

## DIMENSIONS OF 1/2 AREA PLATES-96

1/2Area OptiPlate-96	PN 6002290 / 6002299	white, untreated
1/2Area OptiPlate-96 HB	PN 6002520	white, high protein binding affinity
1/2Area OptiPlate-96 F	PN 6002270 / 6002279	black, untreated
1/2Area AlphaPlate-96	PN 6002350 / 6002359	light-grey, untreated
1/2Area SpectraPlate-96 MB	PN 6002640 / 6002649	clear, medium protein binding affinity

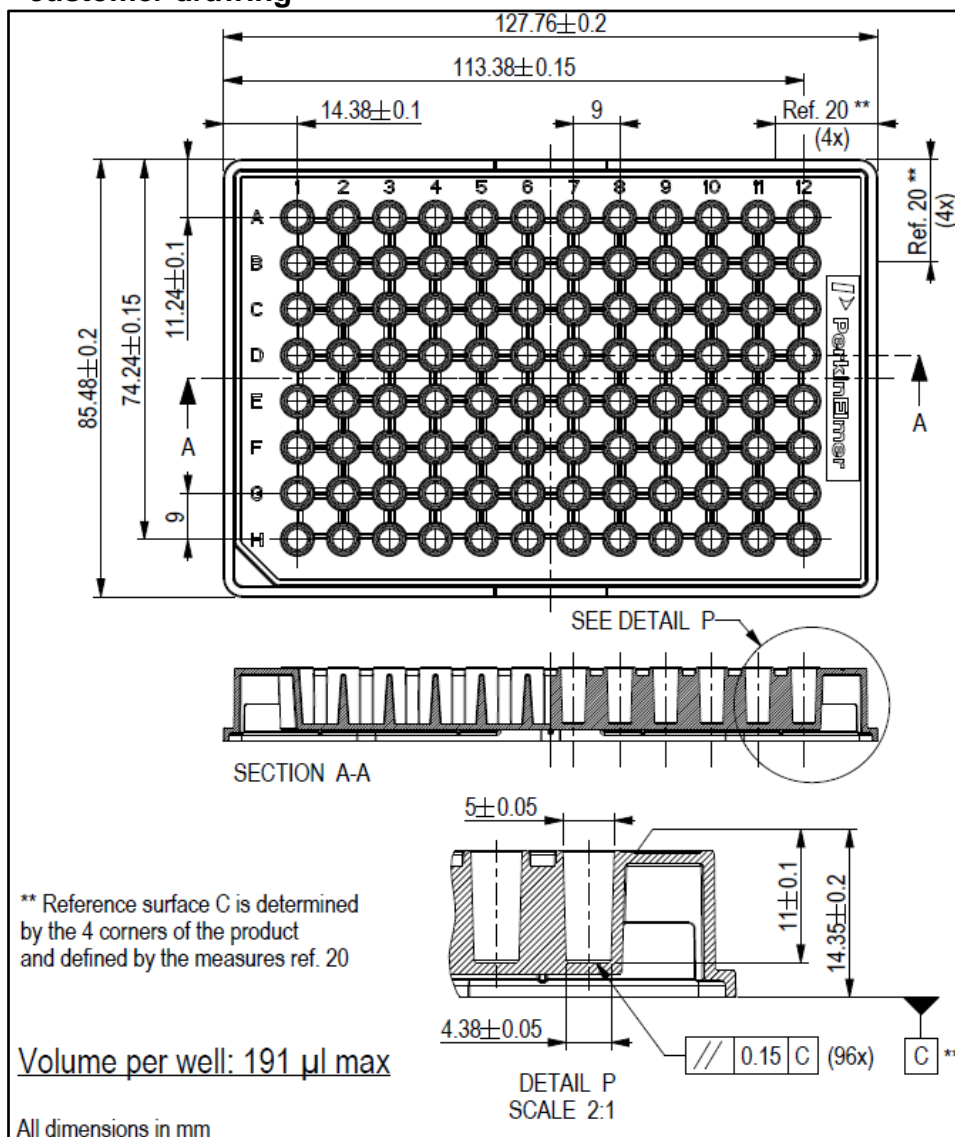
### Material and treatment

Black, white, clear or light-grey polystyrene microplate with 96 flat-bottom wells with reduced diameter / volume. Plate is available untreated, TC-treated or with high protein binding affinity.

### Packaging

6002290/6002520/6002270/ 6002350/6002640 is a case with 50 plates, packed as 2 bags of 25 plates. 6002299/6002279/6002359/6002649 is a box with 200 plates, packed as 8 bags of 25 each.

### Dimensions --- customer drawing



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