

LUMOS MEA 48 DATASHEET

M768-tMEA-48OPT



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

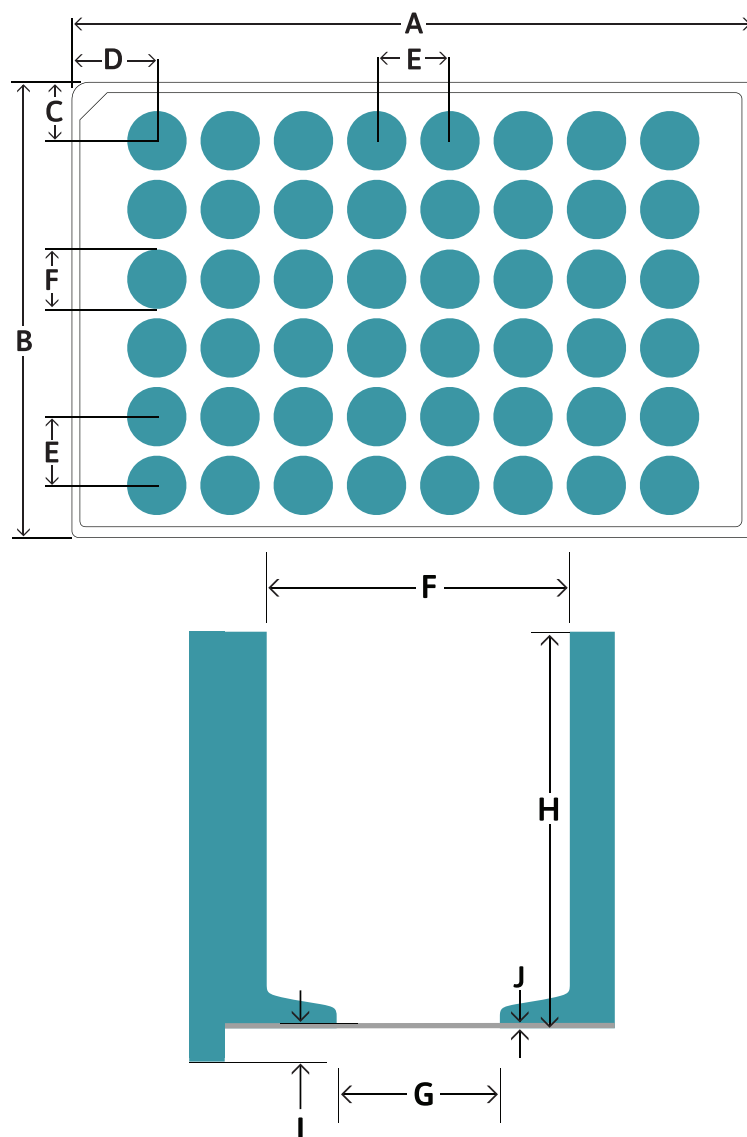
ELECTRODE SPACING 350 μm
ELECTRODE DIAMETER 50 μm
RECORDING AREA 1.1 mm x 1.1 mm

The Lumos MEA 48 combines the robustness and assay flexibility of a CytoView MEA plate with white walls and a custom lensed lid for optimal light delivery in each well.

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

Lumos MEA plates require AxIS software version 2.4 (or later). When recording from Lumos plates, choose "Lumos 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.

