

# Multiplexed structure-function assay for high throughput drug safety testing on human induced pluripotent stem cell-derived cardiomyocytes and neurons

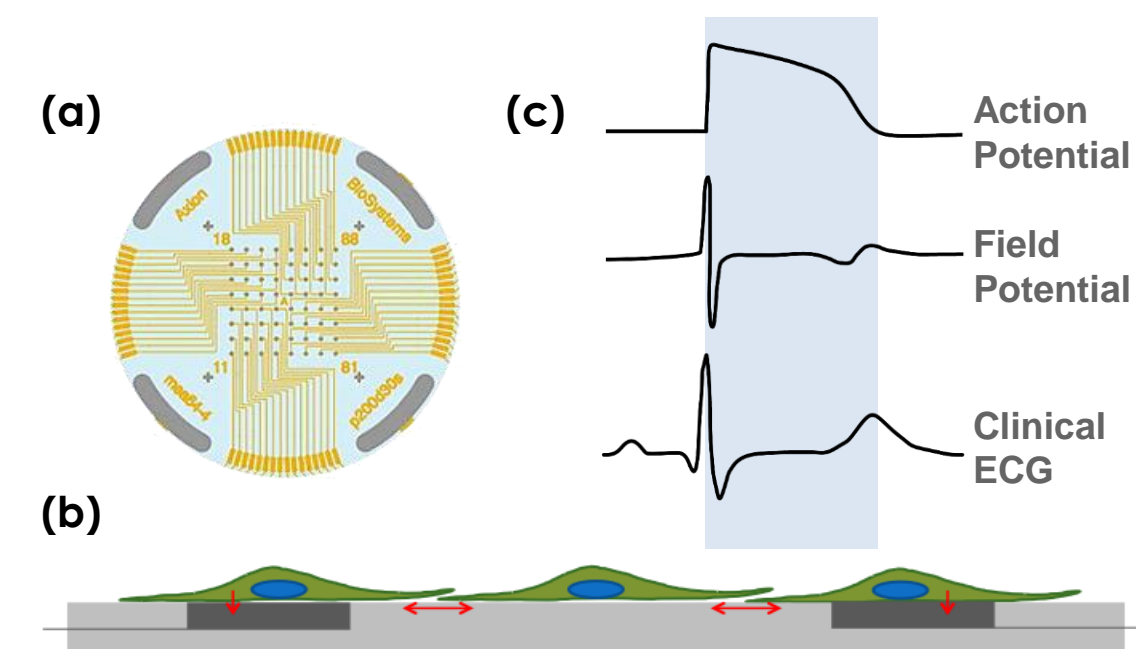
Daniel Millard, Denise Sullivan, Heather Hayes  
Axion BioSystems, Atlanta, GA

## Multiwell MEA Technology

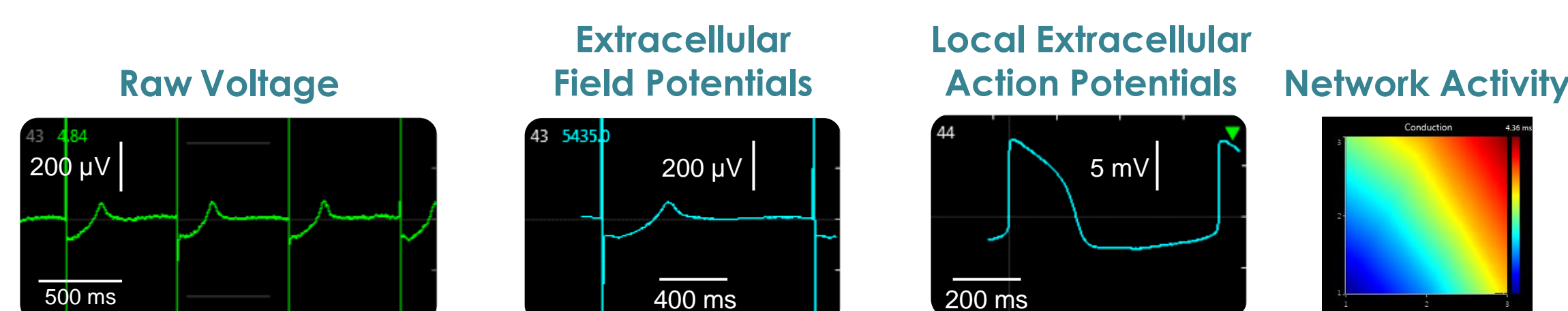
### Microelectrode array technology

The flexibility and accessibility of neural and cardiac *in vitro* models, particularly induced pluripotent stem cell (iPSC) technology, has allowed complex human biology to be recapitulated *in vitro*. Accurate characterization of the function and structure of neurons and cardiomyocytes are required for a comprehensive drug safety assay.

Axion BioSystems' Maestro™ multiwell microelectrode array (MEA) platform is a benchtop system for non-invasive functional and structural characterization of cellular networks cultured on a dense array of extracellular electrodes in each well. The Maestro records functional activity and tracks cell coverage and viability from the same microelectrodes in one simple assay.

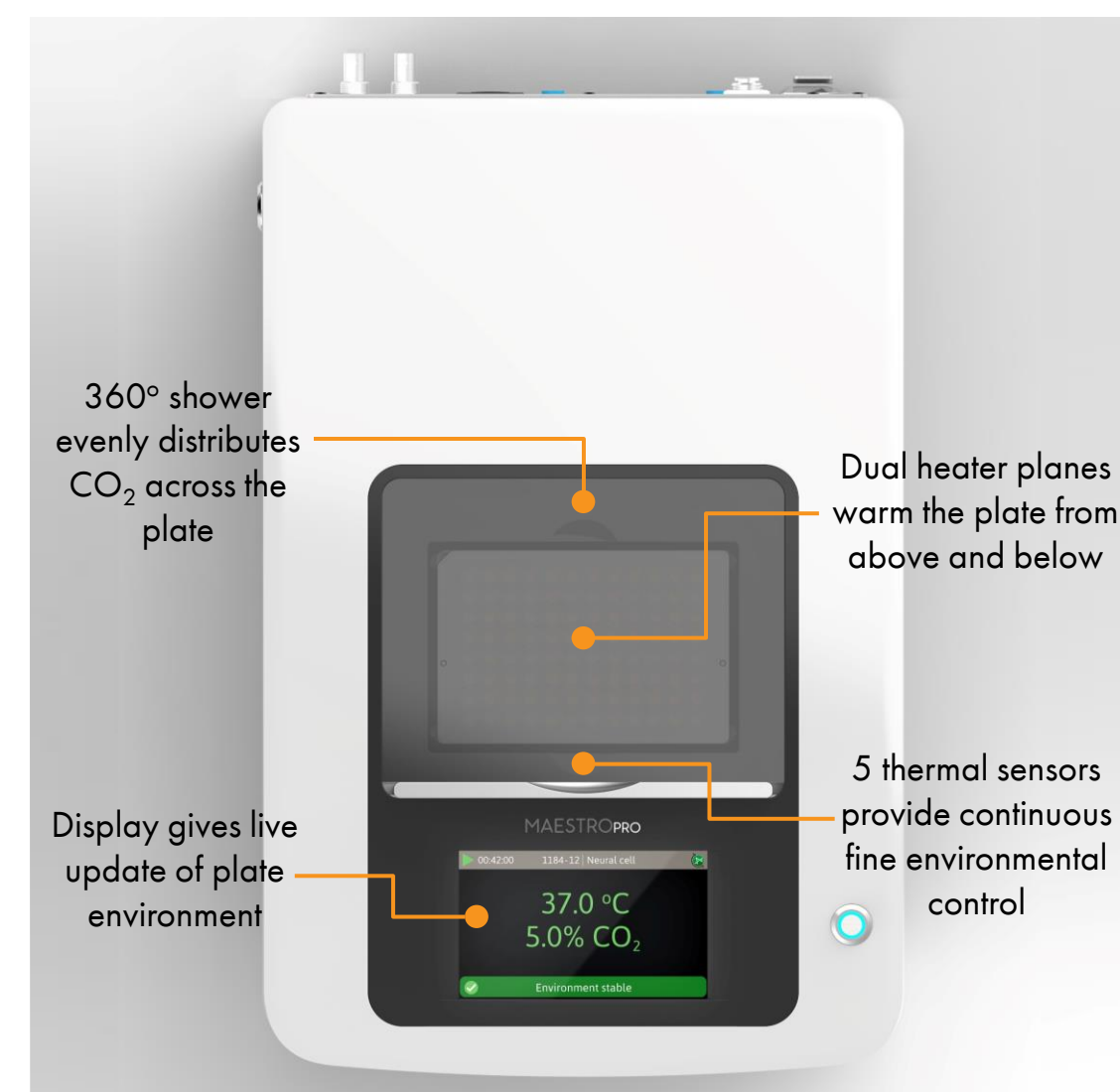


A planar grid of microelectrodes (a) interfaces with cultured neurons or cardiomyocytes (b) to model complex, human systems. Electrodes detect changes in raw voltage and record extracellular field potentials (c). The same microelectrodes can track cell viability using impedance measurements.



Raw voltage signals are processed in real-time to obtain extracellular field potentials from across the network, providing a valuable electrophysiological phenotype for applications in drug discovery, toxicological and safety screening, disease models, and stem cell characterization

## Introducing the Maestro Pro™ and Maestro Edge™



- **Label-free, non-invasive recording** of extracellular voltage from cultured electro-active cells
- **Integrated environmental control** provides a stable benchtop environment for short- and long-term toxicity studies
- **Fast data collection rate (12.5 KHz)** accurately quantifies the depolarization waveform
- **Sensitive voltage resolution** detects subtle extracellular action potential events
- **Industry-leading array density** provides high quality data from across the entire culture
- **Scalable format (6-, 24-, 48- and 96-well plates)** meets all throughput needs on a single system
- **State-of-the-art electrode processing chip (BioCore v4)** offers stronger signals, ultra-low frequency content, and enhanced flexibility



Feature	Maestro Edge	Maestro Pro
Recording Electrodes	384	768
BioCore Chip	6 Chips (v4)	12 Chips (v4)
MEA Plates	24-Well	6-, 24-, 48-, 96-Well
Integrated Hard Drive	0.5 TB	1.0 TB
Touchscreen	No	Yes
Optical Stimulation	No	Yes

The Maestro Pro™ (left) and Maestro Edge™ (right) offer the latest MEA technology for optimal data

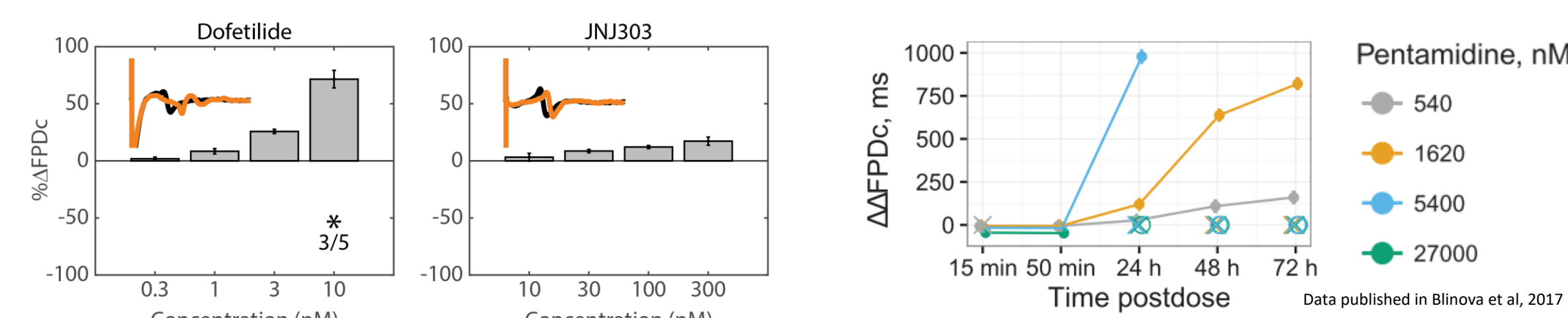
## MEA Assay with Cardiomyocytes

### Structure and Function in One Assay

The Maestro MEA platform enables assessment of *in vitro* cardiomyocyte function and structural integrity with an easy-to-use benchtop system. The Maestro detects electrical signals from cells cultured onto an array of planar electrodes in each well of the MEA plate, with five modes providing critical safety information:

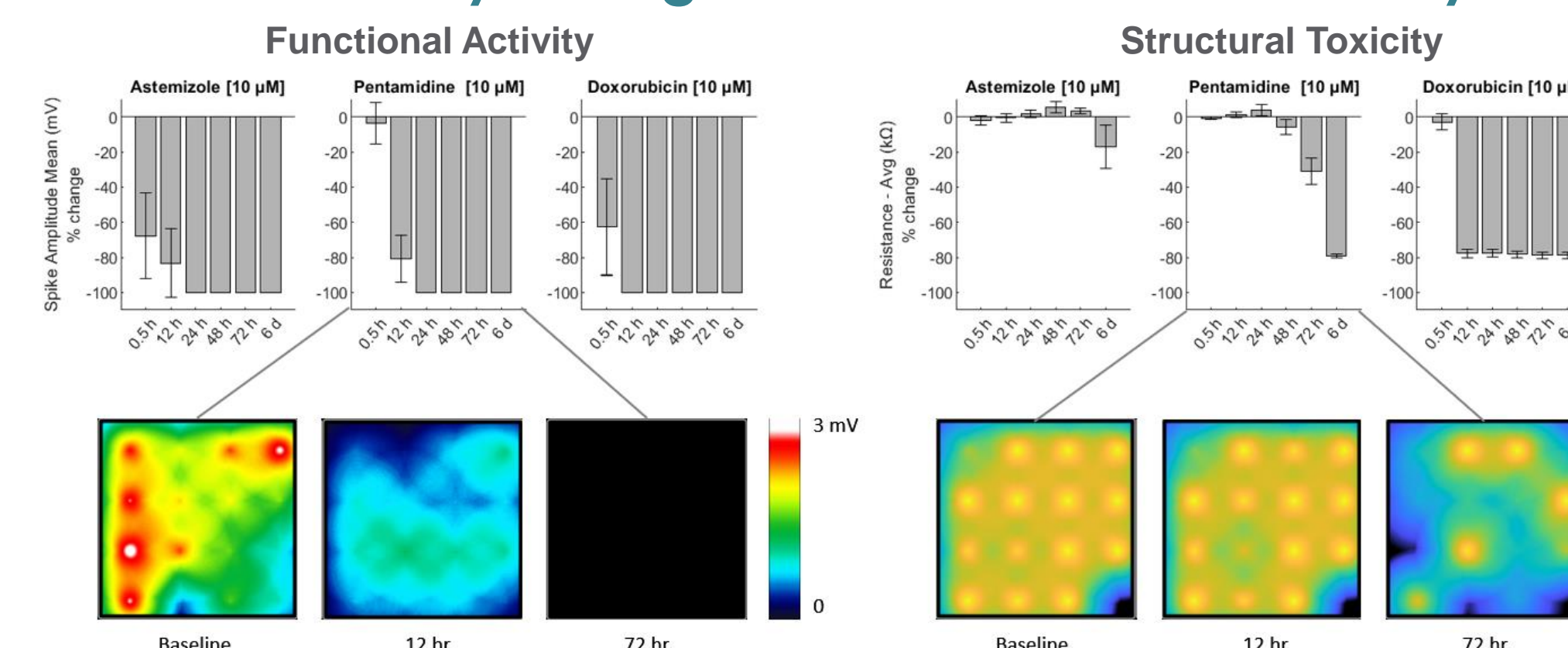
- **Field Potential** – “gold standard” measurement for multiwell cardiac electrophysiology.
- **LEAP** – first truly scalable technique for acquiring action potential signals from intact cardiac monolayers.
- **Conduction** – detect speed and direction of action potential propagation.
- **Contractility** – assess the contractility of cardiac monolayers adhered to the 2D array.
- **MEA Viability** – track cell coverage and viability using impedance technology.

### Field Potential Provides Sensitive, Label-Free Measurements of Acute and Chronic Function



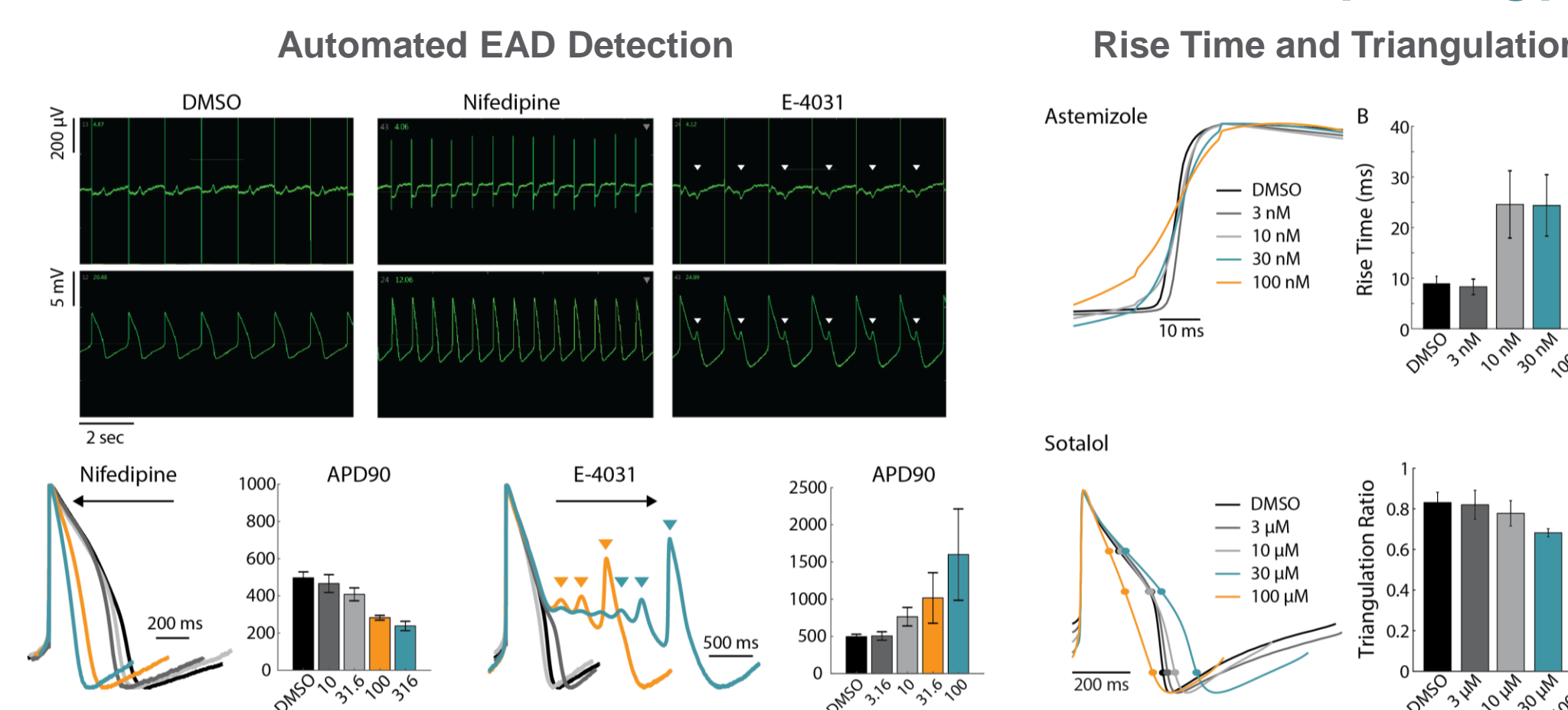
Field potential metrics provide the highest sensitivity for small effects on repolarization across acute and chronic timescales. (Left) Dofetilide caused significant prolongation of FPDc and arrhythmias, whereas JNJ303 produced a more subtle prolongation of FPDc. (Right) A chronic field potential assay detected dose-dependent effects of Pentamidine due to inhibition of hERG trafficking.

### MEA Viability Distinguishes Cardiotoxins from Cytotoxins



Structure and function were tracked over days using field potential and impedance-based viability recordings. All compounds impacted cardiac function by reducing sodium spike amplitude, but only Pentamidine and Doxorubicin induced structural toxicity. Doxorubicin caused cell death within 24 hours, whereas Pentamidine caused cell death over 6 days. In the maps below, brighter colors indicate higher spike amplitudes or cell coverage; darker colors indicate lower spike amplitudes or cell death.

### LEAP Provides Measures of Action Potential Morphology

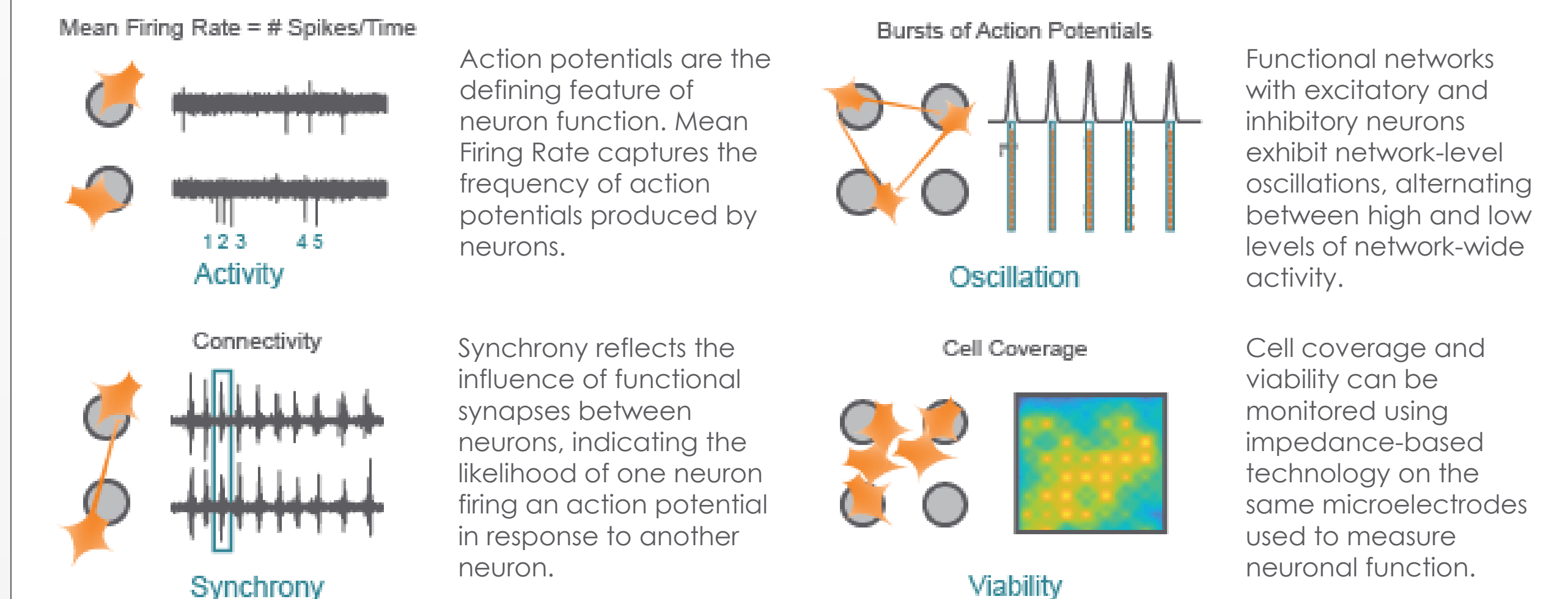


The LEAP signal provides endpoints of action potential morphology, with completely automated analysis of repolarization and EADs. Dose-dependent trends in repolarization timing were detected for Nifedipine (shortening) and E-4031 (prolongation). LEAP also affords measures of rise time and triangulation, providing additional information beyond classical field potential assays.

## MEA Assay with Neurons

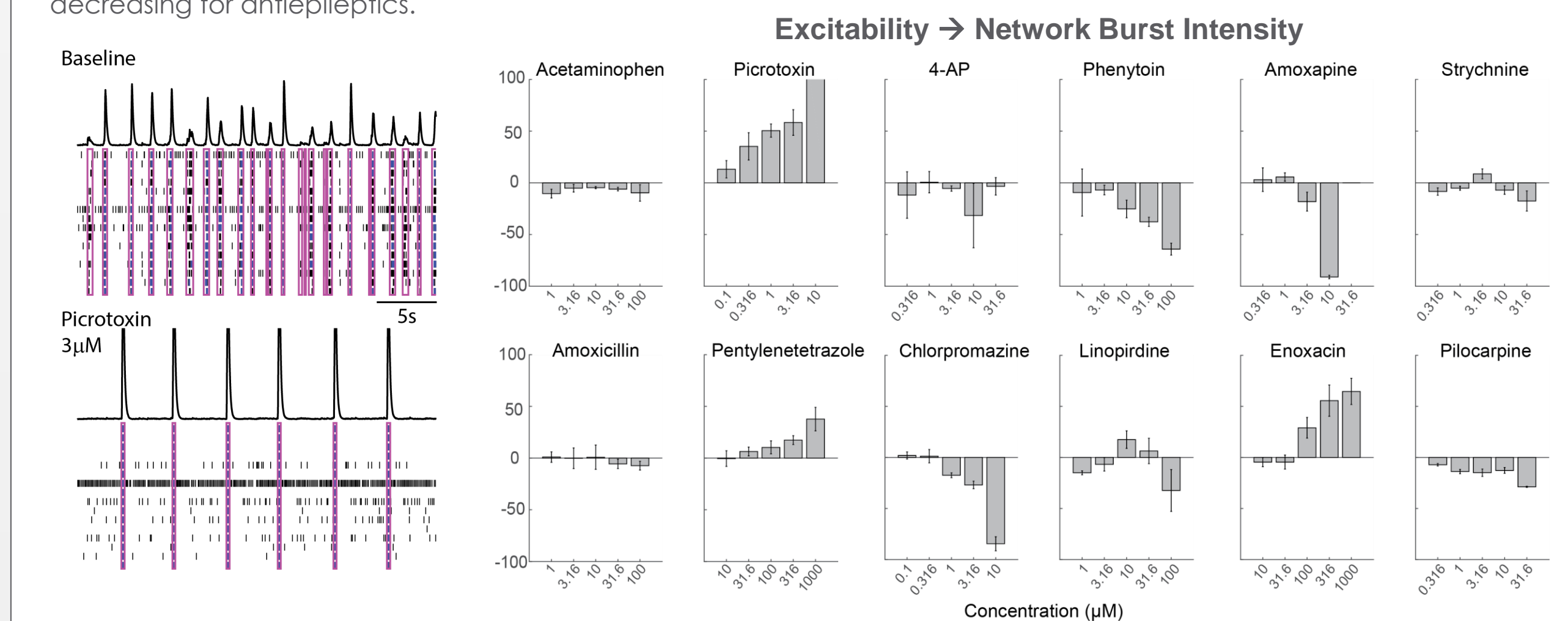
### Structure and Function in One Assay

The Maestro provides a comprehensive assessment of neuronal activity, network connectivity, and structural integrity.



### Network Electrophysiology Assays for Proconvulsant Assessment

As part of the NeuTox consortium (HESI), rodent cortical neurons (Thermo Fisher Scientific) were seeded on CytoView MEA 48 well plates. At DIV28, neurons were dosed with 12 compounds at 5 doses. Network burst intensity, measured as the number of spikes per burst, changed for neuroactive compounds, increasing for most proconvulsants and decreasing for antiepileptics.



### MEA Viability Quantifies Dose-Dependent Cytotoxicity

Many neuroactive compounds, such as antiepileptics and cytotoxins, can cause activity to shutdown, especially at higher doses. Measures of both cell function and viability are required to distinguish compounds that silence neural activity from those that induce cell death. Below, hiPSC-derived neurons (NeuCyte) were dosed with a variety of cytotoxins. Impedance-based MEA Viability was used to monitor cytotoxicity for 72 hrs. Because impedance is non-invasive and label-free, both function and viability can be measured repeatedly without interfering with the biology.

