

Research use only. Not for use in diagnostic procedures.

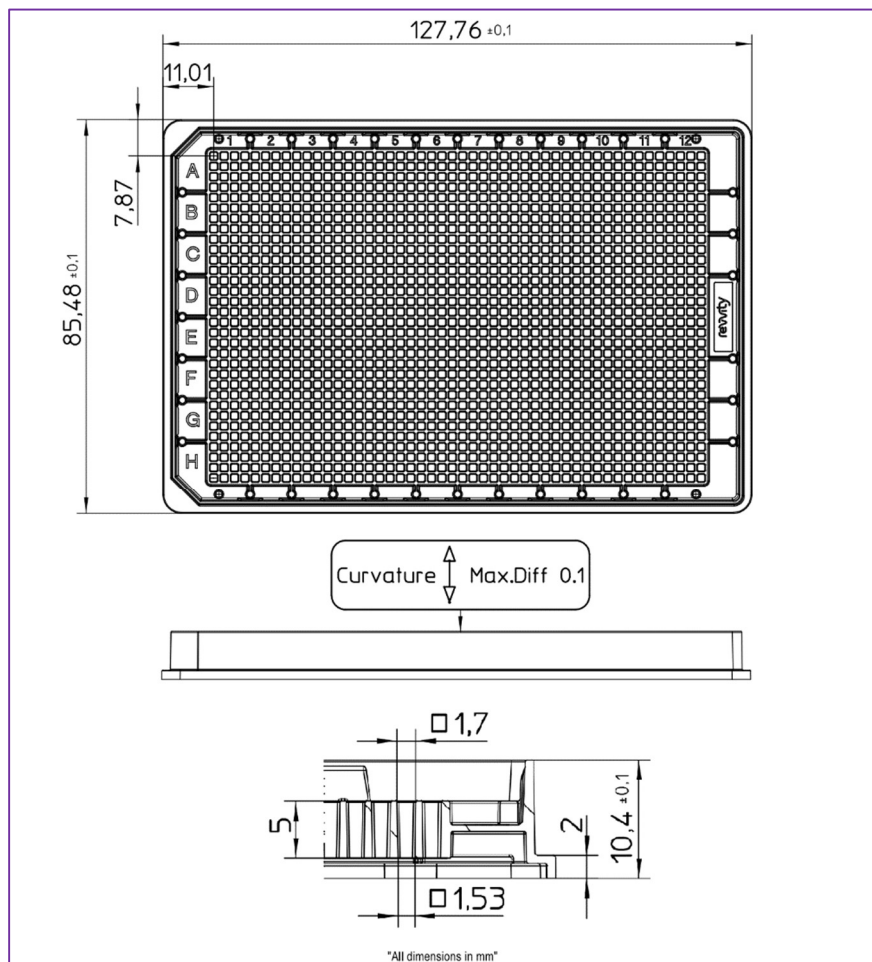
Dimensions of ViewPlate-1536

Part numbers:	PN 6004470	Black, untreated
	PN 6004490	White, untreated
	PN 6004460	Black, TC-treated
	PN 6004480	White, TC-treated
	PN 6004810	Black, Collagen-coated
	PN 6004710	Black, PDL-coated

Material and treatment: Black or white polystyrene plate frame with transparent polystyrene foil bottom (75 µm). Refractive index of plate bottom is 1.58. Plate is available untreated, Tissue-Culture treated (TC), or with PDL or Collagen-coating to facilitate cell adherence and proliferation.

Packaging: 6004470 and 6004490 is a case with 60 untreated plates as 6 bags of 10 plates each. 6004460 and 6004480 is a box with 40 TC-treated plates (4 bags of 10 lidded plates each). 6004810 and 6004710 are boxes with 10 coated and lidded plates, individually packed.

Dimensions --- Customer Drawing:



Dimensions for Harmony® Software Microplate Definition Wizard

Plate Type Definition Wizard

Step 4 / 7

Please enter the dimensions of the plate type.

A: mm C: mm E: mm
 B: mm D: mm F: mm

< Back Next >

Plate Type Definition Wizard

Step 5 / 7

Side:

G: mm
 H: mm
 I: mm

Top: Bottom:

J: mm N: mm

< Back Next >

revvity