

LUMOS MEA 24 DATASHEET

M384-tMEA-24OPT



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

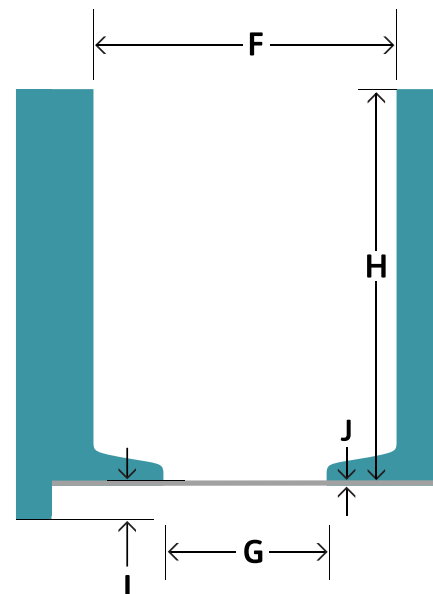
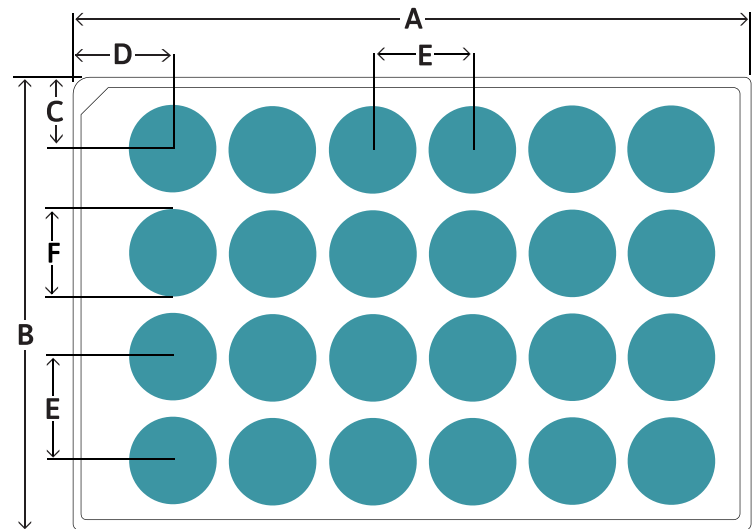
ELECTRODE SPACING: 350 μm
ELECTRODE DIAMETER: 50 μm
RECORDING AREA: 1.1 mm x 1.1 mm

The Lumos MEA 24 combines the robustness and assay flexibility of a CytoView MEA plate with white walls and a custom Lumos lid for optimal light delivery in each well. The dedicated stimulation electrode (300 μm diameter, enlarged blue circle) provides superior stimulation capacity, and 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)	15.69 mm
A1 Offset (Column) (D)	18.83 mm
Well-to-Well Distance (E)	18.00 mm
Well Diameter (Top) (F)	15.00 mm
Well Diameter (Bottom) (G)	5.4 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.15 mm
Well Volume	2,000 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

Lumos MEA 24 plates require AxIS Navigator software version 1.5 (or later) for the Maestro Pro and Edge systems.

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

