

DIMENSIONS OF OPTIPLATE-384 / CULTURPLATE-384 / SPECTRAPLATE-384 / ALPHAPLATE-384

- PN 6007290 / 6007299 --- white, untreated*
PN 6005620 / 6005629 --- white, high protein binding affinity (HB)
PN 6007680 / 6007688 / 6007689 --- white, TC-treated
PN 6007270 / 6007279 --- black, untreated
PN 6005520 / 6005529 --- black, high protein binding affinity (HB)
PN 6007660 / 6007668 / 6007669 --- black, TC-treated
PN 6007640 / 6007649 --- clear, untreated (MB)
PN 6007500 / 6007509 --- clear, high protein binding affinity (HB)
PN 6007650 / 6007658 / 6007659 --- clear, TC-treated
PN 6005350 / 6005359 --- light gray, untreated
PN 6057690 / 6057699 --- light gray, high protein binding affinity (HB)

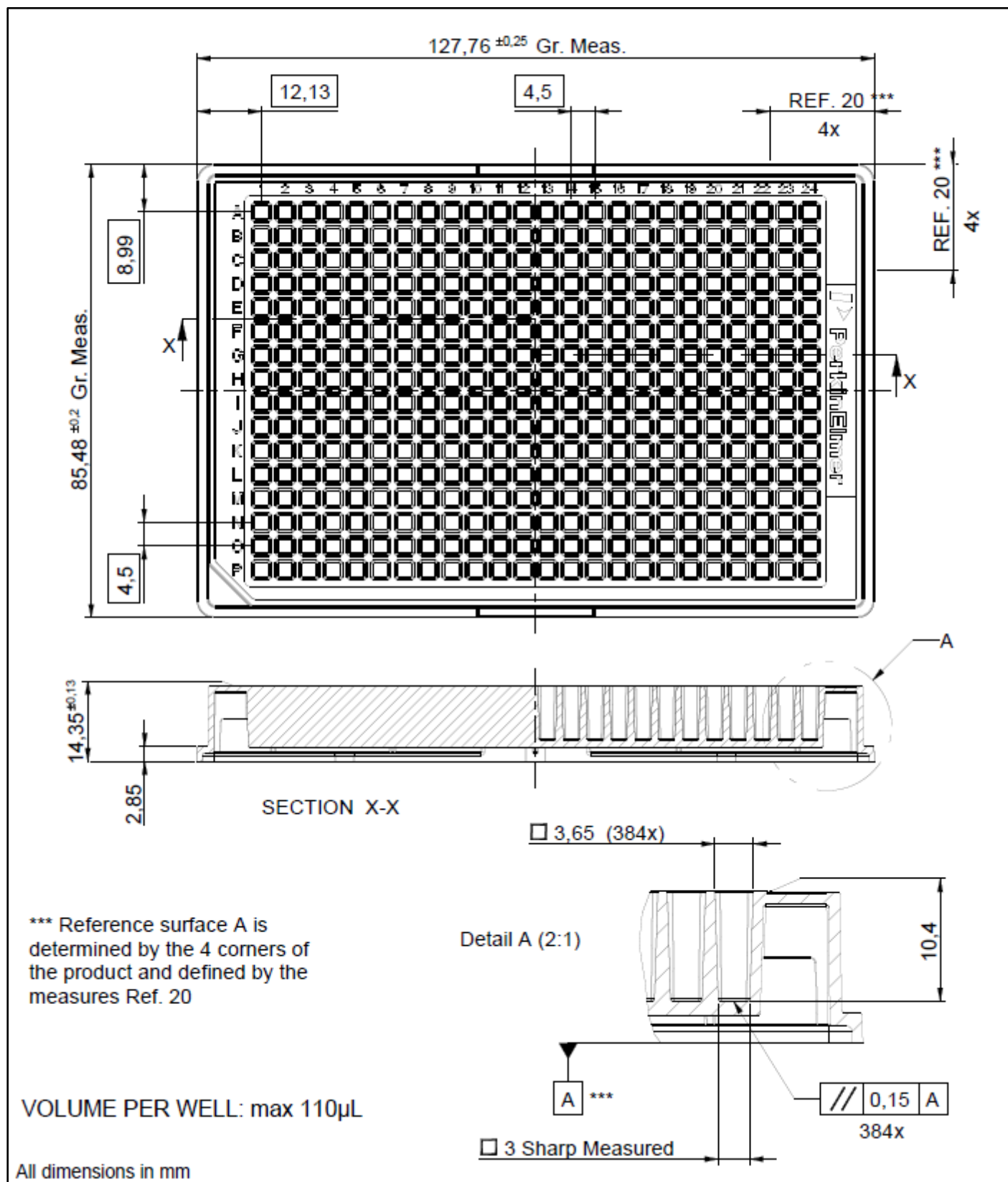
Material and treatment

Black, white, transparent, or light gray polystyrene microplate with 384 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) | 6007680, 6007660, 6007650 |
| Case of 50 plates | 2 bags of 25 plates (no lids) | 6007290, 6005620, 6007270, 6005520, 6007640, 6007500, 6005350, 6057690 |
| Case of 160 plates | 8 bags of 20 plates (with lids) | 6007688, 6007668, 6007658 |
| Case of 200 plates | 8 bags of 25 plates (no lids) | 6007299, 6005629, 6007689, 6007279, 6005529, 6007669, 6007649, 6007509, 6007659, 6005359, 6057699 |

Dimensions --- customer drawing



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