

Dimensions of LBS-coated ProxiPlate-384 / SW SpectraPlate-384 / AlphaPlate-384 SW

Part numbers:	LBS ProxiPlate-384	PN 6059480	white, Low Binding Surface - coated
	LBS ProxiPlate-384 F	PN 6059260	black, Low Binding Surface - coated
	LBS SW AlphaPlate-384	PN 6059340	light-gray, Low Binding Surface - coated
	LBS SW SpectraPlate-384	PN 6059630	clear, Low Binding Surface - coated

Material and treatment: Black, white, clear, and light-grey polystyrene microplate with 384 flat-bottom wells with reduced volume ('Shallow Well'). This range of plates is coated to create a Low Binding Surface (LBS) to reduce non-specific binding of material to the plastic surface. Plates with other surface treatments are available.

Packaging: All plates are available in cases with 50 plates, packed as 2 bags of 25.

Dimensions -- Customer Drawing:

