

## DIMENSIONS OF LBS-COATED PROXIPLATE-384 / SW SPECTRAPLATE-384 / ALPHAPLATE-384 SW

<i>LBS ProxiPlate-384</i>	<i>PN 6059480</i>	<i>white, Low Binding Surface-coated</i>
<i>LBS ProxiPlate-384 F</i>	<i>PN 6059260</i>	<i>black, Low Binding Surface-coated</i>
<i>LBS SW AlphaPlate-384</i>	<i>PN 6059340</i>	<i>light-grey, Low Binding Surface-coated</i>
<i>LBS SW SpectraPlate-384</i>	<i>PN 6059630</i>	<i>clear, Low Binding Surface-coated</i>

### 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 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

