

LUMOS MEA 48 DATASHEET

M768-tMEA-48OPT



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

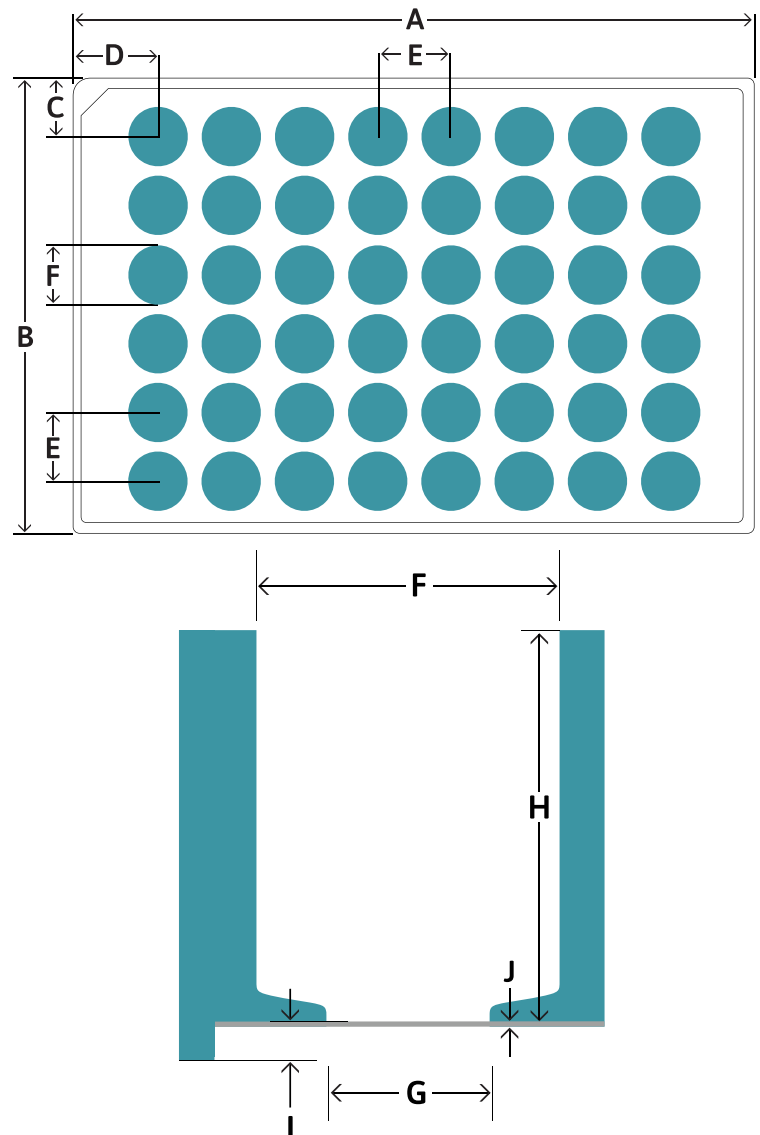
ELECTRODE SPACING: 350 μm
ELECTRODE DIAMETER: 50 μm
RECORDING AREA: 1.1 mm x 1.1 mm

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

Surface 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
- Do not autoclave
- Use clean gloves when handling plates
- Designed for one-time cell plating
- 12-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.

