

Research use only. Not for use in diagnostic procedures.

Dimensions of LBS-coated 1/2Area Plates-96

Part numbers:	LBS 1/2Area OptiPlate-96	PN 6052480	white, Low Binding Surface-coated
	LBS 1/2Area OptiPlate-96 F	PN 6052260	black, Low Binding Surface-coated
	LBS 1/2Area AlphaPlate-96	PN 6052340	light-grey, Low Binding Surface-coated
	LBS 1/2Area SpectraPlate-96	PN 6052630	clear, Low Binding Surface-coated

Material and treatment:

Black, white, clear or light-grey polystyrene microplate with 96 flat-bottom wells with reduced diameter / volume. 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 plates.

Dimensions -- Customer Drawing:

