

CYTOVIEW-Z 96 DATASHEET

Z96-IMP-96B **BLACK**



Each well contains 1 interdigitated electrode.

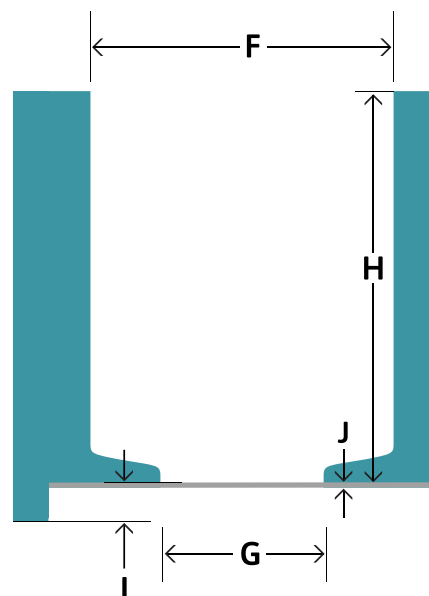
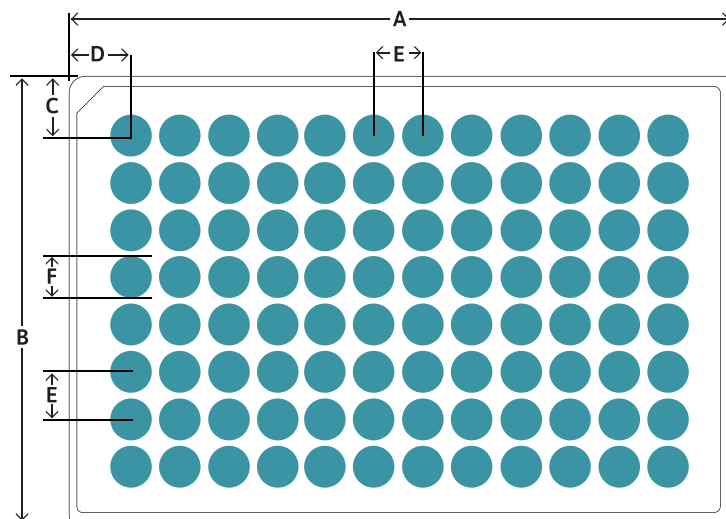
ELECTRODE SPACING: 100 μm
ELECTRODE WIDTH: 100 μm
RECORDING AREA: 20 mm^2

The CytoView-Z 96 delivers high-quality impedance results with a transparent well bottom for cell visualization and assay multiplexing. The on-plate viewing windows (shown in gray, 500 μm x 3000 μm) aid cell visualization and imaging.

PLATE INFORMATION

Electrode material: Gold; Surface material: PET

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)	6.4 mm
Well Diameter (Bottom) (G)	5.0 mm
Well Height (H)	13.2 mm
Well Bottom Elevation (I)	1.15 mm
Base Thickness (J)	0.15 mm
Well Volume	396 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
- 6-month warranty from date of receipt

SOFTWARE REQUIREMENTS

CytoView-Z 96 plates require AxIS Z software version 1.3 (or later).

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

