

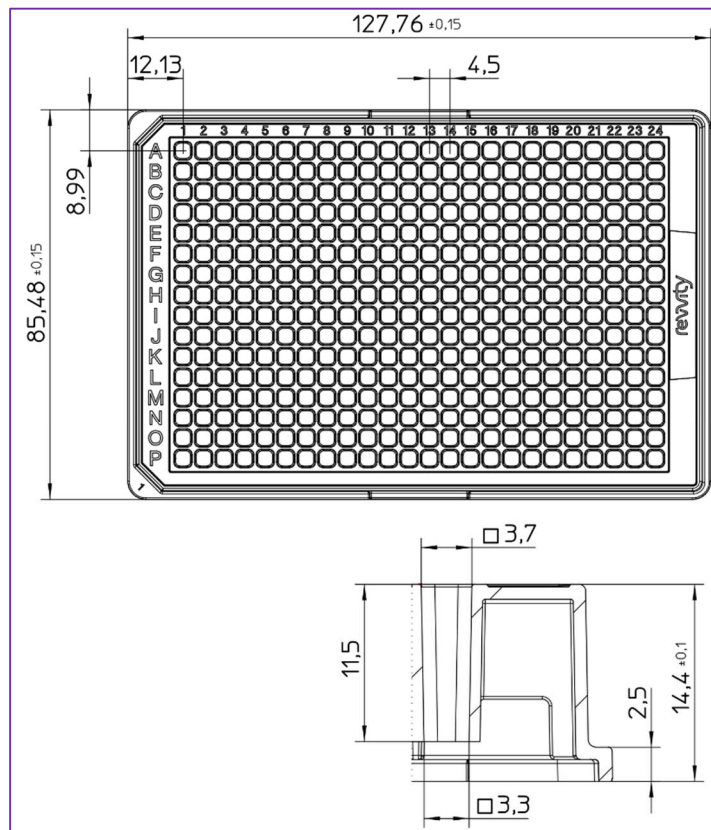
Dimensions of ViewPlate-384

Part numbers:	PN 6007470	Black, untreated
	PN 6007490	White, untreated
	PN 6007460	Black, TC-treated
	PN 6007480	White, TC-treated
	PN 6007810 / 6007818	Black, Collagen-coated
	PN 6007710 / 6007718	Black, PDL-coated

Material and treatment: Black or white polystyrene plate frame with transparent polystyrene foil bottom (190 µm). Refractive index of plate bottom is 1.58. Plate is available untreated, Tissue-Culture (TC) treated, or with PDL or Collagen-coating to facilitate cell adherence and proliferation.

Packaging: 6007470 and 6007490 is a case with 50 untreated plates as 2 bags of 25 plates each. 6007460 and 6007480 is a box with 40 TC-treated plates (4 bags of 10 lidded plates each). 6007810 and 6007710 are boxes with 40 coated and lidded plates, packed as 2 bags of 20 each. 6007818 and 6007718 hold 160 coated and lidded plates as 8 bags of 20.

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:	<input type="text" value="12.13"/>	mm	C:	<input type="text" value="127.76"/>	mm	E:	<input type="text" value="115.63"/>	mm
B:	<input type="text" value="8.99"/>	mm	D:	<input type="text" value="85.48"/>	mm	F:	<input type="text" value="76.49"/>	mm

< Back Next >

Plate Type Definition Wizard

Step 5 / 7

Side:

G:	<input type="text" value="14.40"/>	mm
H:	<input type="text" value="2.71"/>	mm
I:	<input type="text" value="0.19"/>	mm

Top: Bottom:

J:	<input type="text" value="3.70"/>	mm	N:	<input type="text" value="3.30"/>	mm
----	-----------------------------------	----	----	-----------------------------------	----

< Back Next >

