

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of polypropylene StorPlate-96 round well, V-bottom, 450 μ L volume

Part numbers: PN 6008290 / 6008299 Natural polypropylene, untreated

Material and treatment: Round well, polypropylene 96-well microplate with V-shaped bottoms. Well volume 0.45 mL. Free of DNase, and RNase. Untreated, non-sterile.

Packaging: Part number 6008290 is a case with 50 plates, packed as 5 bags of 10 plates each. 6008299 is a case with 200 plates.

Dimensions:

Microplate Dimensions	6008290 / 6008299
Length at base plane	127.76 mm
Width at base plane	85.48 mm
Height overall	14.35 mm
Well depth	12.43 mm
Well diameter at opening	8.50 mm
Well diameter at bottom	NA
Well volume	450 μ l
Well spacing	9.0 mm
Skirt height	2.45 mm
Well offset:	
Left edge to well A1	14.3 mm
Top edge to well A1	11.2 mm