

>> AxIS Analysis

From spikes to insights — instantly

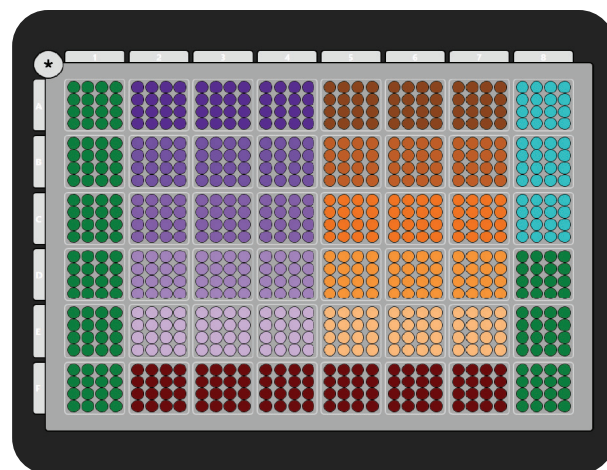
AxIS Analysis transforms neural data into actionable insights. Only on Maestro MEA, the software streamlines complex datasets into clear, publication-ready results.



Organize

Gather all your recordings into one place and export results to the Lab Notebook.

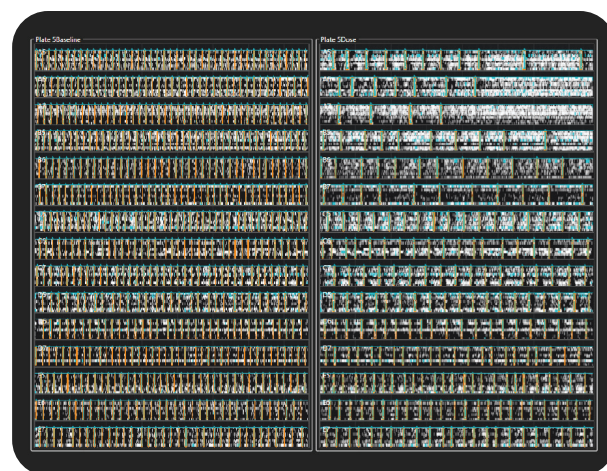
- Import multiple recordings
- Edit platemaps
- Export to single Excel file



Visualize

See your data like never before. Easily optimize detection and compare conditions and time points.

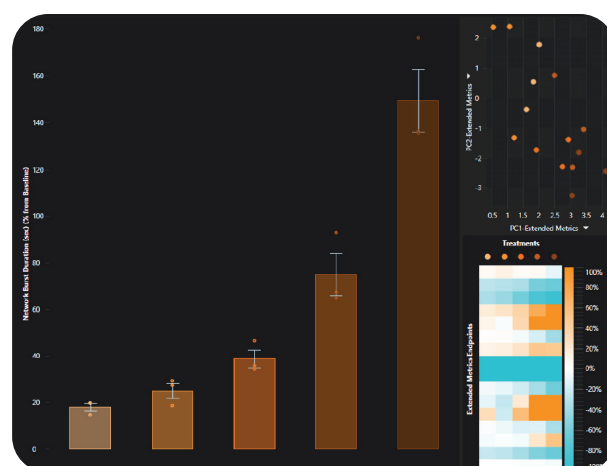
- Side-by-side rasters
- Instant result plotting
- Visual detection verification



Analyze

Gain access to all new multiparameter analysis, apply controls, and export figures with just a few clicks.

- Interactive heat map
- PCA plots
- Customizable metric sets



Learn more:

axionbiosystems.com/products/axis-analysis

Contact us:

axionbiosystems.com/contact

Office locations:

North America - Europe - Asia Pacific

For pricing and ordering:

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, 2025.