

Research use only. Not for use in diagnostic procedures.

Dimensions of LBS-coated OptiPlate-384 / SpectraPlate-384 / AlphaPlate-384

| | | | |
|----------------------|----------------------|------------|--|
| Part numbers: | LBS OptiPlate-384 | PN 6057480 | white, Low Binding Surface - coated |
| | LBS OptiPlate-384 F | PN 6057260 | black, Low Binding Surface - coated |
| | LBS AlphaPlate-384 | PN 6057340 | light-gray, Low Binding Surface - coated |
| | LBS SpectraPlate-384 | PN 6057630 | clear, Low Binding Surface - coated |

Material and treatment: Black, white, clear, and light-grey polystyrene microplate with 384 flat-bottom wells. 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:

