

Research use only. Not for use in diagnostic procedures.

Technical datasheet: Dimensions of 1/2Area Plates-96

Part numbers:	1/2Area OptiPlate-96	PN 6002290 / 6002299	White, untreated
	1/2Area OptiPlate-96 HB	PN 6002520	White, high protein binding
	1/2Area OptiPlate-96 F	PN 6002270 / 6002279	Black, untreated
	1/2Area AlphaPlate-96	PN 6002350 / 6002359	Light-grey, untreated
	1/2Area SpectraPlate-96	PN 6002640 / 6002649	Clear, medium protein binding

Material and treatment: Black, white, clear, or light-grey polystyrene microplate with 96 flat-bottom wells with reduced diameter / volume. Plate is available untreated, TC-treated or with high protein binding affinity.

Packaging: 6002290/6002520/6002270/ 6002350/6002640 is a case with 50 plates, packed as 2 bags of 25 plates. 6002299/6002279/6002359/6002649 is a box with 200 plates, packed as 8 bags of 25 each.

Dimensions -- Customer Drawing:

