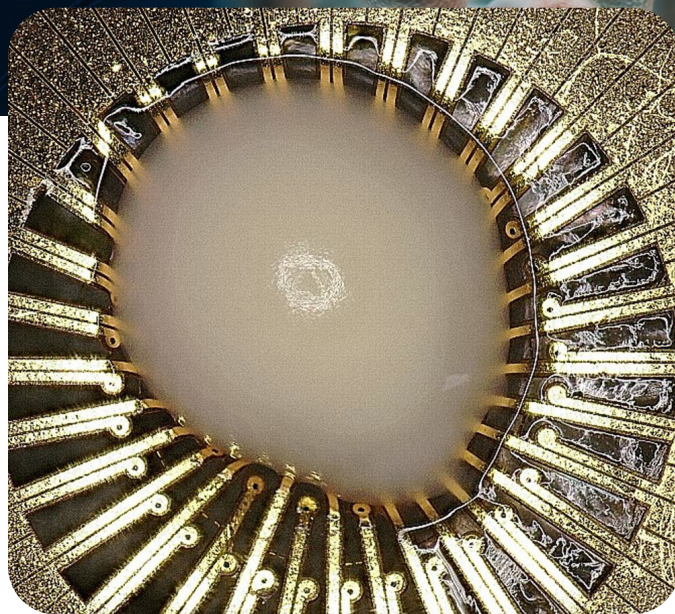




# >> Organoids on Maestro MEA

*The MEA for organoid research*



## The future is 3D

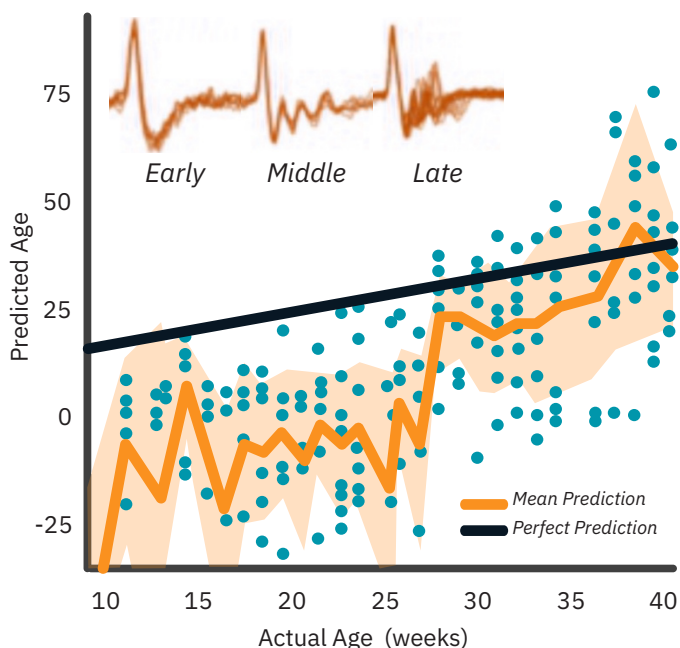
Organoids are transforming brain research, providing new insights into human development, disease modeling, and therapeutic discovery.

Maestro MEA is the **platform built for organoids**, delivering long-term, noninvasive recordings that unlock the full potential of these complex 3D models.

## The MEA of choice

With over 85 peer-reviewed publications, **researchers worldwide choose Maestro MEA** to push the boundaries of neuroscience and accelerate discovery.

 **85+**  
Organoid Publications



## Complete functional readouts

Track how neural organoids grow, mature, and model disease with Maestro MEA's rich, noninvasive electrophysiological recordings.

- **Analyze** over two dozen neural metrics
- **Monitor** the emergence of complex LFP activity
- **Test** multiple conditions simultaneously

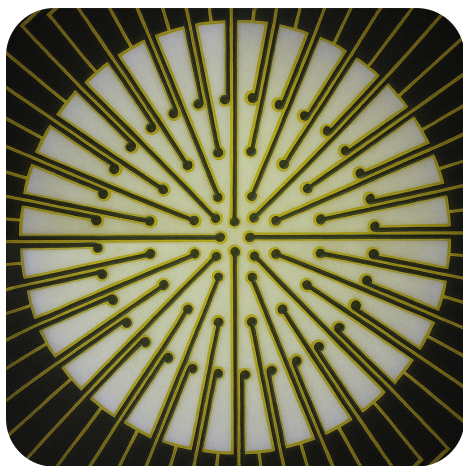
Figure: As cerebral organoids matured, their MEA activity patterns closely mirrored *in vivo* EEG recordings - highlighting the relevance of organoid data to human brain development. Data from: Muotri, Alysson. (2019, Sep. 17). Measuring oscillatory waves in cerebral organoids [Webinar]. Axion BioSystems.

# >> Designed for organoids

To learn more:  
Scan the QR  
code or visit:  
[axionbio.com](http://axionbio.com)

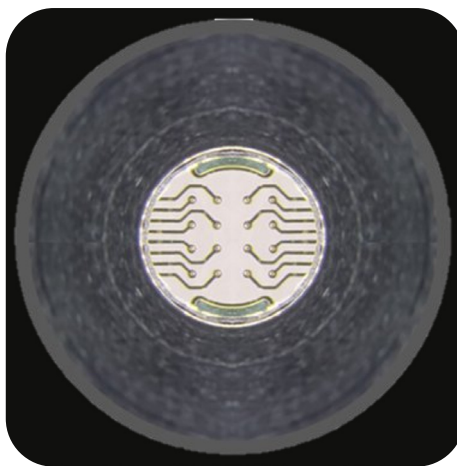


Studying 3D cultures like organoids and spheroids requires specialized tools. Axion BioSystems' **organoid MEA** portfolio is purpose-built to address the unique needs of these complex models.



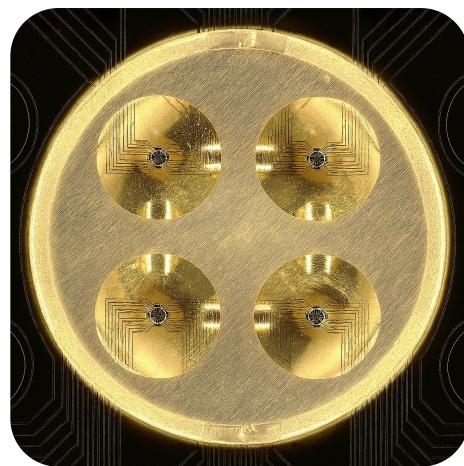
**3DMap™**

Flexible electrodes for tracking activity across the surface



**SpheroGuide™**

High-throughput, targeted placement for large organoids

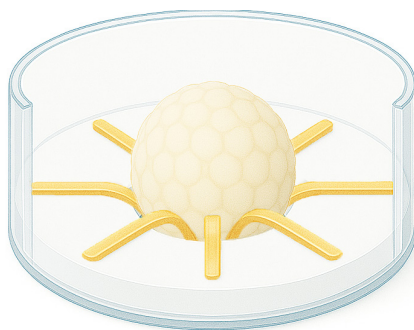


**SpheroHD™**

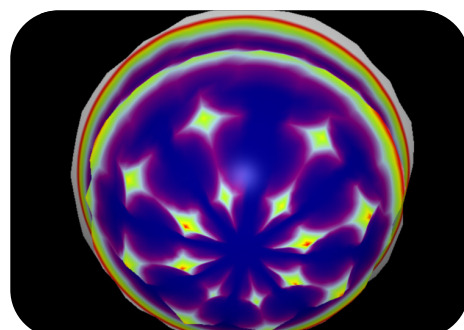
High-density electrodes for small organoids

## Fit to your **biology**

**3DMap** is designed to interface with 3D organoids like never before, using **flexible electrode technology** to capture signals across the surface.



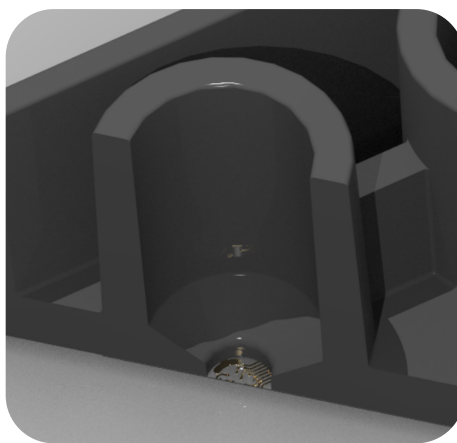
Flexible electrodes wrap around the organoid



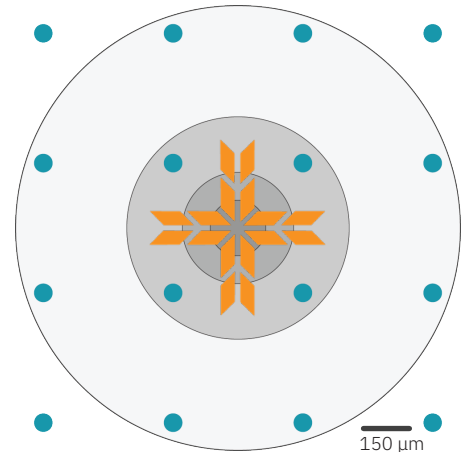
Activity is mapped in 3D in AxIS software

## No matter the **size**

**SpheroGuide** and **SpheroHD** have you covered. Both feature **integrated placement funnels** for fast, accurate positioning. SpheroGuide is ideal for larger organoids, while SpheroHD's **high-density electrode** arrays are perfect for smaller organoids or spheroids.



Well shape guides organoids to the electrodes



**SpheroGuide** and **SpheroHD** electrode layout