

CYTOVIEW MEA 96 DATASHEET

M768-tMEA-96B **BLACK** | M768-tMEA-96W **WHITE**



Each well contains 8 electrodes.

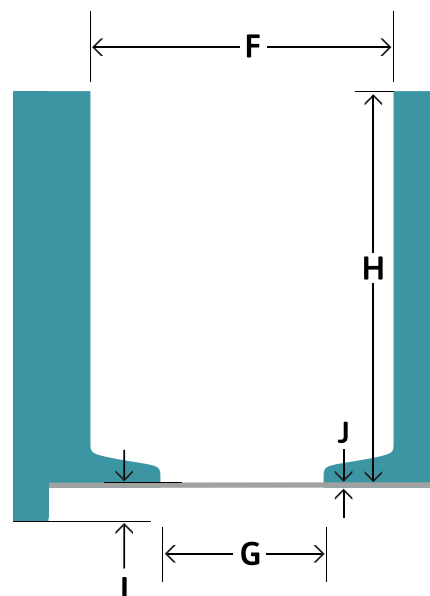
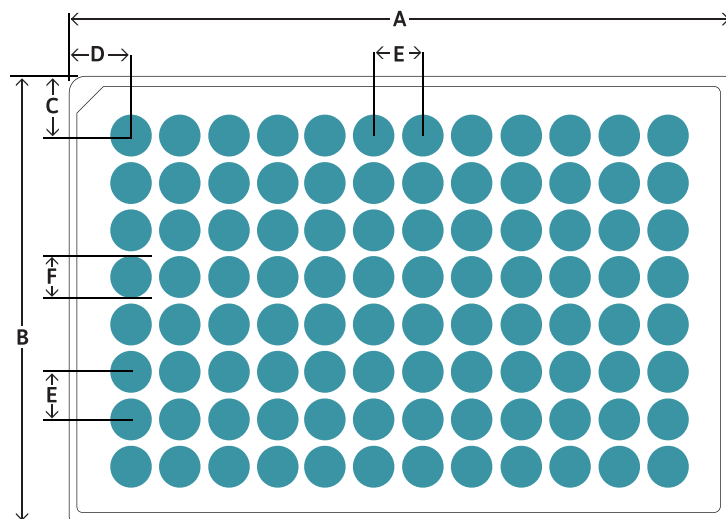
ELECTRODE SPACING: 350 μm
 ELECTRODE DIAMETER: 50 μm
 RECORDING AREA: 0.8 mm x 0.8 mm

The CytoView MEA 96 delivers high-quality MEA results with a transparent well bottom for cell visualization and assay multiplexing. The on-plate spotting guides (shown in gray) aid plate preparation.

PLATE INFORMATION

Electrode material: PEDOT; Insulation/Surface material: SU-8

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)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.15 mm
Well Volume	500 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
- 12-month warranty from date of receipt

SOFTWARE REQUIREMENTS

CytoView MEA 96 plates require AxIS Navigator software version 1.5 (or later) for the Maestro Pro system. For the first-generation Maestro system, CytoView 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.

