

Research use only. Not for use in diagnostic procedures.

Technical datasheet: Dimensions of ProxiPlate-384 / SW SpectraPlate-384 / AlphaPlate-384 SW

Part numbers:	PN 6008280 / 6008289	white, untreated (MB)
	PN 6008230 / 6008238 / 6008239	white, TC-treated
	PN 6008260 / 6008269	black, untreated (MB)
	PN 6008210 / 6008218 / 6008219	black, TC-treated
	PN 6008640 / 6008649	clear, untreated (MB)
	PN 6008650	clear, TC-treated
	PN 6008350 / 6008359	light gray, untreated (MB)

Material and treatment: Black, white, transparent, or light gray polystyrene microplate with 384 flat-bottom, shallow wells (reduced volume). Plate is available untreated (medium binding (MB)), or Tissue-Culture treated (TC).

Packaging: Plates are placed in diverse formats to provide you the right quality and protection during transport.

Format	Type	Part numbers
Case of 50 plates	Individually pouched (with lids)	6008230, 6008210, 6008650
Case of 50 plates	2 bags of 25 plates (no lids)	6008280, 6008260, 6008640, 6008350
Case of 160 plates	8 bags of 20 plates (with lids)	6008238, 6008218
Case of 200 plates	8 bags of 25 plates (no lids)	6008289, 6008239, 6008269, 6008219, 6008649, 6008359

