

# BIOCIRCUIT MEA 96 DATASHEET

M768-BIO-96



Each well contains 8 electrodes.

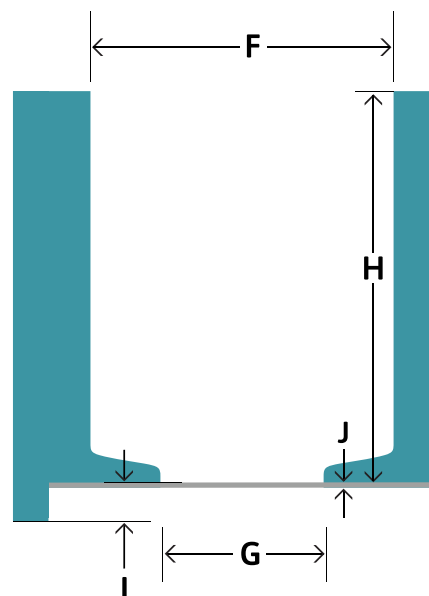
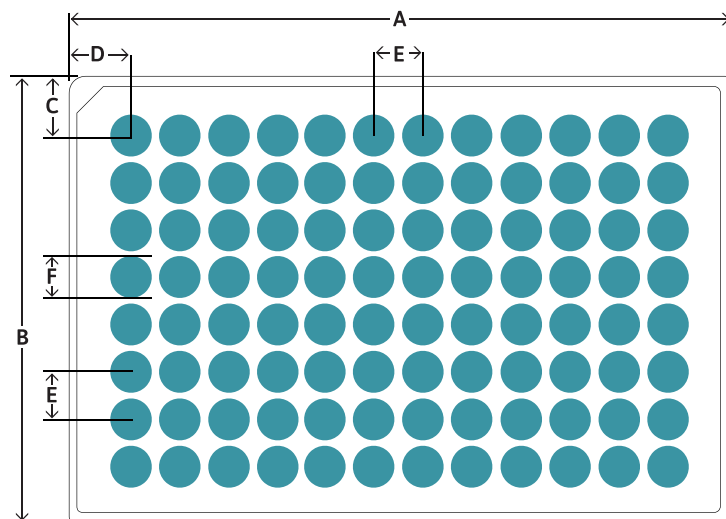
ELECTRODE SPACING: 350  $\mu\text{m}$   
ELECTRODE DIAMETER: 50  $\mu\text{m}$   
RECORDING AREA: 0.8 mm x 0.8 mm

The BioCircuit MEA 96 delivers high-quality MEA results and the benefit of plate preparation ease and simplicity. The AccuSpot feature (shown in gray) acts as an on-plate spotting guide to confine a 5  $\mu\text{l}$  droplet of plated cells to an area of 7.65 mm<sup>2</sup> over the recording electrode region.

## PLATE INFORMATION

Electrode Material: Gold

PLATE DIMENSIONS	
Plate Length (A)	127.76 mm
Plate Width (B)	85.48 mm
A1 Offset (Row) (C)	11.24 mm
A1 Offset (Column) (D)	14.38 mm
Well-to-Well Distance (E)	9.00 mm
Well Diameter (Top) (F)	8.07 mm
Well Diameter (Bottom) (G)	6.40 mm
Well Height (H)	12.75 mm
Well Bottom Elevation (I)	1.60 mm
Base Thickness (J)	0.53 mm
Well Volume	500 $\mu\text{l}$



## 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 96 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 96 plates require AxIS software version 2.5 (or later).

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

