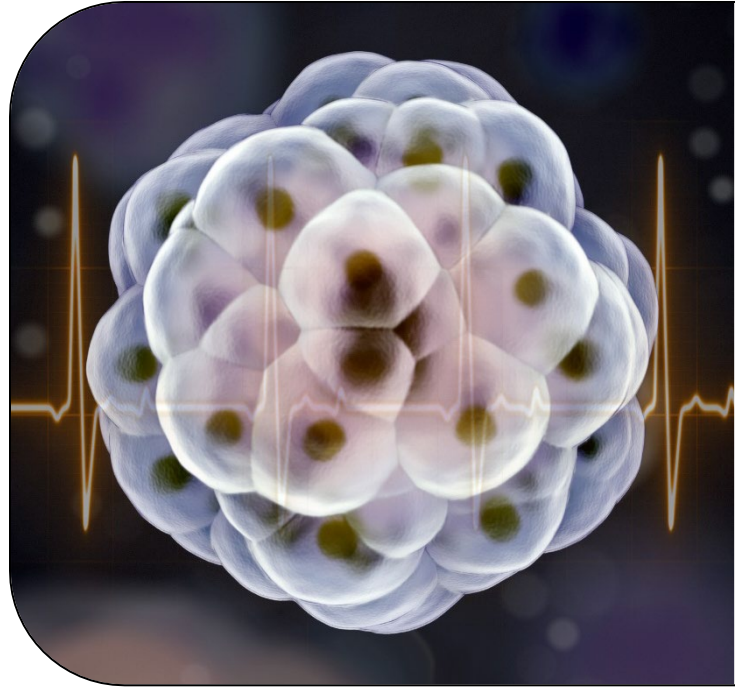




### >> Cardiac Organoids on Maestro MEA

Accurately recapitulating human cardiac activity in the lab is essential for **disease research, therapeutic discovery, and safety testing**—but traditional animal models often fall short. However, new approach methodologies like stem-cell derived cardiac organoids offer a more human-relevant platform for investigating pathological mechanisms and accelerating preclinical testing. Pairing 3D cardiac organoids with sensitive functional analysis provided by next-generation **Maestro MEA** offers high-fidelity replication of disease features and dependable assessments of drug responses *in vitro*.

Learn how Maestro MEA can support your cardiac organoid research with **these selected publications**:



#### Generation of human iPSCs derived heart organoids structurally and functionally similar to heart

Lee S-G, Kim Y-J, et al. *Biomaterials*. (2022)

Increased demand for physiologically relevant *in vitro* models has driven the development of organoids and other complex models to study diseases and screen therapeutics. In this study, researchers present and validate a self-organized heart organoid model from human induced pluripotent stem cells (hiPSCs).

#### Highlights:

- hiPSC-derived heart organoids display physiologically relevant structures, such as atrium- and ventricle-like areas.
- Electrophysiological analysis of hiPSC-derived heart organoids using Maestro MEA indicate functional maturation throughout organoid development.
- Organoids respond as expected to pharmacological treatment with common ion channel blockers, suggesting utility for cardiotoxicology and drug discovery applications.

[Read more >>](#)



## Modeling acute myocardial infarction and cardiac fibrosis using human induced pluripotent stem cell-derived multi-cellular heart organoids

Song M, Choi DB, et al. *Cell Death & Disease*. (2024)

The complexity of the adult heart makes it challenging to recapitulate cardiac diseases and injuries, such as myocardial infarction. Here, the authors present a novel human *in vitro* heart model using self-organizing heart organoids (HOs) containing multiple cardiac cell types. These HOs successfully mimic key pathological features of acute myocardial infarction and ischemia-reperfusion injury, indicating a promising alternative to animal studies for investigating heart diseases and for drug screening.

### Highlights:

- hiPSC-derived heart organoids contain multiple cardiac cell types, including cardiomyocytes, fibroblasts, and endothelial cells.
- An ischemia-reperfusion model using cobalt chloride provides stable hypoxia induction allowing for more accurate modeling of acute myocardial infarction and reperfusion.
- Organoids exposed to ischemia-reperfusion conditions exhibit tachycardia, prolonged field potential duration, and decreased spike amplitude and conduction velocity, indicative of significant functional deficits detected by the Maestro.

[Read more >>](#)

## Evaluation of the cardiotoxicity of Echinchrome A using human induced pluripotent stem cell-derived cardiac organoids

Lee S-J, Kim E, et al. *Ecotoxicology and Environmental Safety*. (2025)

Despite its therapeutic potential for cardiovascular disease, the cardiac safety of Echinchrome A (EchA) requires further investigation. To address this gap, the authors employed human iPSC-derived cardiac organoids (hCOs) to assess whether EchA causes heart damage or irregular heart rhythms.

### Highlights:

- hiPSC-derived cardiac organoids express key ion channels and markers indicative of functional maturity.
- The Maestro MEA platform was used to evaluate hCO function as an indicator of cardiotoxicity, finding low cardiotoxic potential with little change in beat period and field potential duration.



- In addition to supporting the safety and therapeutic use of EchA, this study highlights the translational relevance of cardiac organoids for toxicity assessment.

[Read more >>](#)

### Frameshift variants in C10orf71 cause dilated cardiomyopathy in human, mouse, and organoid models

Li Y, Ma K, et al. *The Journal of Clinical Investigation*. (2024)

Dilated cardiomyopathy (DCM) has widely been associated with various genetic mutations. Researchers identified a candidate causal gene, C10orf71, and used hiPSC-derived cardiomyocytes and organoids to assess the effects of C10orf71 mutation on cardiomyocyte electrophysiology and DCM.

#### Highlights:

- C10orf71 frameshift mutations are associated with DCM, and C10orf71 is specifically expressed in cardiomyocytes.
- Contractility analysis of hiPSC-derived heart organoids using Maestro MEA indicates impaired contractile function – evident via decreased amplitude and increased excitation-contraction delay.
- Contractile protein activation via OM treatment rescues contractile dysfunction *in vivo*, suggesting therapeutic potential for DCM treatment.

[Read more >>](#)

*See how easy it is to get functional data for your next cardiac organoid publication.*

**Find out more at:**  
[axionbiosystems.com/maestro-pro](https://axionbiosystems.com/maestro-pro)

