

BIOCIRCUIT MEA 24 DATASHEET

M384-BIO-24



Each well contains 16 electrodes arranged in a 4 x 4 grid.

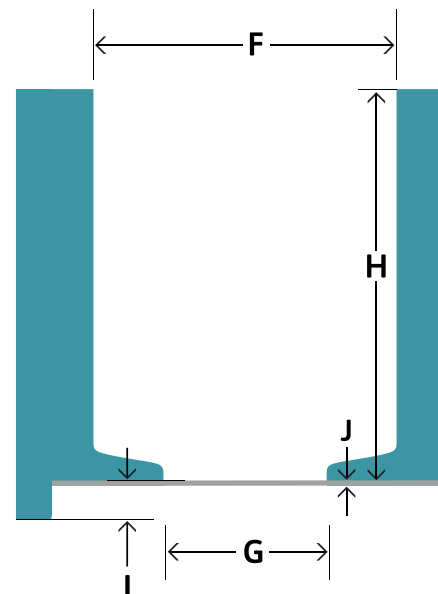
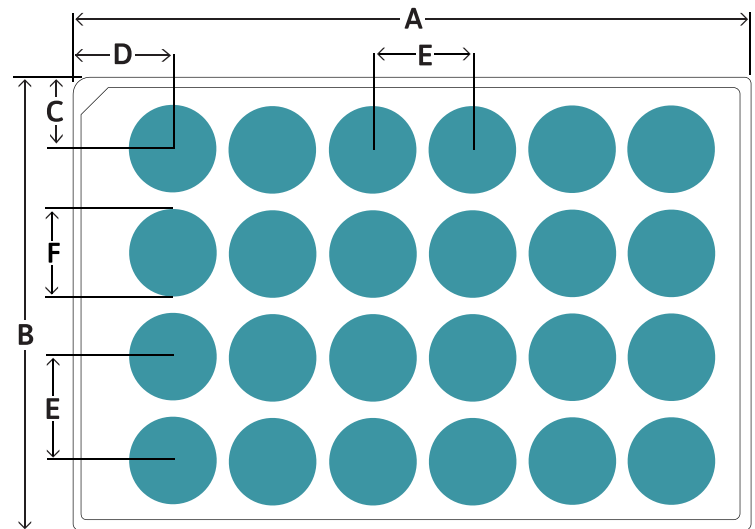
ELECTRODE SPACING: 350 μm
ELECTRODE DIAMETER: 50 μm
RECORDING AREA: 1.1 mm x 1.1 mm

The BioCircuit MEA 24 delivers high-quality MEA results and additional stimulation capacity with a dedicated stimulation electrode (enlarged blue circle). The AccuSpot feature (shown in gray) acts as an on-plate spotting guide to confine a 5 μl droplet of plated cells to an area of 7.65 mm^2 over the recording electrode region, to improve plate preparation ease and simplicity.

PLATE INFORMATION

Electrode Material: Gold

PLATE DIMENSIONS	
Plate Length (A)	127.76 mm
Plate Width (B)	85.48 mm
A1 Offset (Row) (C)	15.69 mm
A1 Offset (Column) (D)	18.83 mm
Well-to-Well Distance (E)	18.00 mm
Well Diameter (Top) (F)	15.00 mm
Well Diameter (Bottom) (G)	5.4 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.33 mm
Well Volume	2,000 μl



STORING AND HANDLING

- Store in clean, dry environment at room temperature
- Do not expose to ultraviolet light
- Do not autoclave
- Use clean gloves when handling plates
- Designed for one-time cell plating
- 6-month warranty from date of receipt

SOFTWARE REQUIREMENTS

BioCircuit MEA 24 plates require AxIS Navigator software version 1.5 (or later) for the Maestro Pro and Edge systems. For the first-generation Maestro system, BioCircuit MEA 24 plates require AxIS software version 2.5 (or later).

Please contact us if you have any questions or need assistance with a software upgrade.

