

# SPHEROGUIDE MEA 48 DATASHEET

M768-SPH-48B **BLACK**



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

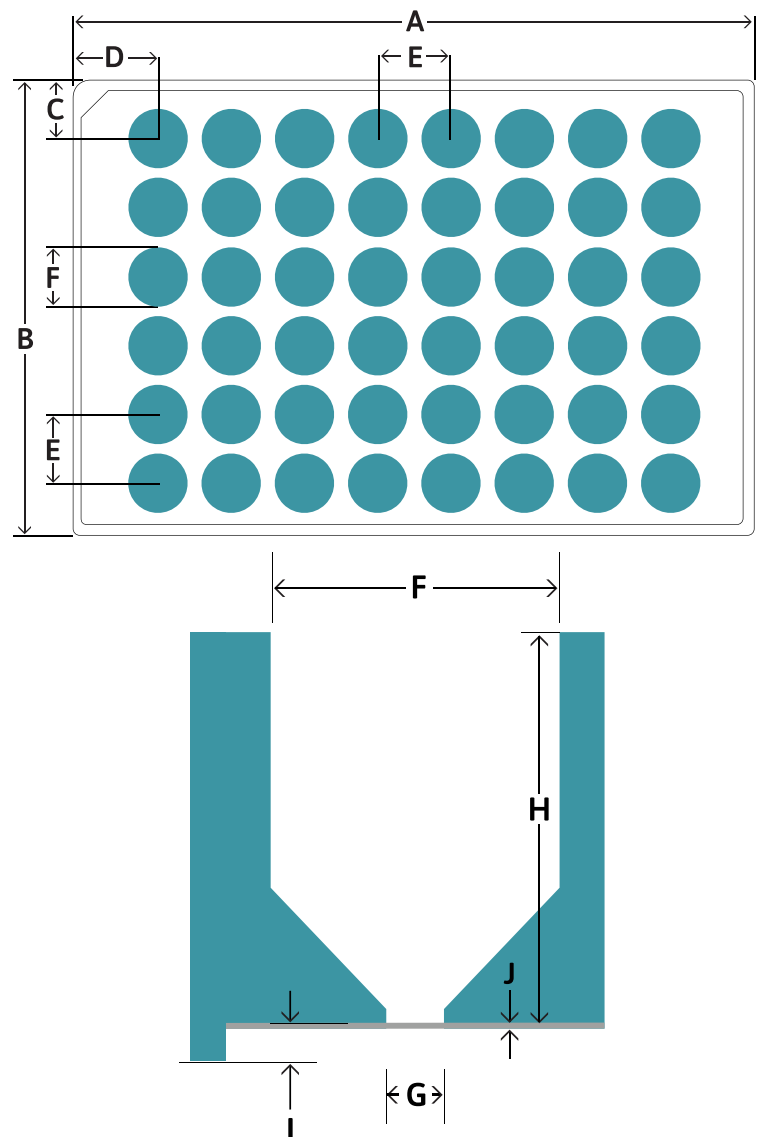
ELECTRODE SPACING: 350  $\mu\text{m}$   
ELECTRODE DIAMETER: 50  $\mu\text{m}$   
RECORDING AREA: 1.1 mm x 1.1 mm

The Spheroguide MEA 48 provides consistency and accuracy to spheroid and organoid MEA assays with an integrated placement funnel for reliable spheroid positioning and transparent bottom for spheroid visualization and multiplexing

## 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)	7.0 mm
Well Diameter (Bottom) (G)	2.5 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	150 $\mu\text{m}$
Well Volume	435 $\mu\text{l}$
Working Volume	300 $\mu\text{l}$



## 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
- 12-month warranty from date of receipt

## SOFTWARE REQUIREMENTS

Spheroguide plates require AxIS Navigator version 3.11 (or later). When recording from Spheroguide plates, choose "Spheroguide MEA 48" from the drop-down plate selector, if not auto-selected upon plate docking.

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

