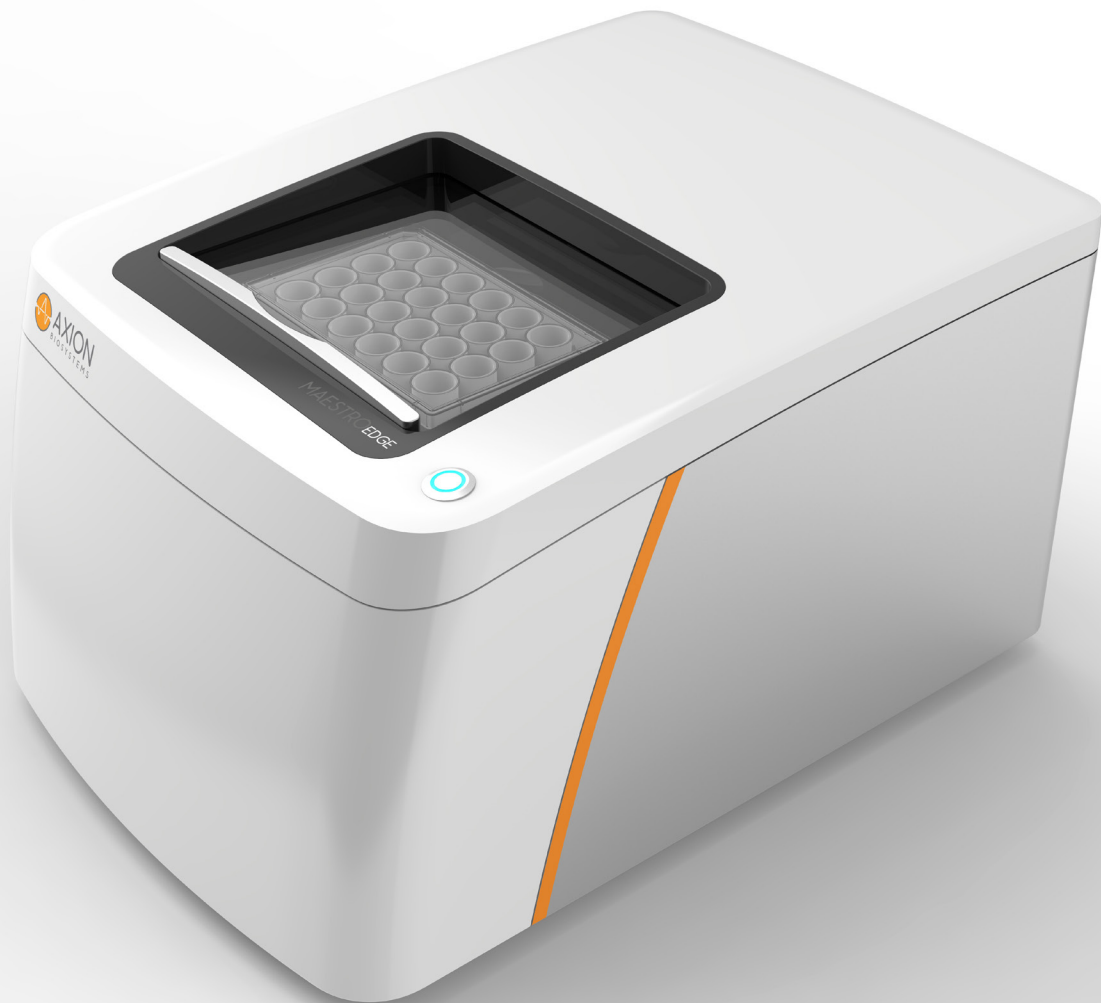


# MAESTRO EDGE

A versatile bioelectronic assay system



# IMAGINE

## A simple assay...

Bioelectronic assays monitor *in vitro* cell health and behavior. Sensitive electrodes track cellular activity, while the culture remains undisturbed.



## Flexible

Noninvasively measure activity anytime, as often as needed, while cells remain in an optimally controlled environment.

## Easy

Use basic cell culture techniques to perform quantitative assays that deliver high-resolution, functional data – with no need for dyes, labels, or complicated steps.

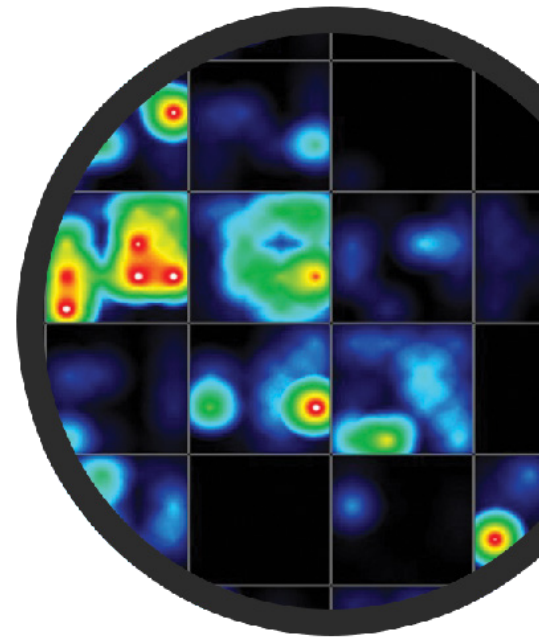
## Time-saving

Measure with the push of a button or schedule recordings ahead of time. No incubation steps or complex handling required with a mostly hands-free assay.

# EXPLORE

## Complex biology...

The versatility of the Maestro Edge makes it **ideal for any lab**. Monitor dynamic cellular activity and perform complex functional experiments with only basic cell culture.



## Electrical *Activity*

Create detailed functional profiles of electrically active cells using microelectrode array (MEA) technology. Repeatedly measure from the same culture and track network development and activity.

- Neurons
- Cardiomyocytes
- Skeletal muscles
- Retinal cells
- Primary or stem cells
- 2D or 3D cultures

## Cellular *Kinetics*

Monitor the health and behavior of any cell type using impedance-based technology. Bioelectrodes measure complex dynamic responses, capturing every minute to reveal detailed cellular kinetics.

- Immuno-oncology
- Cytotoxicity
- Cell proliferation
- Barrier function
- Migration/invasion
- Cell signaling

# DISCOVER

## The Maestro Edge

The latest technological advances to facilitate your research. **The Maestro Edge features:**

### *Dynamic responses*

Detect key parameters of neural activity, network dynamics, and cardiac functionality, measure cell growth and cytolysis, barrier function, and signaling – all label-free and in real time.

### *Push-button acquisition*

Simplify processes with Maestro's automatically adjusted temperature and CO<sub>2</sub> levels and integrated barcode scanner that conveniently tracks plate usage.

### *Precise cellular control*

Recreate specific patterns of cellular activity using electrical stimulation or light pulses (with the Lumos™ optical stimulation system).

### *Integrated environment*

Easily control temperature and CO<sub>2</sub> levels while suppressing electrical noise and mechanical vibrations with Maestro's smart environmental chamber. No need to take up incubator space.

### *Efficient design*

The Maestro Edge is designed to save time, cost, and space. Record from 6 or 24 wells of MEA data, or 96 wells of impedance data.

### *On-the-go connectivity*

Use the Impedance module to track changes in cell proliferation, viability, and cell death from any mobile device. No need to be in the lab.



# Customize your MAESTRO EDGE



## Standard Equipment

- Environmental control
- Push-button start
- Computer
- Analysis software
- Warranty

## Software



Cardiac



Neural



MEA Viability



MEA Automation



Impedance



GxP Impedance



Impedance Automation

## Throughput

### MEA:

6- or 24-well

### Impedance:

96-well

## Lumos Optogenetic Stimulation



24-well, 4 LED colors

## Consumables



BioCircuit MEA



CytoView MEA



Lumos MEA



CytoView-Z

# SOFTWARE MODULES

## To expand your assay

The Maestro Edge platform is available with seven software modules.  
Select the software modules to **match your assay needs**:



**Neural** - Measure electrical network behavior of neurons, including: activity, synchrony, and network oscillations, label-free.



**Cardiac** - Record the four key measures of functional cardiac performance: action potential, field potential, propagation, and contractility.



**MEA Viability** - Measure cell viability and coverage on MEA plates for a complete structure-function assay.



**MEA Automation** - Automate Cardiac and Neural MEA assays with this API for interfacing with liquid handling platforms.



**Impedance** - Track cell proliferation, viability, barrier function, immune cell-mediated killing, viral cytopathic effects, and more.



**GxP Impedance** - Achieve FDA 21 CFR Part 11 compliance in GMP/GLP labs with this version of the Impedance Software Module.



**Impedance Automation** - Automate impedance assays with this API for interfacing with liquid handling platforms.

# MULTIWELL PLATES

## In a range of formats

All of Axion's MEA and impedance assay multiwell plates can be used with the Maestro Edge system. Select the multiwell plates to **match your assay needs**:

Plate Technology	Assay requirements							
	Field Potential	Action Potential (LEAP)	Contractility	Propagation	MEA viability	Electrical Stimulation	Optical Stimulation	Impedance
<b>Biocircuit MEA</b> For lowest cost per well MEA assays	●	●		●		●		
<b>Cytoview MEA</b> For MEA & cell imaging assays	●	●	●	●	●	●		
<b>Cytoview-Z</b> For impedance assays								●
<b>Lumos MEA</b> For optical stimulation MEA assays	●	●	●	●	●	●	●	



**BioCircuit MEA** - Maestro MEA plates with an opaque well bottom delivering high-quality results at the lowest cost per well. Available in 24-well format.



**CytoView MEA** - The premium Maestro MEA plate with a transparent well bottom for cell visualization and assay multiplexing. Available in 24-, and 6-well formats.



**Lumos MEA** - Maestro MEA plates designed for use with the Lumos system, featuring a transparent well bottom and light-focusing lid. Available in 24-well plate format.



**CytoView-Z** - The Maestro impedance plate with a transparent well bottom for cell visualization and assay multiplexing. Available in 96-well plate format.

**Learn more:**

*[axionbiosystems.com/maestro-edge](https://axionbiosystems.com/maestro-edge)*

**Contact us:**

*[axionbiosystems.com/contact](https://axionbiosystems.com/contact)*

**Office locations:**

North America - Europe - Asia Pacific

**For pricing and ordering:**

*[sales@axionbio.com](mailto:sales@axionbio.com)* or scan the QR code



For research use only. Not for use in diagnostic procedures.  
This information is subject to change without notice. © Axion BioSystems, Inc, 2022.