

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

(PN 6007290, 6007299, 6007270, 6007279, 6007680, 6007688, 6007689, 6007660, 6007688, 6007689, 6005350, 6005359, 6007640, 6007649, 6007500, 6007509, 6007650, 6007658, 6007559)

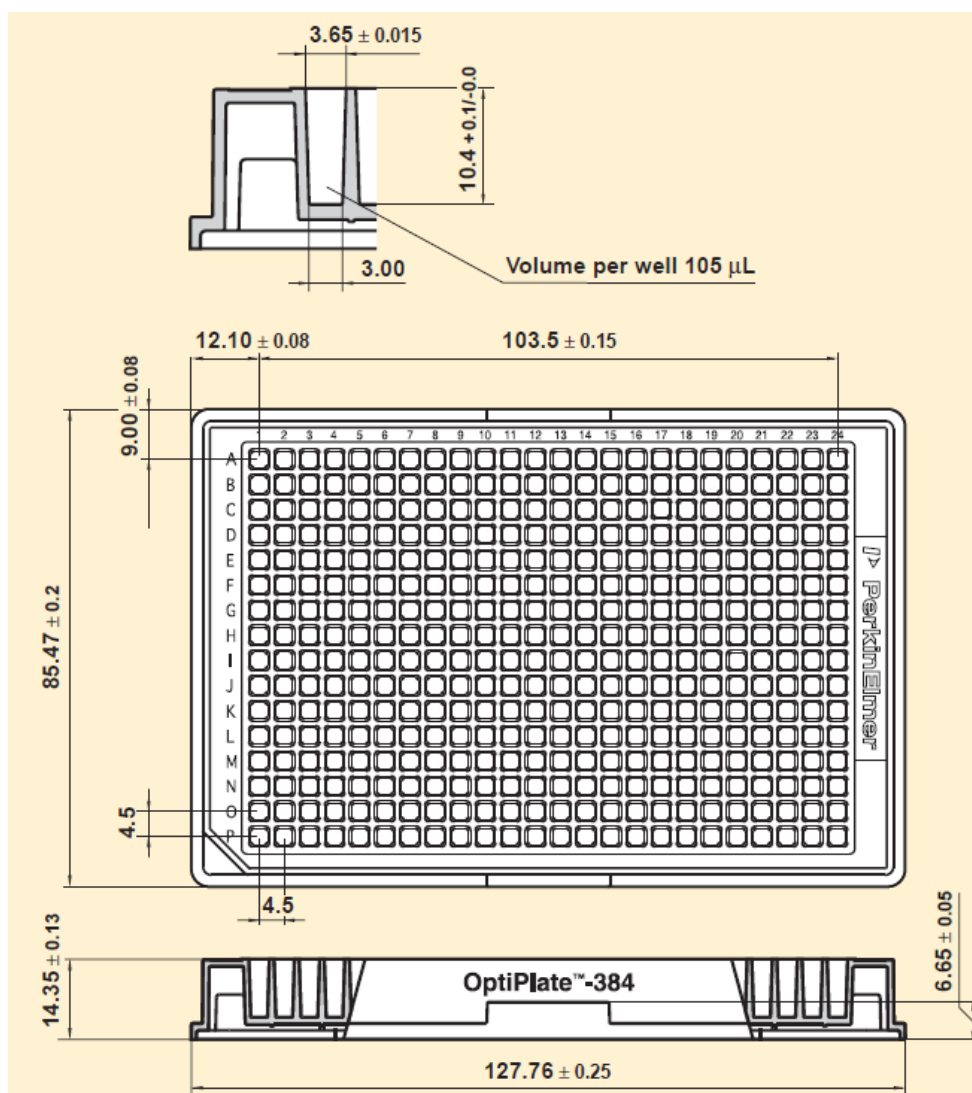
Material and treatment

White, clear, black or grey microplate with 384 flat-bottom wells. Plates are available in untreated version, with high protein binding affinity (HB) and as tissue culture treated (TC), lidded and irradiated.

Packaging

Untreated and HB plates are packed in cases of 50 plates (2 bags of 25 plates each) or 200 plates (8 x 25). TC-treated plates are available in cases of 50 plates (individually pouched) or cases of 160 plates (8 bags of 20 lidded plates each).

Dimensions --- customer drawing



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