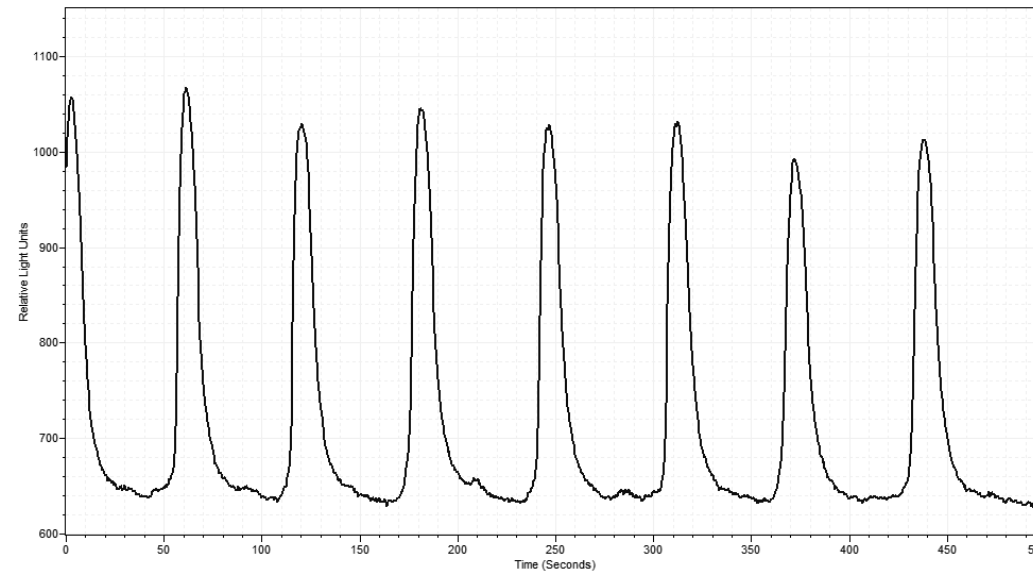
- Neuronal activity is monitored as calcium oscillations detected by high throughput kinetic fluorescence (FLIPR® Tetra System)

Spontaneous neuronal activity



FLIPR Tetra System

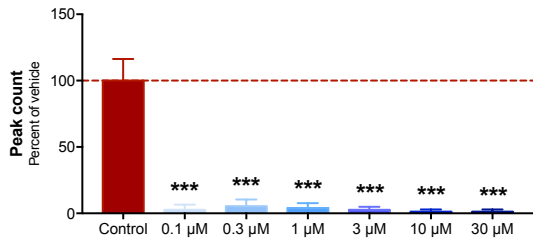
- Spontaneous activity on one well
- Detected oscillations correspond to synchronized calcium oscillation occurring on the sphere



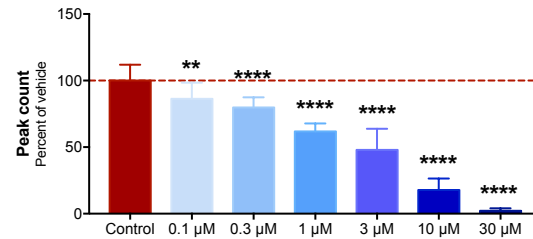
microBrain 3D activity respond to known neuromodulators

- Spontaneous neural activity can be monitored by high throughput calcium flux analysis (FLIPR Tetra System)
- Graphs show modulation on spontaneous oscillations 30min after adding compounds

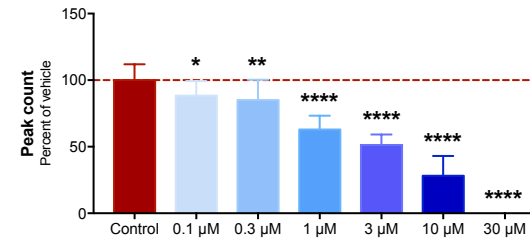
Tetrodotoxin



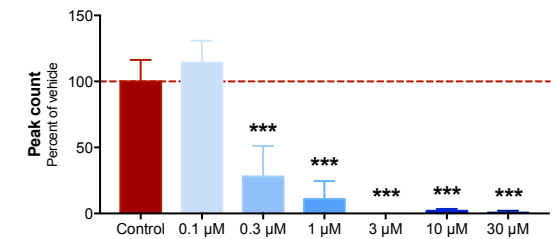
CNQX



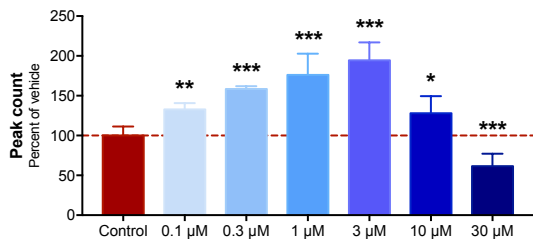
MK-801



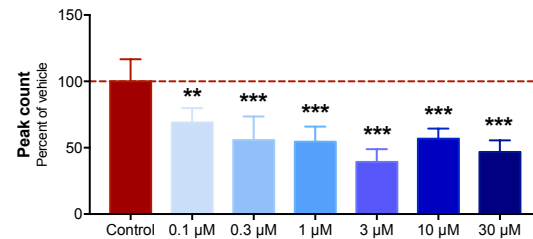
Muscimol



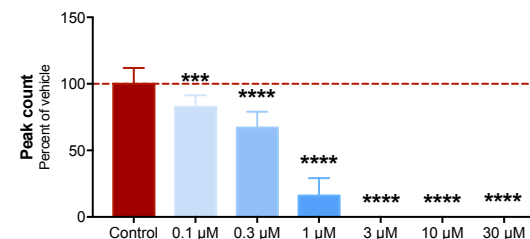
Kainic Acid



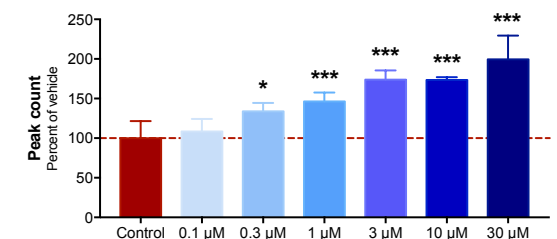
NBQX



MK-801 + CNQX



4-Aminopyridine



Tetrodotoxin: sodium channel blocker (potent neurotoxin)

Kainic Acid: Kainate receptor agonist

MK-801: NMDA receptor antagonist

CNQX: competitive AMPA/kainate receptor antagonist

NBQX: AMPA receptor antagonist

Muscimol: potent GABA agonist

4-Aminopyridine: Non-selective K_v channel blocker

microBrain 3D showcase

Drug Discovery / Screening:

- Epilepsy Drug Discovery
- High Throughput Screening of LOPAC®1280
- Thousands of compounds screened as a service with high reproducibility among replicates

Safety Pharmacology and Investigative Toxicology:

- Neurotoxicity Screen of Drug Candidates to Treat Zika Infection
- Screening of Environmental Toxins for Neurotoxic Effects

Disease Modeling:

- Neurodevelopmental disorders (Rett Syndrome)

Screening of Environmental Toxins for Neurotoxic Effects

- Neurotoxicity is a major reason for attrition in the final stages of a drug development pipeline.
- Current neurotoxic evaluation of chemicals rely mostly on animal models.
- Among hiPSC neurotoxicity assays, neurite outgrowth and multi-electrode arrays (MEAs) have been used as tool for evaluation of toxicity on the Central Nervous System (CNS).
- There is a need to more predictive and scalable human *in vitro* models to test compounds for their toxicological effects on the CNS.

Screening of Environmental Toxins for Neurotoxic Effects

- Goal:

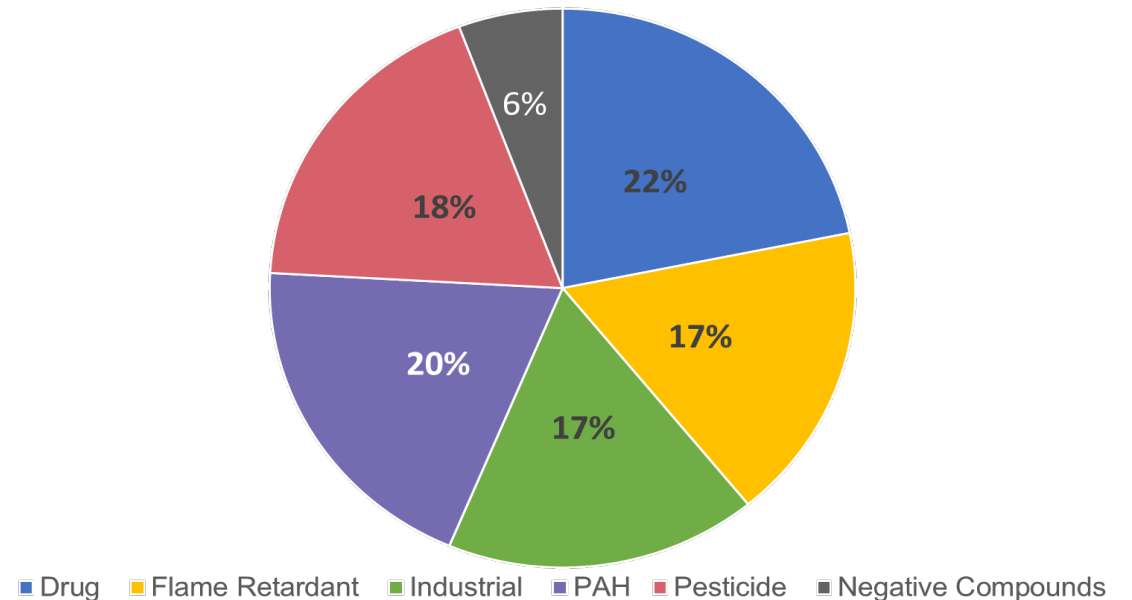
We used microBrain 3D platform to characterize the phenotypic responses to various compounds by monitoring the impact on the frequency and pattern of the spontaneous calcium oscillations.

Screening of Environmental Toxins for Neurotoxic Effects

- Method:

We screened a diverse library of 91 compounds comprised of representative examples of compounds from various environmentally relevant and potentially neurotoxic groups using an HTS (FLIPR Tetra System) and HCS (ImageXpress Micro Confocal system) workflow.

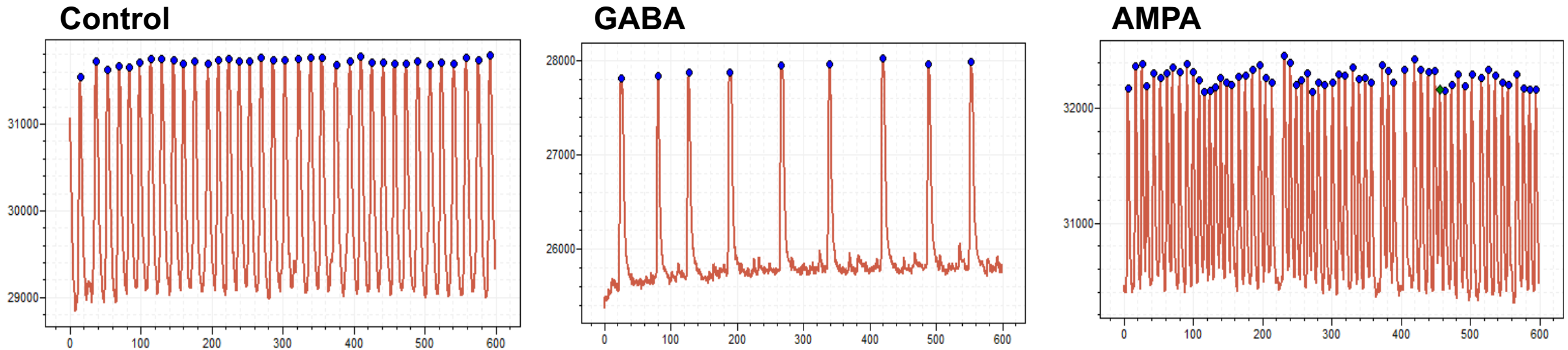
Compound Library Categories



microBrain 3D activity modulation

- Representative FLIPR Tetra System recordings of spontaneous neuronal activity

Calcium oscillation traces recorded by the FLIPR Tetra System:

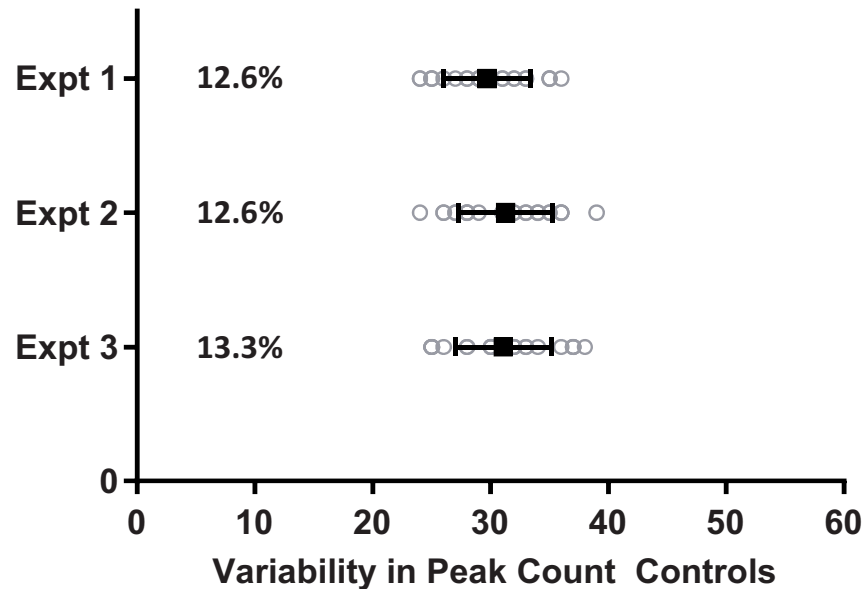


Sirenko et al. 2018

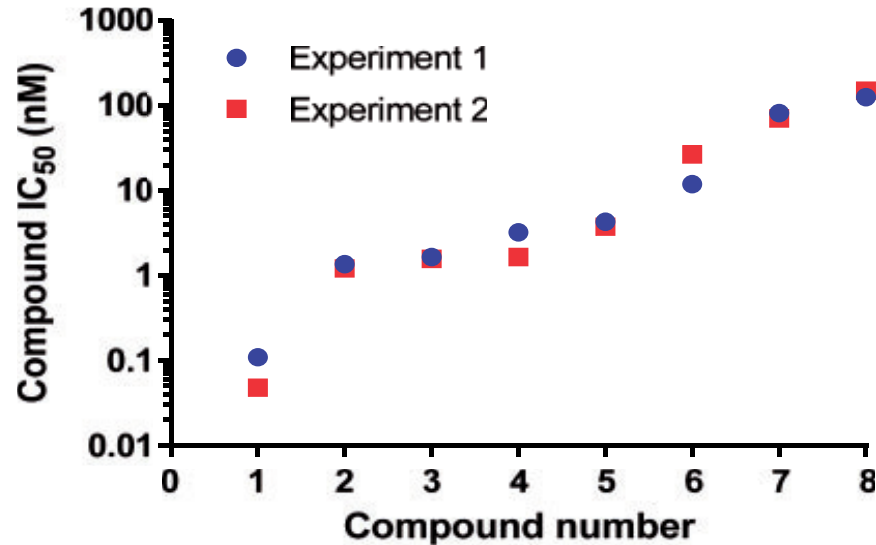
microBrain 3D activity is reproducible

- Inter-plate variability of spontaneous neuronal activity
- Each experiment number correspond to a different microBrain 3D plate

Inter-plate variability in vehicle controls



Inter-plate compound response comparison



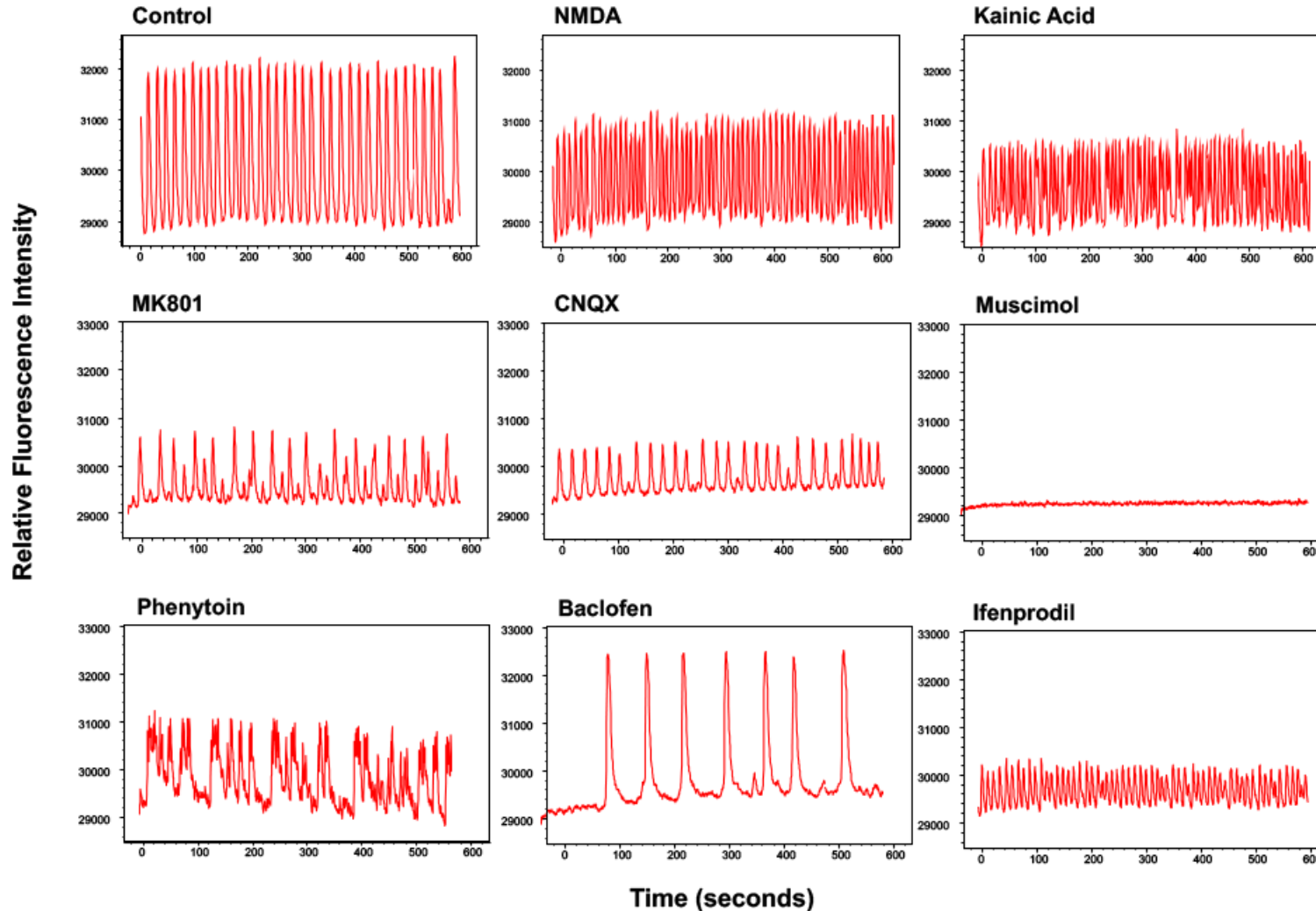
Compound Legend:

- ¹Rotenone
- ²2-Ethylhexyl diphenyl phosphate
- ³Methyl mercuric (II) chloride
- ⁴Isodecyl diphenyl phosphate
- ⁵2,2',4,4',5-Pentabromodiphenyl ether
- ⁶2,2',4,4'-Tetrabromodiphenyl ether
- ⁷Lead (II) acetate trihydrate
- ⁸1-Ethyl-3-methylimidazolium diethylphosphate

Sirenko et al. 2018

microBrain 3D activity modulation

- Representative FLIPR Tetra System recordings of spontaneous neuronal activity



Legend:

- Control: Vehicle DMSO control
- NMDA: NMDA receptor agonist
- Kainic Acid: Kainate receptor agonist
- MK-801: NMDA receptor antagonist
- CNQX: AMPA/Kainate receptor antagonist
- Muscimol: GABA_A agonist
- Phenytoin: Blocker of voltage gated Na⁺ channels
- Baclofen: GABA_B agonist
- Ifenprodil: non-competitive NMDA antagonist

Sirenko et al. 2018

microBrain 3D activity modulation

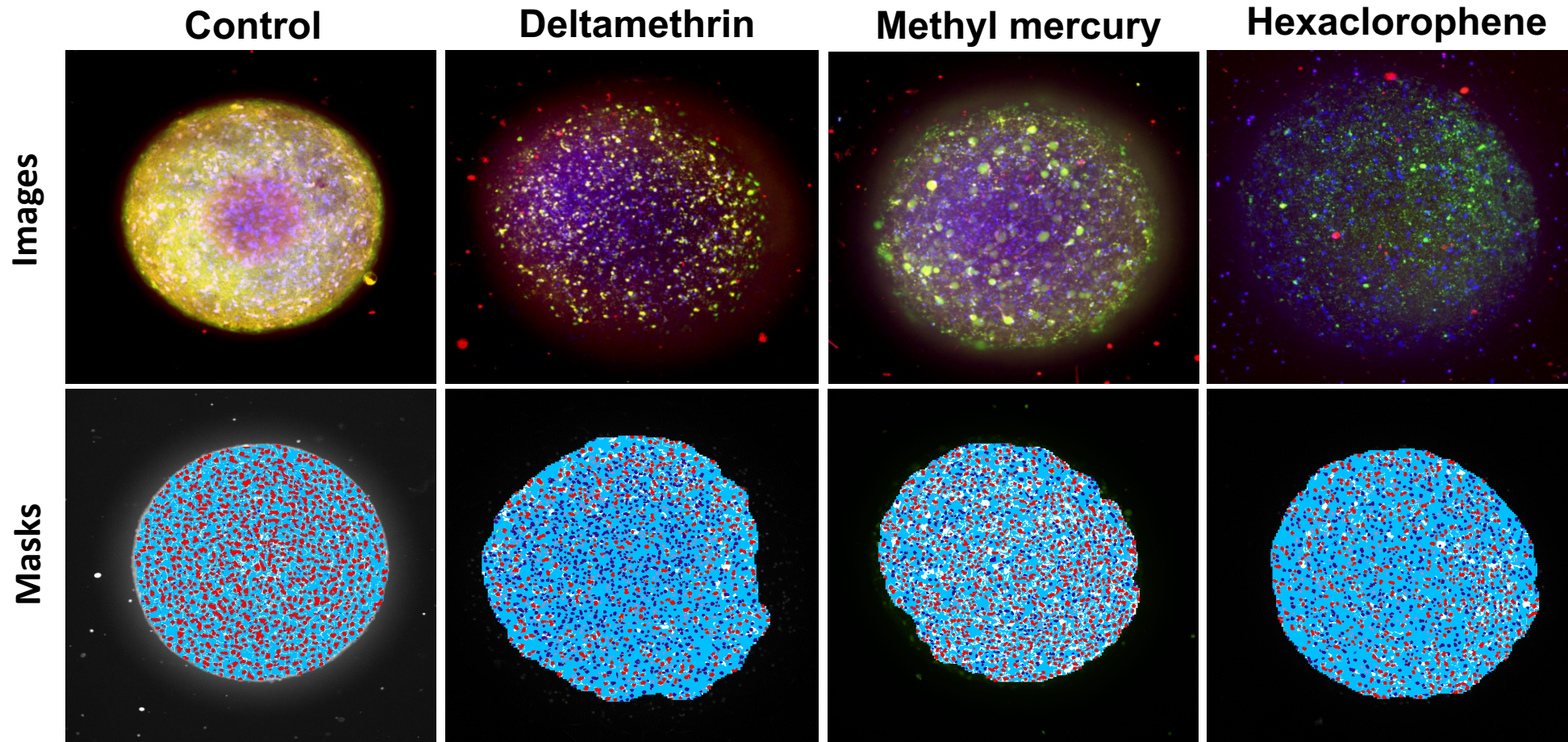
- IC₅₀ determination and comparison with the literature of some control compounds

Compound	Mechanism of Action	IC ₅₀ (μM), This Study	Reference IC ₅₀ (μM)
Kainic acid	Agonist of kainate receptor	2.66	3; 20
MK-801	Antagonist of NMDA receptor	0.033	0.037
CNQX	Antagonist of AMPA/kainate receptor	2.05	0.3–1.5
Muscimol	Agonist of GABA _A receptor	0.021	0.47
(R)-Baclofen	Agonist of GABA _B receptor	0.45	0.70
GABA	Endogenous agonist of GABA receptor	5.93	7; 27
Haloperidol	Antagonist of D2 dopamine receptor	0.13	0.037
Lidocaine	Voltage-gated Na ⁺ channel blocker	9.47	5
Phenytoin (dilantin)	Voltage-gated Na ⁺ channel blocker	9.41	16
Lamotrigine isothionate	Voltage-gated Na ⁺ channel blocker	34.1	66

Sirenko et al. 2018

microBrain 3D is amenable to HCS

- Representative images of Cell Viability quantification using ImageXpress Micro Confocal system



Nuclei (Hoechst): Blue / Viability (Calcein AM): Green / Mitochondria (MitoTracker Orange): Red

Sirenko et al. 2018

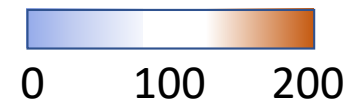
microBrain 3D is amenable to HCS



Legend:

- Drugs: Pharmaceutical compounds (n=17)
- FR: Flame retardants (n=19)
- Industrial (n=15)
- PAH: Polycyclic Aromatic Hydrocarbons (n=20)
- Pesticides (n=17)
- Controls (n=11)

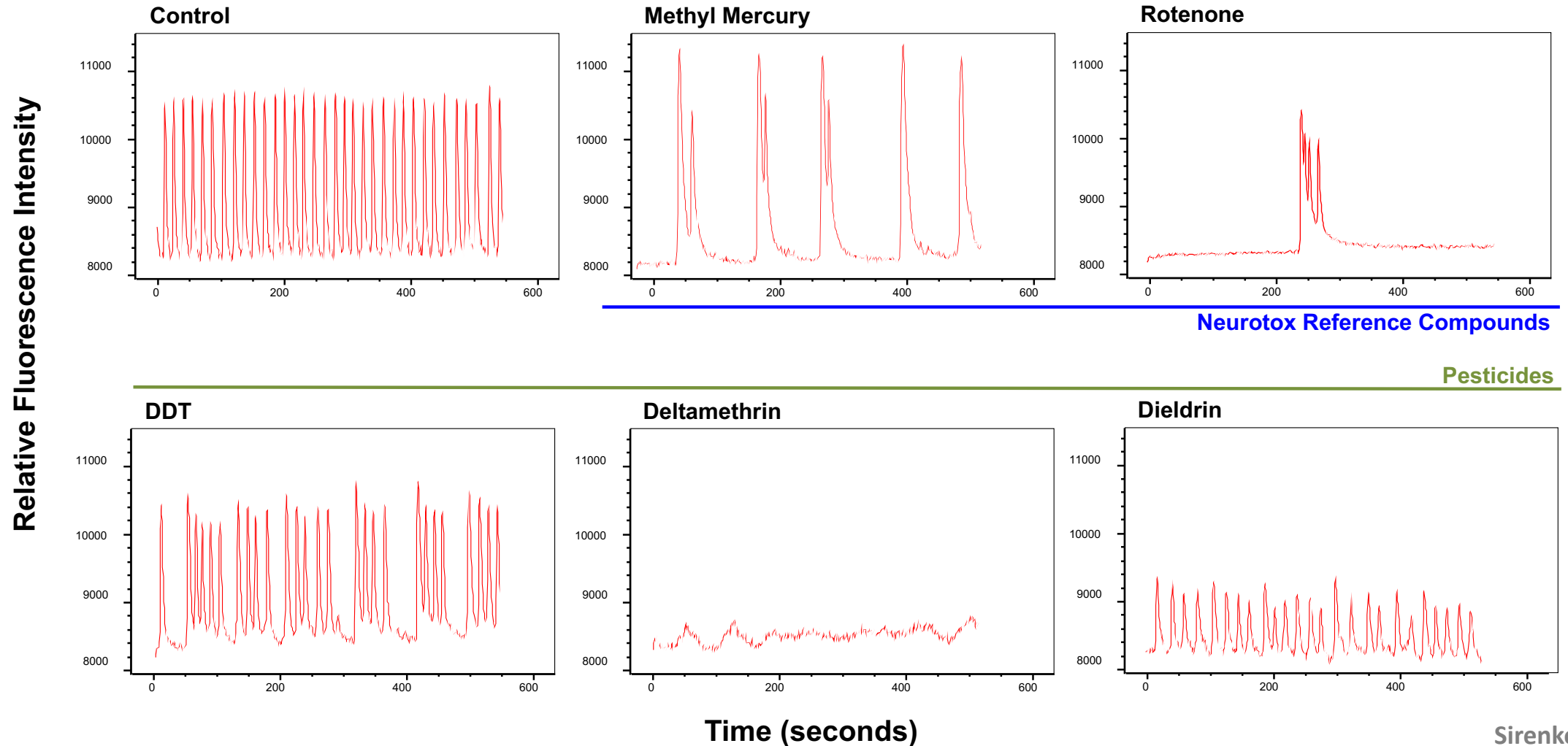
Percentage of Normalized Response



Sirenko et al. 2018

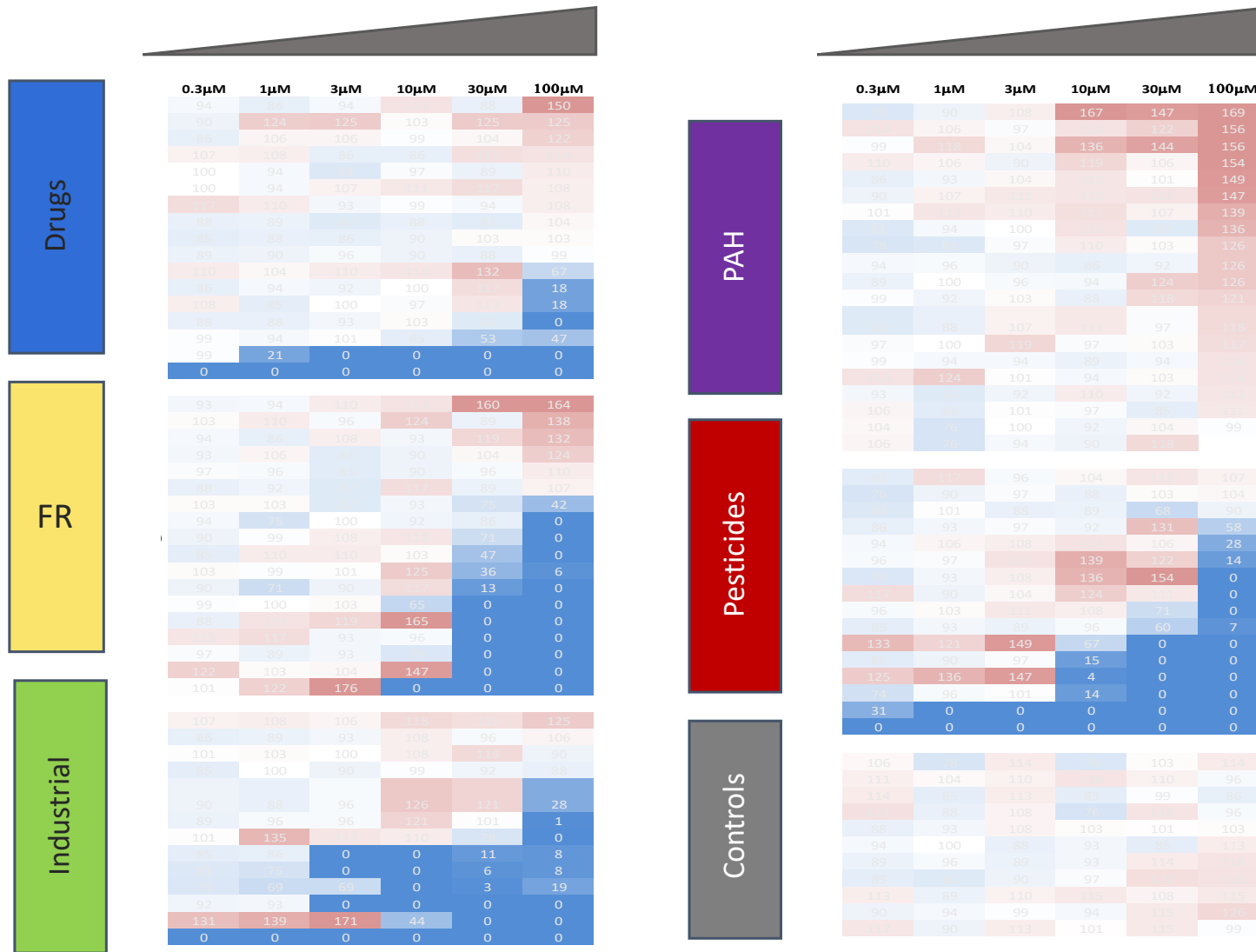
Neurotoxic compounds interfere with microBrain 3D activity

- Representative traces of spontaneous activity recorded on FLIPR Tetra System



Sirenko et al. 2018

Neurotoxic compounds interfere with microBrain 3D activity



Legend:

- Drugs: Pharmaceutical compounds (n=17)
- FR: Flame retardants (n=19)
- Industrial (n=15)
- PAH: Polycyclic Aromatic Hydrocarbons (n=20)
- Pesticides (n=17)
- Controls (n=11)

Percentage of Normalized Response



0 100 200

Sirenko et al. 2018

Conclusions

- microBrain 3D is a very homogenous human-based tool for CNS interrogation *in vitro*.
- The platform was successfully used for HCS and HTS when paired with FLIPR Tetra System and ImageXpress Micro Confocal system.
- A functional phenotype (Ca²⁺ spontaneous activity) was more sensitive to capture toxicity of compounds than cell viability.
- Many compounds investigated on this study presented a detrimental effect on the neuronal activity of microBrain 3D, which could be used as a phenotype to investigate CNS toxicity.

Acknowledgements

Oksana Sirenko, PhD

Carole Crittenden

Krithika Sridhar

Sarah Vargas-Hurlston, PhD

Kristen Ryan, PhD

Oivin Guicherit, PhD

Ryan Gordon, PhD

Fabian Zanella, PhD

Poster:

*Multiplexed Automated Imaging Assays for
Compound Testing Using Induced Pluripotent
Stem Cell-Derived Cells*

Poster W-3210

Date: 06/28/2019

Time: 7pm



<https://bit.ly/2WXNS3W>

For additional information:

StemoniX contact:

David Buché

David.Buche@stemonix.com

612-325-4273

Molecular Devices contact:

info@moldev.com

Read More:

Sirenko, O. et al. Functional and Mechanistic Neurotoxicity Profiling Using Human iPSC-Derived Neural 3D Cultures, *Toxicological Sciences*, Volume 167, Issue 1, January 2019, Pages 58–76,

<https://doi.org/10.1093/toxsci/kfy218>

microBrain® 3D Neurotoxicity Profiling Application Note

<http://bit.ly/2YstGc1>