

DIMENSIONS OF OPTIPLATE-1536 / CULTURPLATE-1536 / SPECTRAPLATE-1536 / ALPHAPLATE-1536

- PN 6004290 --- white, untreated*
- PN 6004620 --- white, high protein binding affinity (HB)*
- PN 6004680 / 6004684 --- white, TC-treated*
- PN 6004270 --- black, untreated*
- PN 6004520 --- black, high protein binding affinity (HB)*
- PN 6004660 / 6004664 --- black, TC-treated*
- PN 6004640 --- clear, untreated (MB)*
- PN 6004500 --- clear, high protein binding affinity (HB)*
- PN 6004650 / 6004654 --- clear, TC-treated*
- PN 6004350 --- light gray, untreated*
- PN 6004360 --- dark gray (HS), untreated*

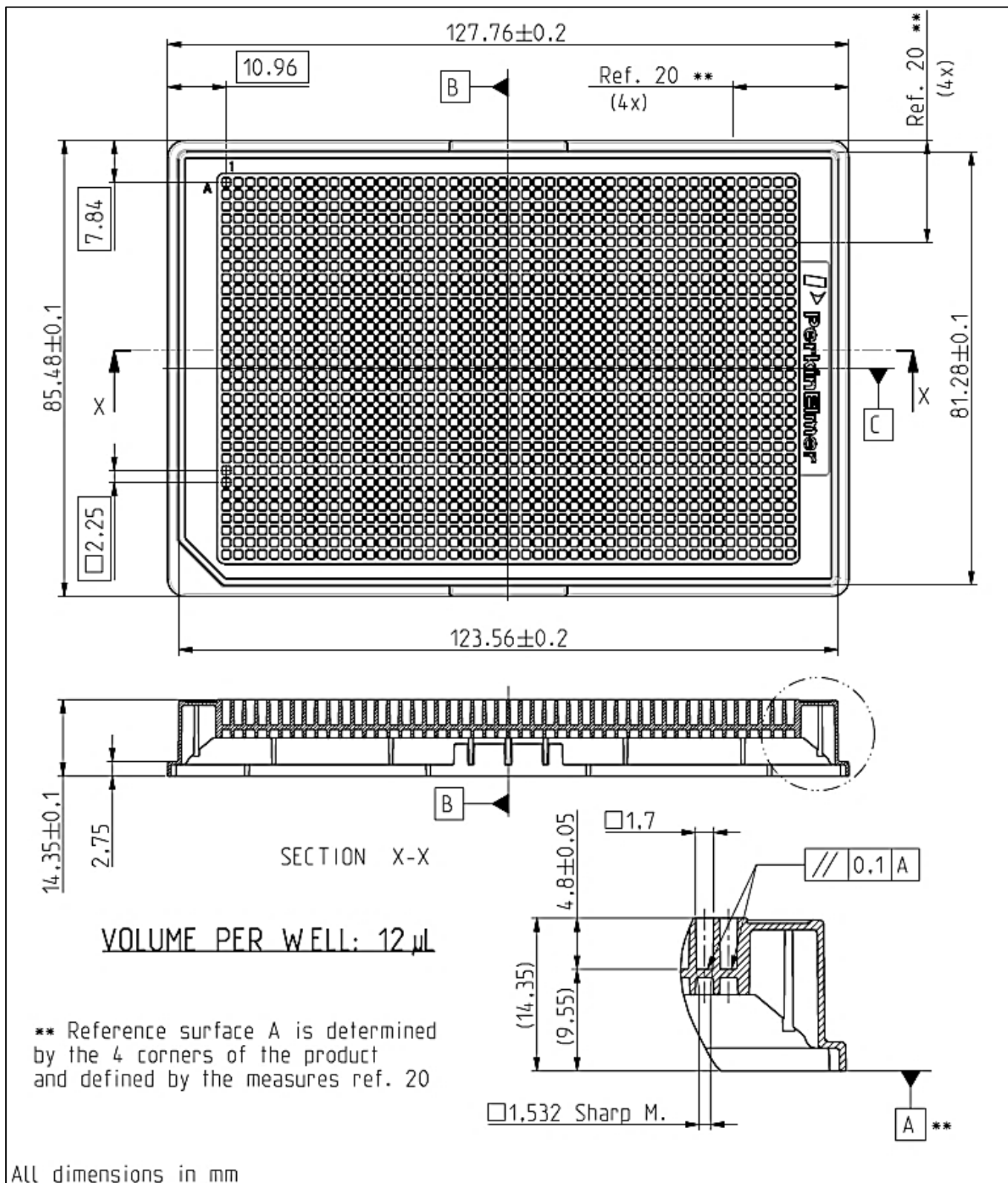
Material and treatment

Black, white, transparent, light gray or dark gray polystyrene microplate with 1536 flat-bottom wells. Plate is available untreated, with high protein binding affinity or Tissue-Culture treated.

Packaging

format	type	part numbers
Case of 50 plates	Individually pouched (with lids)	6004680, 6004660, 6004650
Case of 50 plates	2 bags of 25 plates (no lids), packed in styrofoam protection	6004290, 6004620, 6004270, 6004520, 6004640, 6004500, 6004350, 6004360
Case of 80 plates	4 bags of 20 plates (with lids), packed in Styrofoam protection	6004684, 6004664, 6004654

Dimensions --- customer drawing



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