

A research product for research purposes only.

## DIMENSIONS OF LBS-COATED 1/2 AREA PLATES-96

<b>LBS 1/2Area OptiPlate-96</b>	<b>PN 6052480</b>	<i>white, Low Binding Surface-coated</i>
<b>LBS 1/2Area OptiPlate-96 F</b>	<b>PN 6052260</b>	<i>black, Low Binding Surface-coated</i>
<b>LBS 1/2Area AlphaPlate-96</b>	<b>PN 6052340</b>	<i>light-grey, Low Binding Surface-coated</i>
<b>LBS 1/2Area SpectraPlate-96</b>	<b>PN 6052630</b>	<i>clear, Low Binding Surface-coated</i>

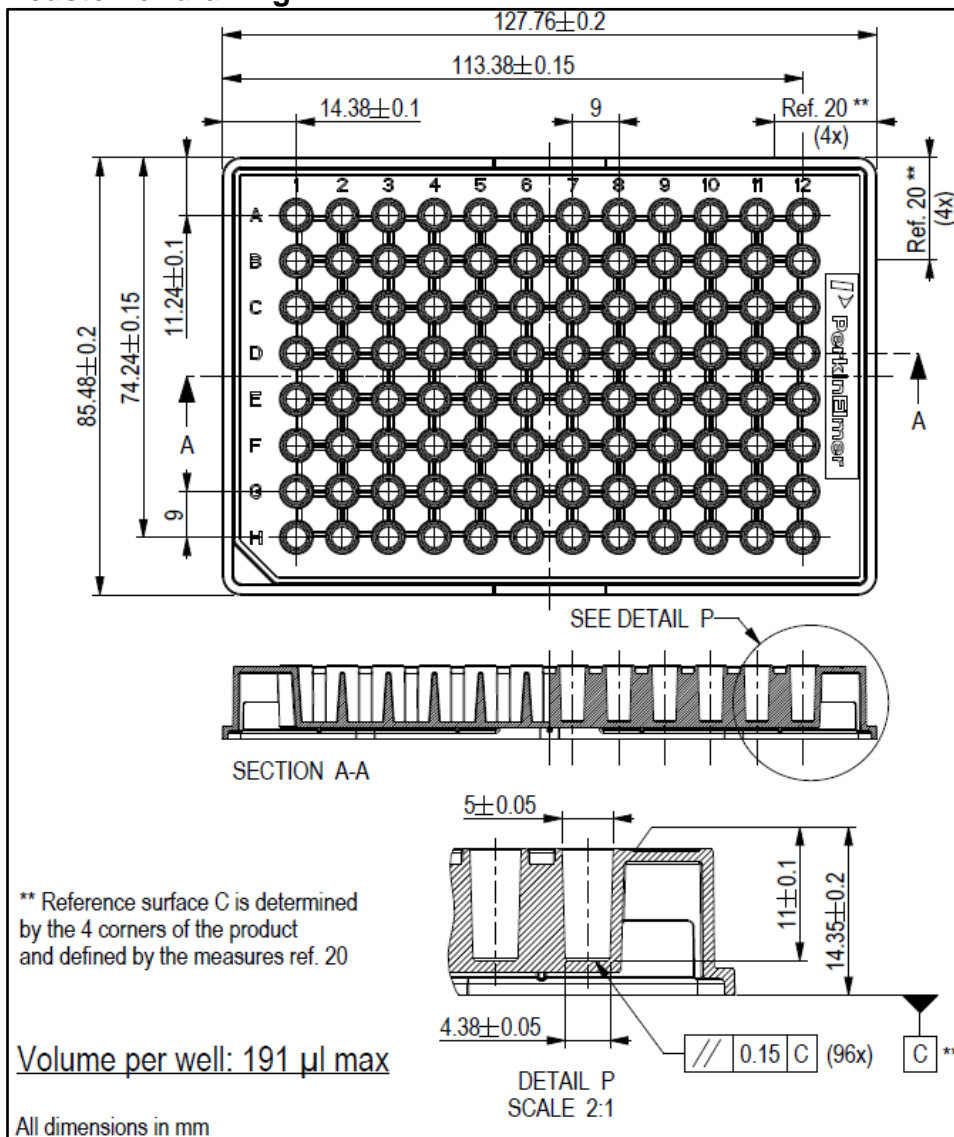
### Material and treatment

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



Product manufactured and distributed by:  
 PerkinElmer Health Sciences B.V.  
 The Netherlands