

# LUMOS MEA 96 DATASHEET

M768-tMEA-96OPT



Each well contains 8 electrodes.

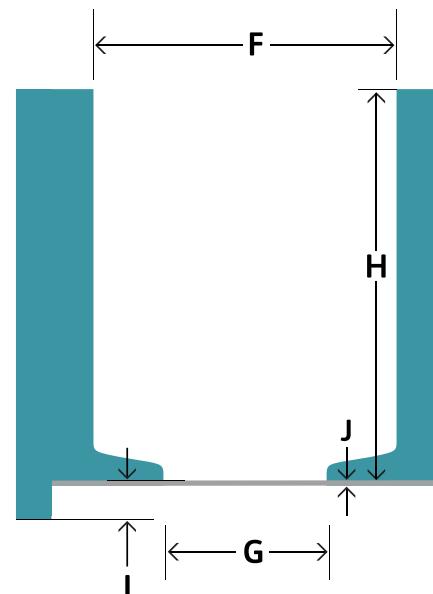
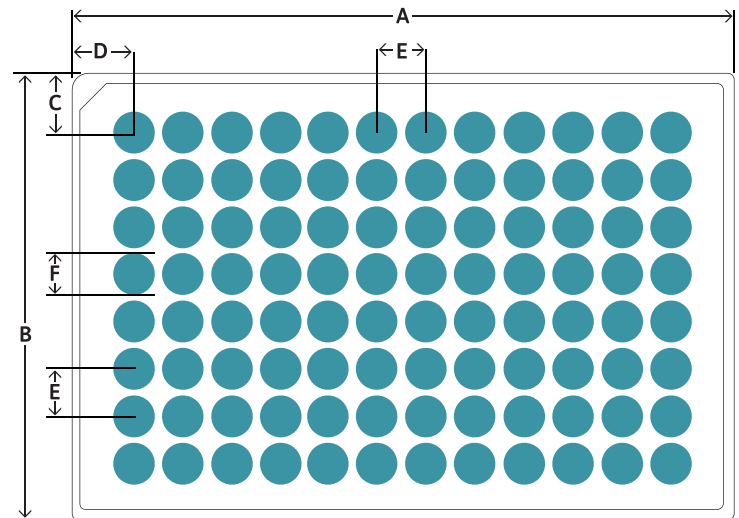
ELECTRODE SPACING: 350  $\mu\text{m}$   
ELECTRODE DIAMETER: 50  $\mu\text{m}$   
RECORDING AREA: 0.8 mm x 0.8 mm

The Lumos MEA 96 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. On-plate spotting guides (shown in gray) aid plate preparation.

## 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)	11.24 mm
A1 Offset (Column) (D)	14.38 mm
Well-to-Well Distance (E)	9.00 mm
Well Diameter (Top) (F)	8.07 mm
Well Diameter (Bottom) (G)	6.4 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.15 mm
Well Volume	500 $\mu\text{l}$



## 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
- 6-month warranty from date of receipt

## SOFTWARE REQUIREMENTS

Lumos MEA 96 plates require AxIS Navigator software version 1.5 (or later) for the Maestro Pro system.

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

