

CYTOVIEW MEA 12 DATASHEET

M768-tMEA-12W WHITE



Each well contains 64 electrodes arranged in an 8 x 8 grid.

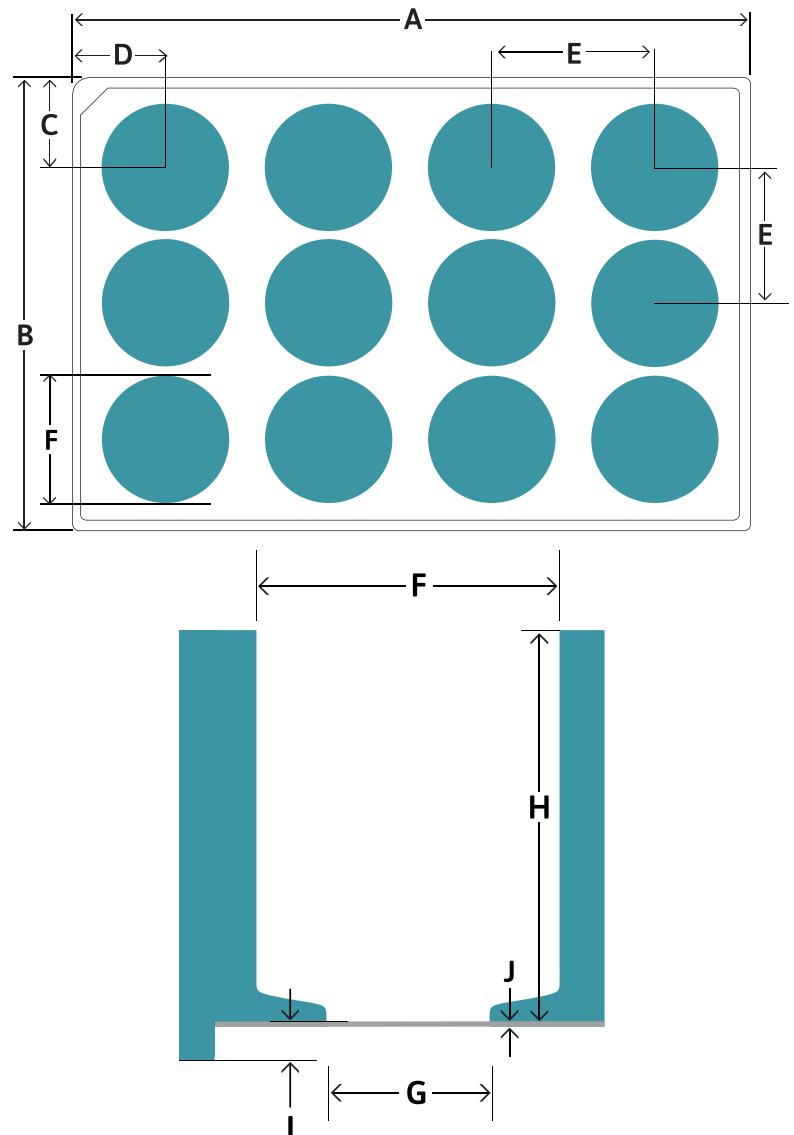
ELECTRODE SPACING: 300 μm
ELECTRODE DIAMETER: 50 μm
RECORDING AREA: 2.1 mm x 2.1 mm

The CytoView MEA 12 delivers high-quality MEA results with a transparent well 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)	16.69 mm
A1 Offset (Column) (D)	24.83 mm
Well-to-Well Distance (E)	26.00 mm
Well Diameter (Top) (F)	22.00 mm
Well Diameter (Bottom) (G)	11.5 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.15 mm
Well Volume	3,000 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
- 12-month warranty from date of receipt

SOFTWARE REQUIREMENTS

CytoView MEA 12 plates require AxIS Navigator software version 3.12.4 (or later) for the Maestro Pro system.

Note: These plates are only compatible with the Maestro Pro. CytoView MEA 12 plates sold prior to 2021 are now designated as CytoView MEA 12 (Legacy) in AxIS Navigator and have a different layout.

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

