

# CYTOVIEW MEA 24 DATASHEET

M384-tMEA-24W WHITE



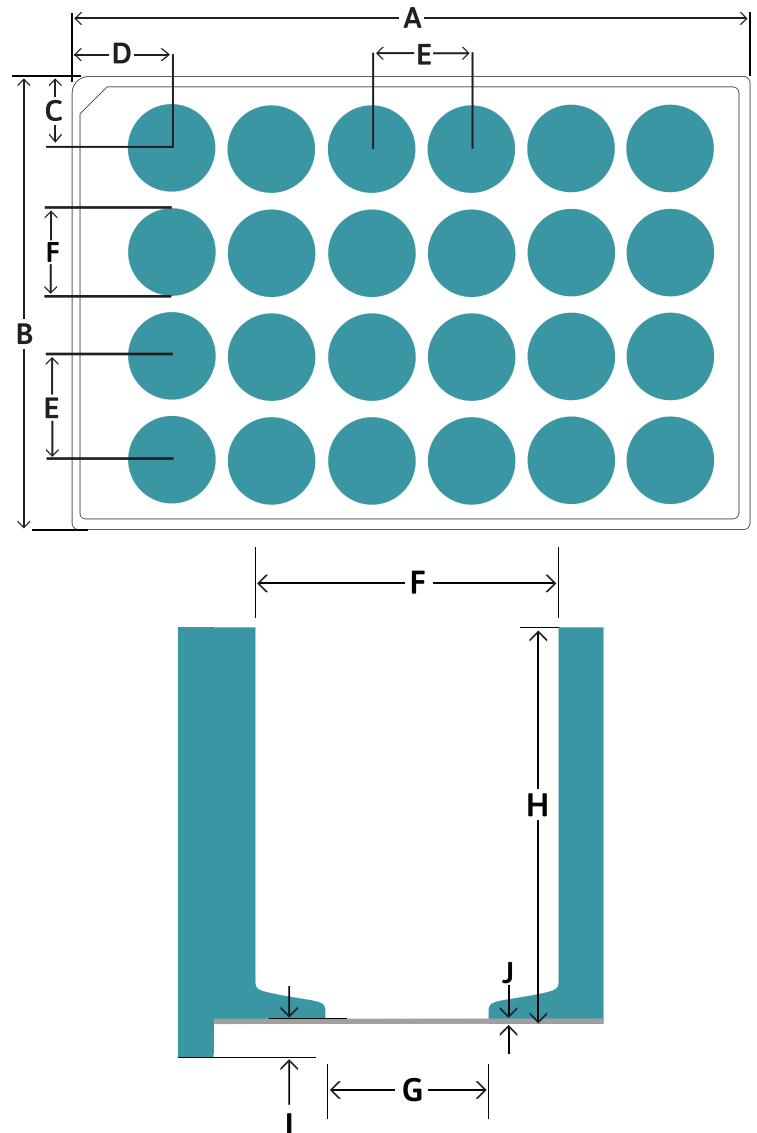
Each well contains 16 electrodes arranged in a 4 x 4 grid.  
**ELECTRODE SPACING** 350  $\mu\text{m}$   
**ELECTRODE DIAMETER** 50  $\mu\text{m}$   
**RECORDING AREA** 1.1 mm x 1.1 mm

The CytoView MEA 24 delivers high-quality MEA results with a transparent well bottom for cell visualization. The dedicated stimulation electrode (300  $\mu\text{m}$  diameter, enlarged blue circle) provides superior stimulation capacity, and on-plate spotting guides (shown in gray) aid plate preparation.

## PLATE INFORMATION

Microelectrode 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)	150 $\mu\text{m}$
Well Volume	2,000 $\mu\text{L}$



## 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
- 12-month warranty from date of receipt

