

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of polypropylene StorPlate-384 square well, V-bottom, 120 µL volume

Part numbers: PN 6008590 / 6008598 Natural polypropylene, untreated

Material and treatment: Square well, polypropylene 384-well microplate with V-shaped bottoms. Well volume 120µL. Free of DNase, and RNase. Untreated, non-sterile.

Packaging: Part number 6008590 is a case with 50 plates, packed as 5 bags of 10 plates each. 6008598 is a case with 200 plates.

Dimensions:

Microplate Dimensions	6008590 / 6008598
Length at base plane	127.76 mm
Width at base plane	85.48 mm
Height overall	14.35 mm
Well depth	11.85 mm
Well diameter at opening	3.50 mm
Well diameter at bottom	NA
Well volume	120 µl
Well spacing	4.5 mm
Skirt height	2.5 mm
Well offset:	
Left edge to well A1	12.13 mm
Top edge to well A1	8.99 mm