

Research use only. Not for use in diagnostic procedures.

Technical datasheet:

Dimensions of OptiPlate-1536 / CulturPlate-1536 / SpectraPlate-1536 / AlphaPlate-1536

Part numbers:	OptiPlate-1536	PN 6004290	White, untreated
	OptiPlate-1536 HB	PN 6004620	White, high protein binding
	CulturPlate-1536	PN 6004680 / 6004684	White, TC-treated
	OptiPlate-1536 F	PN 6004270	Black, untreated
	OptiPlate-1536 F HB	PN 6004520	Black, high protein binding
	CulturPlate-1536 F	PN 6004660 / 6004664	Black, TC-treated
	SpectraPlate-1536	PN 6004640	Clear, medium protein binding
	SpectraPlate-1536 HB	PN 6004500	Clear, high protein binding
	SpectraPlate-1536 TC	PN 6004650 / 6004654	Clear, TC-treated
	AlphaPlate-1536	PN 6004350	Light-grey, untreated

Material and treatment: Black, white, transparent or light gray polystyrene microplate with 1536 flat-bottom wells. Plate is available untreated, with high protein binding affinity (HB) or Tissue-Culture (TC) treated.

Packaging:

Format	Type	Part numbers
Case of 50 plates	Individually pouched (with lids)	6004680, 6004660, 6004650
Case of 50 plates	2 bags of 25 plates (no lids), packed in Styrofoam protection	6004290, 6004620, 6004270, 6004520, 6004640, 6004500, 6004350
Case of 80 plates	4 bags of 20 plates (with lids), packed in Styrofoam protection	6004684, 6004664, 6004654

Dimensions -- Customer Drawing:

