

Research use only. Not for use in diagnostic procedures.

Dimensions of Isoplate-96, ScintiPlate-96 and B&W Isoplate-96

Part numbers:

WHITE ISOPLATE-96	PN 6005040 / 6005049 PN 6005590 / 6005599 PN 6005070 / 6005078	White, untreated White, high protein binding affinity (HB) White, Tissue Culture-treated (TC)
BLACK ISOPLATE-96	PN 6005020 / 6005029 PN 6005570 / 6005579 PN 6005050	Black, untreated Black, HB affinity Black, TC-treated
SCINTIPLATE-96	PN 6005340 / 6005349 PN 6005390	White with scintillating wells, HB affinity White with scintillating wells, TC-treated
B&W ISOPLATE-96	PN 6005030 / 6005039 PN 6005580 / 6005589 PN 6005060 / 6005068	Black with white wells (B&W), untreated Black with white wells (B&W), HB affinity Black with white wells (B&W), TC-treated

Material and treatment or coating:

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 (TC) to facilitate cell adherence and proliferation.

Packaging:

Format	Type	Part numbers
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, 6005068
Case of 200 plates	8 bags of 25 plates (no lids)	6005049, 6005599, 6005029, 6005579, 6005349, 6005039, 6005589

Dimensions --- Customer Drawing:

