

## DIMENSIONS OF ISOPLATE-96, SCINTIPLATE-96 and B&W ISOPLATE-96

### WHITE ISOPLATE-96

*PN 6005040 / 6005049 --- white, untreated*  
*PN 6005590 / 6005599--- white, high protein binding affinity (HB)*  
*PN 6005070 / 6005078 --- white, TC-treated*

### BLACK ISOPLATE-96

*PN 6005020 / 6005029 --- black, untreated*  
*PN 6005570 / 6005579--- black, high protein binding affinity (HB)*  
*PN 6005050 / 6005058 --- black, TC-treated*

### SCINTIPLATE-96

*PN 6005340 / 6005349 --- white with scintillating wells, high protein binding affinity (HB)*  
*PN 6005390 / 6005398 --- white with scintillating wells, TC-treated*  
*PN 1450-551 --- white with scintillating wells, streptavidin-coated*

### B&W ISOPLATE-96

*PN 6005030 / 6005039 --- black with white wells (B&W), untreated*  
*PN 6005580 / 6005589--- black with white wells (B&W), high protein binding affinity (HB)*  
*PN 6005060 / 6005068 --- black with white wells (B&W), TC-treated*

### Material and treatment

Black or white polystyrene plate frame with the inside of the wells made from different material:

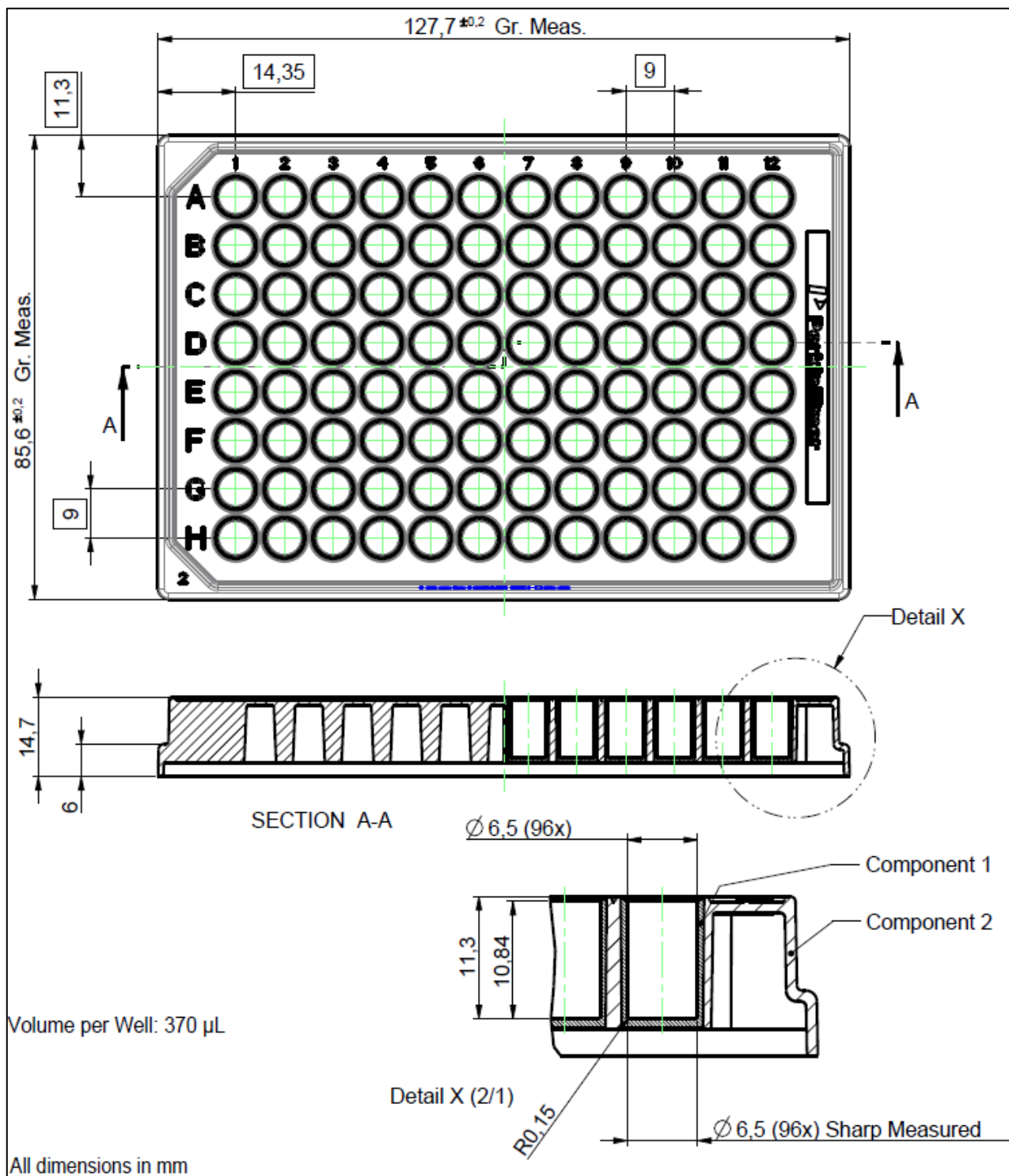
Clear wells = Isoplate-96  
White wells = B&W Isoplate-96  
Scintillating wells = Scintiplate-96

The wells have a flat bottom. Refractive index of clear plate bottom is 1.58. Plate is available untreated, with high protein binding affinity (HB) or Tissue-Culture treated to facilitate cell adherence and proliferation. The Scintiplate-96 is also available with a streptavidin coating.

### Packaging

| format             | type                             | part numbers  |
|--------------------|----------------------------------|---|
| Box of 10 plates   | Individually pouched             | 1450-551  |
| Case of 50 plates  | Individually pouched (with lids) | 6005070, 6005050, 6005390, 6005060                            |
| Case of 50 plates  | 2 bags of 25 plates (no lids)    | 6005040, 6005590, 6005020, 6005570, 6005340, 6005030, 6005580 |
| Case of 160 plates | 8 bags of 20 plates (with lids)  | 6005078, 6005058, 6005398, 6005068                            |
| Case of 200 plates | 8 bags of 25 plates (no lids)    | 6005049, 6005599, 6005029, 6005579, 6005349, 6005039, 6005589 |

# Dimensions --- customer drawing



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