

CYTOVIEW MEA 48 DATASHEET

M768-tMEA-48B BLACK | M768-tMEA-48W WHITE



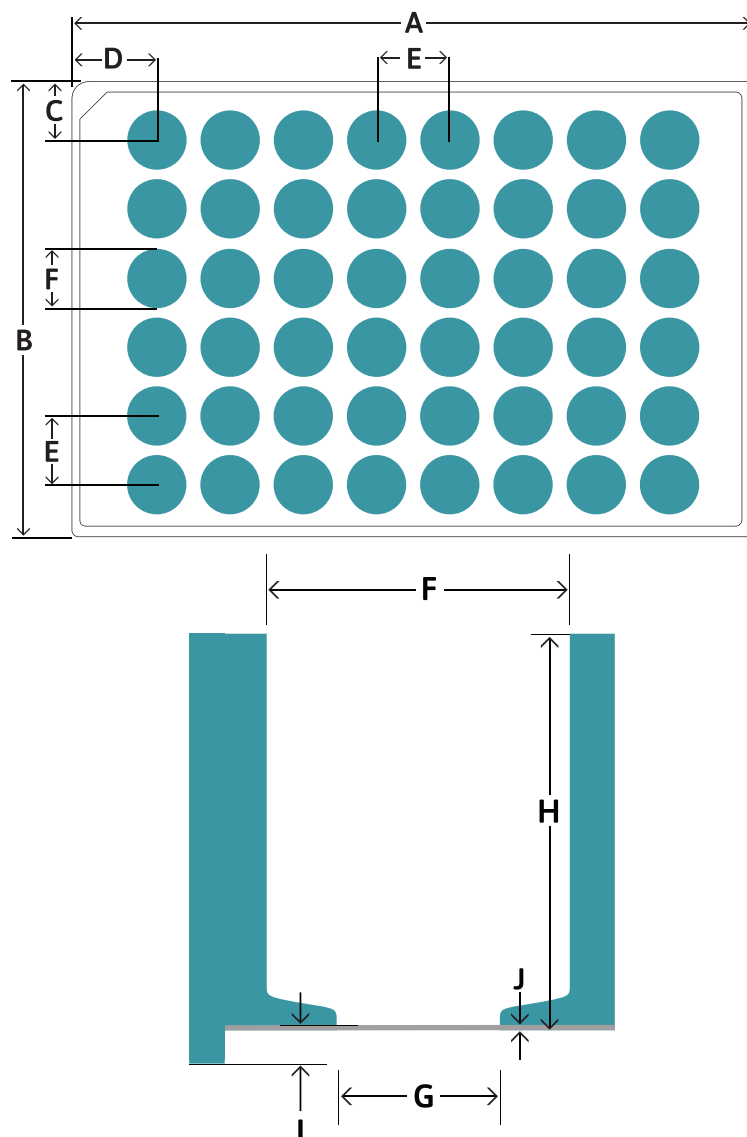
Each well contains 16 electrodes arranged in a 4 x 4 grid.
ELECTRODE SPACING 350 μm
ELECTRODE DIAMETER 50 μm
RECORDING AREA 1.1 mm x 1.1 mm

The CytoView MEA 48 provides confidence and versatility to 48-well MEA assays with a transparent bottom for cell visualization and assay multiplexing.

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)	10.04 mm
A1 Offset (Column) (D)	18.10 mm
Well-to-Well Distance (E)	13.08 mm
Well Diameter (Top) (F)	10.35 mm
Well Diameter (Bottom) (G)	5.30 mm
Well Height (H)	12.15 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	150 μm
Well Volume	950 μl



STORING AND HANDLING

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

SOFTWARE REQUIREMENTS

CytoView MEA plates require AxIS software version 2.4 (or later). When recording from CytoView plates, choose "CytoView MEA 48" from the drop-down plate selector, and remember to change your selection back to "Classic MEA 48" for traditional/opaque plates.
Please contact us if you have any questions or need assistance with a software upgrade.

