

Multiplexed automated assays for neurotoxicity evaluation using induced pluripotent stem cell-derived neural 3D cell models

O. SIRENKO¹, C. CRITTENDEN¹, R. DELLE BOVI⁴, E.F. CROMWELL³, B. ANSON², C. CARROMEU²;
¹Molecular Devices, San Jose, CA; ²Stemonix, San Diego, CA; ³Protein Fluidics, Burlingame, CA; ⁴Enzo Life Sciences, Farmingdale, NY

Introduction

Cell-based phenotypic assays have become an increasingly attractive alternative to traditional in vitro and in vivo testing in pharmaceutical drug development and toxicological safety assessment. The effectiveness of automated imaging assays combined with the organotypic nature of human induced pluripotent stem cell (iPSC)-derived cells opens new opportunities to employ physiologically relevant in vitro model systems to improve screening for new drugs or potential chemical toxicities. In our studies, we used human iPSC-derived neural cultures to test functional and morphological end points for toxicity evaluation in a multi-parametric assay format.

For neurotoxicity assessment we employed iPSC-based 3D neural platform composed of mature cortical neurons and astrocytes (Stemonix Co.). The neuronal cells in neurospheroids generated spontaneous synchronized calcium oscillations. We used fast kinetic fluorescence imaging to measure the patterns and frequencies of the Ca²⁺ oscillations. Characterization of oscillation profiles in neurospheroids was performed through multi-parametric analysis presenting more than 15 read-outs including frequency, amplitude, characterized primary and secondary peaks, peak width, and waveform irregularities. In addition, cellular and mitochondrial toxicities were assessed by high-content imaging. The assay was optimized for high throughput screening in 384-well plates and displayed a highly consistent performance. We evaluated neuroactive and neurotoxic profiles of more than 120 compounds, including set of neuromodulators with known mechanisms of action, set of various pharmaceutical drugs, sample library of cannabinoids, and a library of neurotoxic chemicals including flame retardants, pesticides, and poly-aromatic hydrocarbons.

Our results show that the iPSC-derived 3D neurospheroid assay platform is a promising biologically-relevant tool to assess the neuroactive and neurotoxic potential of pharmaceutical drugs and environmental toxicants.

Stemonix microBrain 3D spheroids

The Stemonix® microBrain® 3D Assay Ready Platform is a high-throughput 3D culture platform that more closely resembles the tissue development and constitution of native human brain tissue. In this platform, human iPSC-derived neuronal spheroids, approx. 600 µm in diameter, are composed of a physiologically relevant co-culture of functionally active cortical glutamatergic and GABAergic neurons (identified by MAP2; green) and astrocytes (identified by GFAP; red). This balanced cellular mix allows the development of a neural network enriched in synapses, creating a highly functional neuronal circuitry. The neuronal cells in the microBrain 3D spheroids are physiologically active, with spontaneous synchronized, readily detectable calcium oscillations.

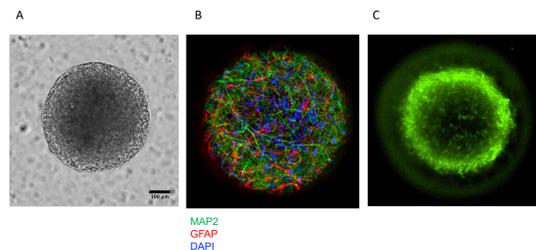


Figure 1. A. Human iPSC-derived neural spheroids composed of a co-culture of active cortical neurons B. (identified by MAP2; green) and astrocytes (identified by GFAP; red), approximately 600 µm diameter, imaged with ImageXpress® Micro Confocal system, 20X magnification. C. Stained with Calcium dye.

Instrument

- The FLIPR® Penta System is powered by a new high-speed EMCCD camera and the new ScreenWorks® Peak Pro™ 2 software.
- The system allows measurement and analysis of complex patterns of calcium oscillations in human iPSC-derived cardiomyocytes and neurons.



We used the new high speed EMCCD camera on the FLIPR Penta system to measure the fast kinetic patterns and frequencies of Ca²⁺ oscillations of neuro-spheroids as monitored by changes in intracellular Ca²⁺ levels with FLIPR® Calcium 6 Assay Kit (Molecular Devices). The instrument equipped with the new ScreenWorks Peak Pro 2 peak analysis software module allows analysis and characterization of the primary and secondary peaks and complex oscillation patterns.

Methods

3D neural cultures

microBrain 3D Assay Ready 384-Well plates were obtained from Stemonix, Inc. Plates were shipped pre-plated under ambient conditions. Each well contained a single, uniformly sized human iPSC-derived cortical neural spheroid matured 8-9 weeks. The human spheroids were exposed to compounds for 24 hours, or as indicated in the figures.

Calcium Flux Assay

The intracellular Ca²⁺ oscillations were assessed using the FLIPR Calcium 6 Assay Kit according to the regular protocol; spheroids were loaded with dye for 2 hours before measurements.

Cell Staining

To assess phenotypic changes, cells were stained live using a mixture of three dyes: the viability dye Calcein AM (1 µM), the mitochondria potential dye MitoTracker Orange (0.2 µM), and the Hoechst nuclear dye (2 µM). Hoechst 33258 (ultra pure) (ENZ-52402) Calcein AM (ultra pure) (ENZ-52002), MITO-ID® Membrane potential detection kit (ENZ-51018-K100)

Results

Calcium Oscillations Evaluated by the FLIPR Penta System

The neuronal cells in the microBrain 3D spheroids generate spontaneous synchronized calcium oscillations. We used fast kinetic fluorescence imaging on the FLIPR Penta system to measure the patterns and frequencies of the Ca²⁺ oscillations of neuro-spheroids as monitored by changes in intracellular Ca²⁺ levels with FLIPR Calcium 6 Assay Kit. A set of known neuromodulators was tested, including agonists and antagonists of NMDA, GABA and AMPA receptors; kainic acid, analgesic, and anti-epileptic drugs.

Analysis of Kinetic Patterns

Advanced analysis methods implemented in the ScreenWorks Peak Pro 2 software module provide multi-parametric characterization of the Ca²⁺ flux oscillation patterns. This phenotypic assay allows for the characterization of readouts such as oscillation frequency, amplitude, peak width, secondary peaks, and amplitude and spacing irregularity. The effects of modulators of neuronal activity were evaluated by measuring changes in multiple read-outs.

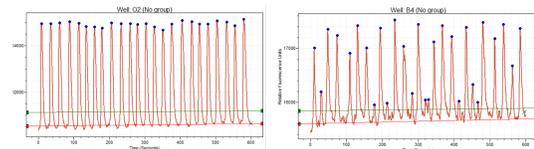


Figure 2. The assay was used for high-throughput assessment of neurotoxicity effects of chemicals in vitro which would help to evaluate and prioritize different drugs and substances for further testing. ScreenWorks PeakPro2 software allows more than 20 descriptors of waveforms. We have evaluated multi-parametric approach for description of changes in oscillations.

Effects of Neuromodulators on Calcium Oscillations

A number of known modulators of neuronal activity were assayed at different time-points and the calculated EC₅₀ values for compound effects. Changes were observed as inhibitions or activations of the peak frequency or other measurements, corresponding to the expected effect of the tested neuromodulator.

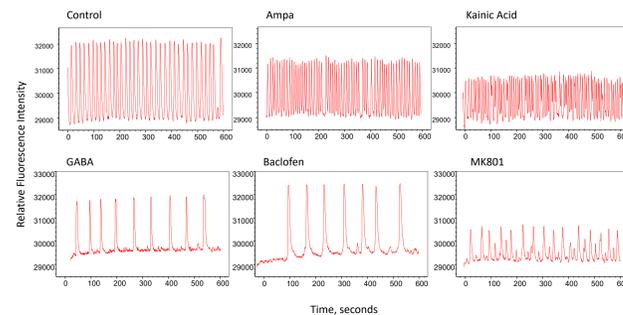


Figure 3. Using high speed EMCCD fluorescence imaging (FLIPR Penta system), we have tested the impact of various compounds on the rates and patterns of intracellular Ca²⁺ oscillations using calcium-sensitive dye (FLIPR Calcium 6 Assay Kit). High throughput recording and analysis of the entire 384-well plate was simultaneously performed using the FLIPR Penta system. Representative traces of calcium oscillations are shown for the control and compound treated spheroids. Spheroids were loaded with dye for 2h and then treated with compounds for 30min. Concentration-dependent changes in the patterns were observed. Oscillation traces were recorded by the FLIPR system for 10 min. Patterns were characterized using ScreenWorks Peak Pro 2 software. Representative traces of calcium oscillations are shown for 1µM concentrations, 30min treatment. ¹

Effects of Environmental Neurotoxicants

A library of environmental compounds containing known and unknown neurotoxicants was tested in the assay. Large number of compounds (57 out of 80) demonstrated deviations from the control pattern and were flagged in the assay. Effective concentrations for compounds effect were determined for various read-outs demonstrating sensitivity of the method for HTS in vitro assay for potential neurotoxicity.

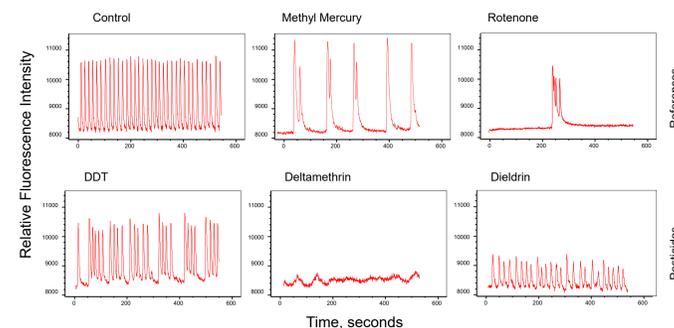


Figure 4. Modulations of calcium oscillation patterns for selected neurotoxic compounds. Detailed information about screen and effective concentrations can be found at ToxSci. ¹ Sirenko et al., Toxicol Sci 2019 Jan 1;167(1):58-76.

Results, continued

Effects of Neuro-active compounds

As a next step, we tested a selected set of neuro-active drugs that have expected effects on the nervous system, including established drugs used for treatment of neurological disorders as well as substances known to cause psychotic effects.

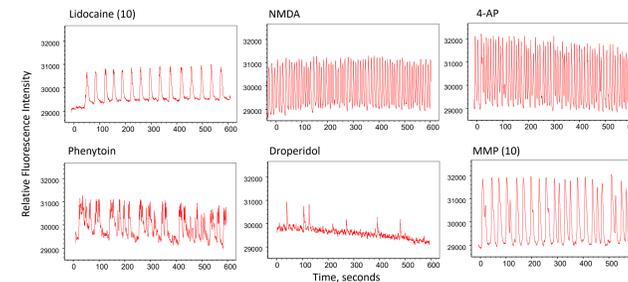


Figure 5. Representative traces of calcium oscillations are shown for 3µM concentrations, 30min treatment.

Effects of Neurotoxic Drugs

To test if the assay can identify chemicals for their potential neurotoxic effects, we tested a panel of 23 compounds that represent different classes of drugs, known to cause peripheral neuropathy (cancer drugs), confusion (selected antibiotics), hearing loss (gentamycin), or other types of disorders. For neurotoxicity test neurospheroids were treated with compounds for 24h. Selected drugs caused mostly moderate perturbations of calcium oscillation patterns.

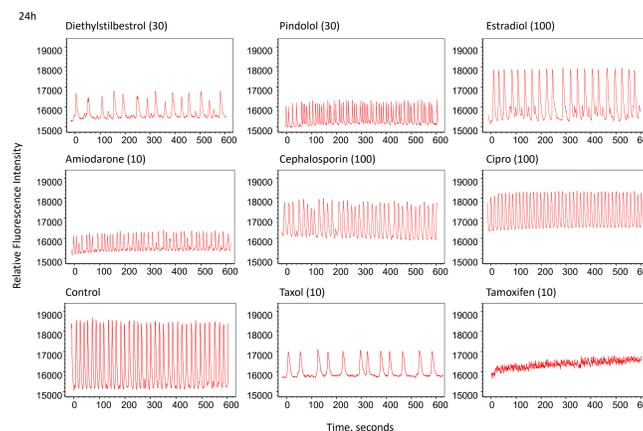


Figure 6. Representative traces of calcium oscillations are shown for selected anti-cancer drugs and antibiotics. Concentrations (µM), 24h treatment.

Multi-parametric assessment of compound effects

The assay shows promise for high-throughput assessment of neurotoxicity effects of chemicals in vitro. ScreenWorks PeakPro2 software allows more than 20 descriptors of waveforms. We have evaluated multi-parametric approach for description of changes in oscillations.

Compounds	Peak Count	Amplitude	Oscillation Frequency	Peak-width	Oscillation Amplitude	Peak-spacing	Secondary Peaks	Irregularity	Peak* ¹	Secondary Peak*	Cycloactivity*	Mitochondria Intensity	Max Intensity	Drug Action
1. Amla	0.2	0.48	up	3	10	1	0.3	0.3	0.3			100	100	AMPA agonist
2. Kainic acid	2.2	2.2	up/down	3	10	1	0.3	0.3	0.3			100	100	Kainate receptor agonist
3. NBQX	0.2	0.38	down	3	10	1	0.3	0.3	0.3			100	100	NMDA antagonist, Channel Blocker
4. CNQX	10.2	down		15	15	15	15	15	15			100	100	AMPA/kainate antagonist
5. Gabazine	0.3	0.3	down	0.3	0.3	0.3	0.3	0.3	0.3			100	100	GABA A agonist
6. Bicuculline	0.2	down		3	3	3	3	3	3			100	100	GABA B agonist
7. GABA	-15	3.56	down									100	100	Engorgement agonist
8. Gabazine acid	1.0	up										1000	1000	Non-specific agonist
9. Flunitrazepam	-2	-1	up	3	10	3	3	3	3			100	100	NMDA 2B agonist
10. MK801	3.56	-5	up	10	3							20	20	NMDA agonist, Non-Specific
11. DMSO	-1	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	100	100	Control
12. DMSO 1%	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	100	100	Control

Compound	Peak Count	Amplitude	Oscillation Frequency	Peak-width	Oscillation Amplitude	Peak-spacing	Secondary Peaks	Irregularity	Peak* ¹	Secondary Peak*	Cycloactivity*	Mitochondria Intensity	Max Intensity	Drug Action
1. Lidocaine	2.5	0.71	down	3	3	3	3	3	3			100	100	Anti-epileptic
2. Lidocaine	0.8	5.7	down	30	30	30	30	30	30			100	100	Anti-epileptic
3. Chlorpromazine	30.7	30.3	down	10	30	10	10	30	30			100	100	Anti-epileptic
4. Phenytoin	124.8	-100	down	100	10	30	10	100	100			100	100	Anti-epileptic
5. Lamotrigine	-50	no effect	up	30	30	30	30	30	30			100	100	Anti-epileptic
6. Carbamazepine	1.4	no effect	up	30	30	30	30	30	30			100	100	Stimulant
7. Valproic acid	0.128	-1	down	1	0.03	0.03	0.03	3	3			100	100	Anticonvulsant, anti-epileptic
8. MMP	59.5	no effect	down	100	100	100	100	100	100			100	100	Stimulant
9. Nitroic Acid Na Salt	1.0	no effect	down	100	100	100	100	100	100			100	100	Anti-epileptic
10. Propofol	29.5	1.28	up	none	none	none	none	none	none			100	100	Stimulant, antiseptic
11. Propofol	49.2	4.6	up	10	10	10	10	10	10			100	100	Stimulant
12. Acetanilopam	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	100	100	Negative control

Compounds	Peak Count	Amplitude	Oscillation Frequency	Peak-width	Oscillation Amplitude	Peak-spacing	Secondary Peaks	Irregularity	Peak* ¹	Secondary Peak*	Cycloactivity*	Mitochondria Intensity	Max Intensity	Drug Action
1. Pindolol	57.3	30	down	100	100	100	100	100	100			100	100	Beta-blocker
2. Diethylstilbestrol	>100	39.3	down	100	100	100	100	100	100			100	100	Estrogen agonist
3. Tamoxifen	4.48	3.58	down	100	100	100	100	100	100			100	100	Estrogen receptor modulator
4. Taxol	9.15	100	down	30	10	3	3	100	100			100	100	Anti-cancer
5. Chlorpromazine	100	100	down	100	100	100	100	100	100			100	100	Antipsychotic
6. Ciprofloxacin	100	100	down	100	100	100	100	100	100			100	100	Antibiotic
7. Cephalosporin	100	100	down	100	100	100	100	100	100			100	100	Antibiotic
8. Gentamicin sulfate	100	100	down	100	100	100	100	100	100			100	100	Antibiotic
9. Ibuprofen	100	100	down	100	100	100	100	100	100			100	100	Anti-inflammatory
10. Ibuprofen	100	100	down	100	100	100	100	100	100			100	100	Anti-inflammatory
11. Estradiol	202	down		30	10	30	30	30	30			100	100	Hormone, oestrogen
12. Amitriptyline	8	1.48	up/down	30	3	10	10	100	100			100	100	Anti-arrhythmic drug
13. Oxcarbazepine	100	100	down	100	100	100	100	100	100			100	100	Anti-epileptic
14. Propylthiouracil	100	100	down	100	100	100	100	100	100			1000	1000	Regulatory stimulant
15. Propylthiouracil	100	100	down	100	100	100	100	100	100			100	100	Glucocorticoid
16. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
17. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
18. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
19. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
20. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
21. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
22. Hydroxypropyl-beta-cyclodextrin	100	100	down	100	100	100	100	100	100			100	100	Neurostimulant
23. Carbamazepine	100	100	up/down	100	100	100	100	100	100			100	100	Anti-epileptic drug
24. Amoxicillin	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	no effect	100	100	Negative control

Table 1. EC₅₀ values or lowest effective concentrations (µM) that cause specific changes (*) are indicated for different read-outs. No changes indicated with blank cells.

Effects of Cannabinoids

The assay was used for testing neuroactive effects of the panel of cannabinoid substances (from Enzo Life Sciences). Several substances were active in the assay.

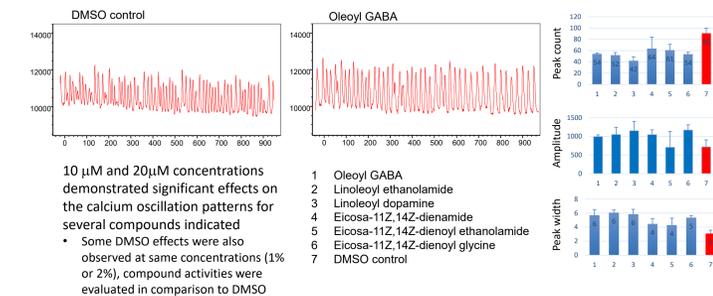


Figure 7. Calcium oscillation traces and measurements are shown for 10µM concentrations, 30min treatment. SCREEN-WELL® Endocannabinoid library (BML-2801-0100).
 10 µM and 20µM concentrations demonstrated significant effects on the calcium oscillation patterns for several compounds indicated
 • Some DMSO effects were also observed at same concentrations (1% or 2%), compound activities were evaluated in comparison to DMSO

Assessment of Viability by High Content Imaging

Confocal imaging and 3D image analysis methods were used to characterize compound effects on the morphology and viability of 3D neural spheroids. To evaluate cytotoxicity effects, cells were treated with various compounds for 24h, and then, live cells were stained with Hoechst nuclear stain, Calcein AM, and MitoTracker Orange dyes. Images were acquired using the ImageXpress Micro Confocal system, using the confocal option and 3D imaging. Projection images were then analyzed using the Custom Module Editor and Cell Scoring algorithms for detection of cell numbers for all cells, live cells (Calcein AM positive cells), and cells with intact mitochondria (MitoTracker positive cells). The analysis methods provide efficient tools for characterization of cell and spheroid morphology.

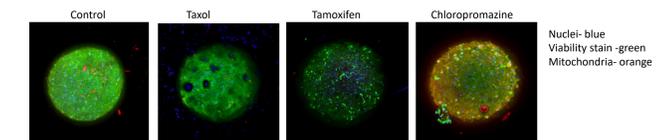
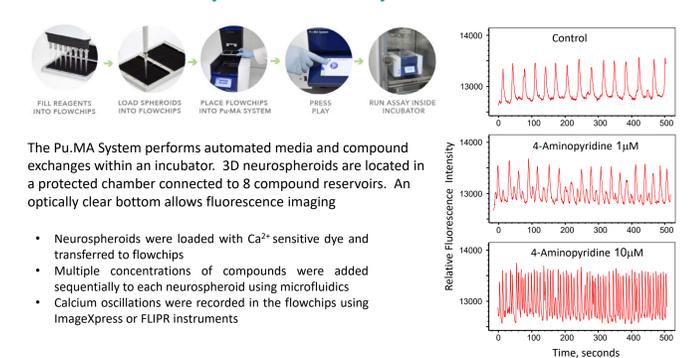


Figure 8. Composite projection images of neural spheroids. ImageXpress Micro Confocal system. Color balance reflects decreases in Calcein AM or MitoTracker signal intensities.

Automated Assay with Pu-MA System



The Pu-MA System performs automated media and compound exchanges within an incubator. 3D neurospheroids are located in a protected chamber connected to 8 compound reservoirs. An optically clear bottom allows fluorescence imaging

- Neurospheroids were loaded with Ca²⁺ sensitive dye and transferred to flowchips
- Multiple concentrations of compounds were added sequentially to each neurospheroid using microfluidics
- Calcium oscillations were