

# LUMOS OPTICLEAR 24 DATASHEET

OPT-24



Designed for optical stimulation with the Lumos system.

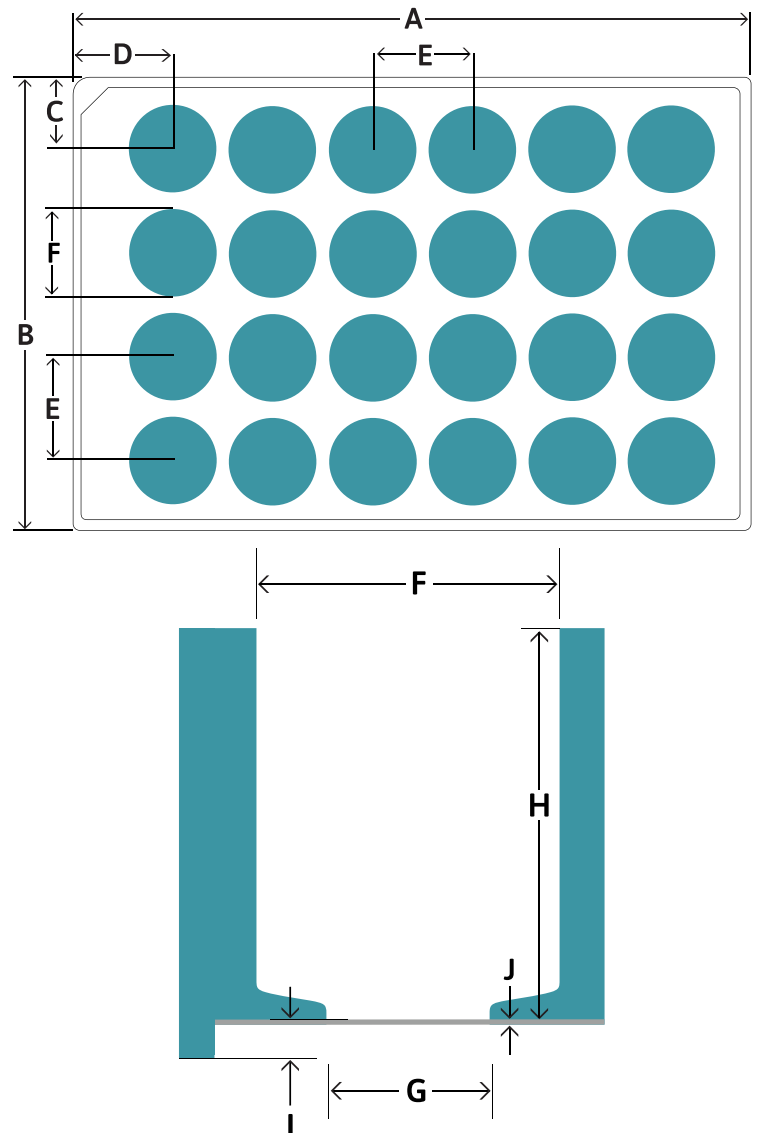
**WAVELENGTHS/WELL: 4**

The Lumos OptiClear 24 plate has a high clarity bottom culture surface for microscopy and high content imaging. The specially formulated reflective white walls and a custom Lumos lid facilitate optimal light delivery within each well, with minimum crosstalk between wells.

## PLATE INFORMATION

**Surface material:** Cyclic Olefin Copolymer (COC)

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)	0.15 mm
Well Volume	2,000 $\mu$ 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

For optical stimulation, use the Lumos 24 head and Lumos Studio software, or AxIS Navigator software version 1.5 (or later), or AxIS software version 2.5 (or later).

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

