

DIMENSIONS OF PROXIPLATE-384 / SW SPECTRAPLATE-384 / ALPHAPLATE-384 SW

PN 6008280 / 6008289 --- white, untreated
PN 6008230 / 6008238 / 6008239 --- white, TC-treated
PN 6008260 / 6008269 --- black, untreated
PN 6008210 / 6008218 / 6008219 --- black, TC-treated
PN 6008270 / 6008279 --- dark gray (HS), untreated
PN 6008640 / 6008649 --- clear, untreated (MB)
PN 6008650 --- clear, TC-treated
PN 6008350 / 6008359 --- light gray, untreated

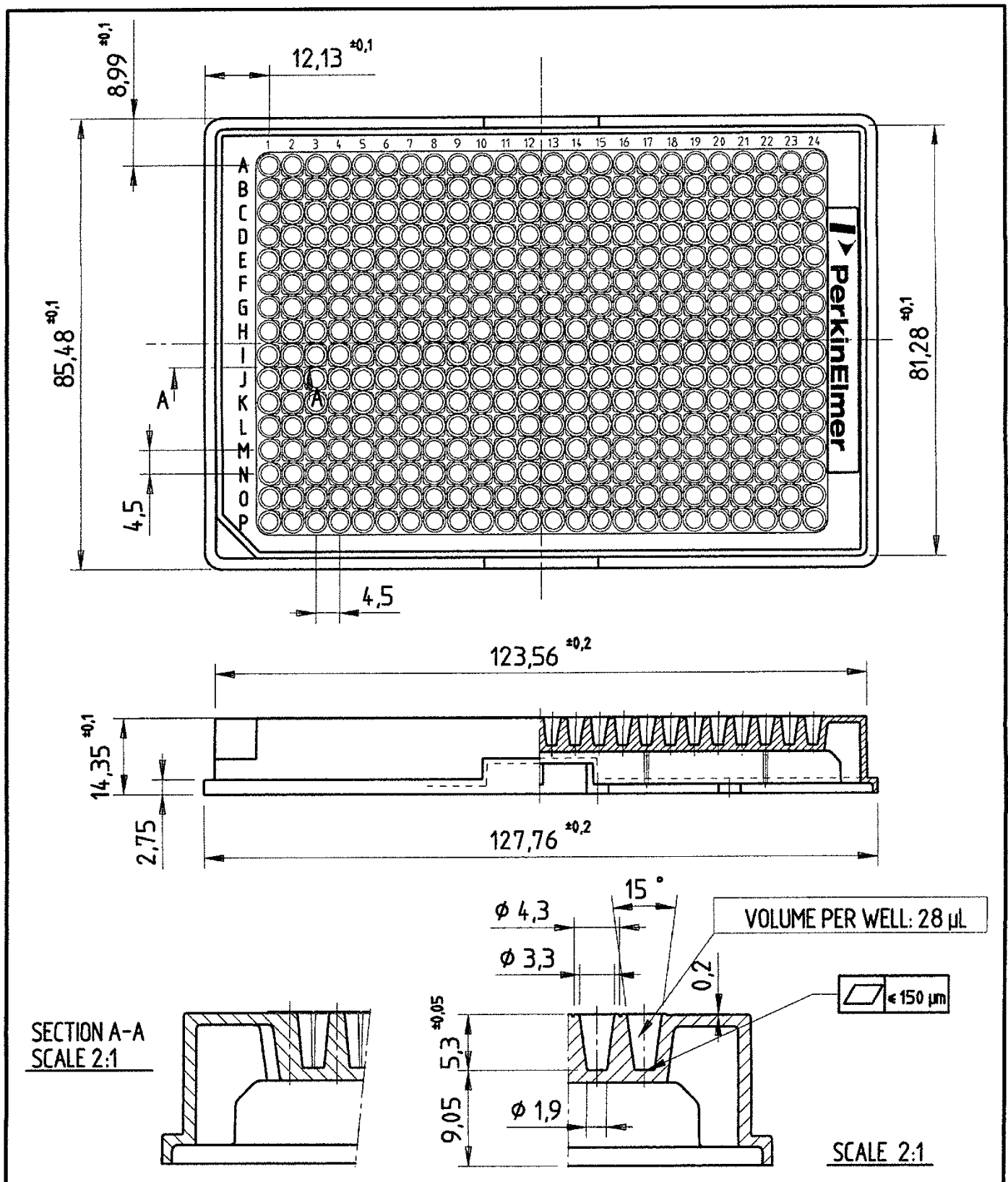
Material and treatment

Black, white, transparent, light gray or dark gray polystyrene microplate with 384 flat-bottom, shallow wells (reduced volume). Plate is available untreated or Tissue-Culture treated.

Packaging

format	type	part numbers
Case of 50 plates	Individually pouched (with lids)	6008230, 6008210, 6008650
Case of 50 plates	2 bags of 25 plates (no lids)	6008280, 6008260, 6008270, 6008640, 6008350
Case of 160 plates	8 bags of 20 plates (with lids)	6008238, 6008218
Case of 200 plates	8 bags of 25 plates (no lids)	6008289, 6008239, 6008269, 6008219, 6008279, 6008649, 6008359

Dimensions --- customer drawing



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