

# CYTOVIEW MEA 96 DATASHEET

M768 - tMEA - 96B **BLACK** | M768 - tMEA - 96W **WHITE**



Each well contains 8 electrodes.

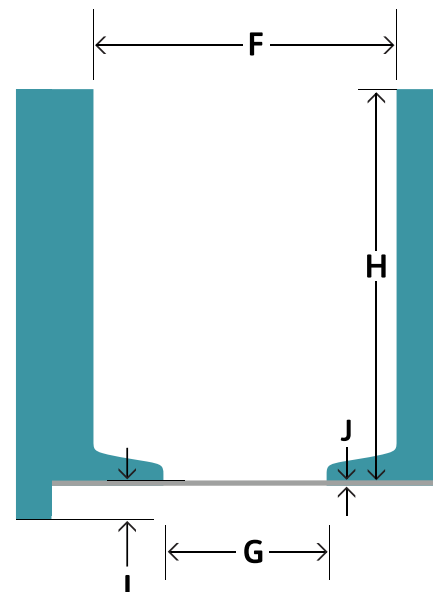
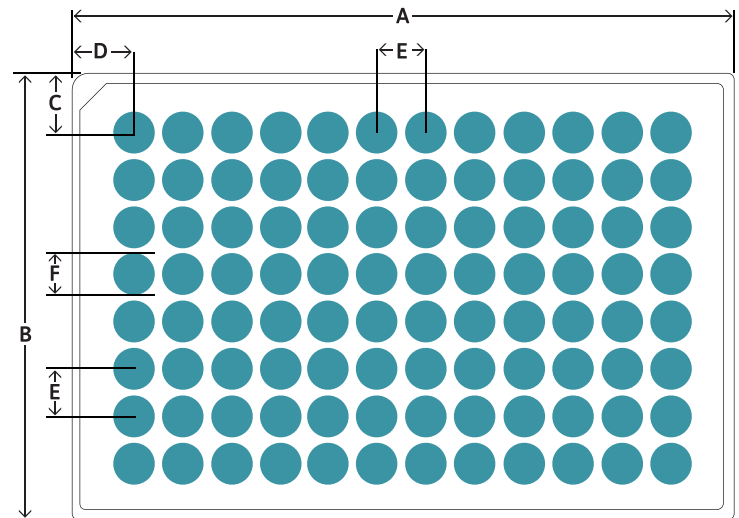
**ELECTRODE SPACING:** 350  $\mu\text{m}$   
**ELECTRODE DIAMETER:** 50  $\mu\text{m}$   
**RECORDING AREA:** 0.8 mm x 0.8 mm

The CytoView MEA 96 delivers high-quality MEA results with a transparent well bottom for cell visualization and assay multiplexing. The 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)        | 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) | 6.40 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

CytoView MEA 96 plates require AxIS Navigator software version 1.5 (or later) for the Maestro Pro system. For the first-generation Maestro system, CytoView MEA 96 plates require AxIS software version 2.5 (or later).

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

