

Research use only. Not for use in diagnostic procedures.

Technical datasheet: Dimensions of ViewPlate-96

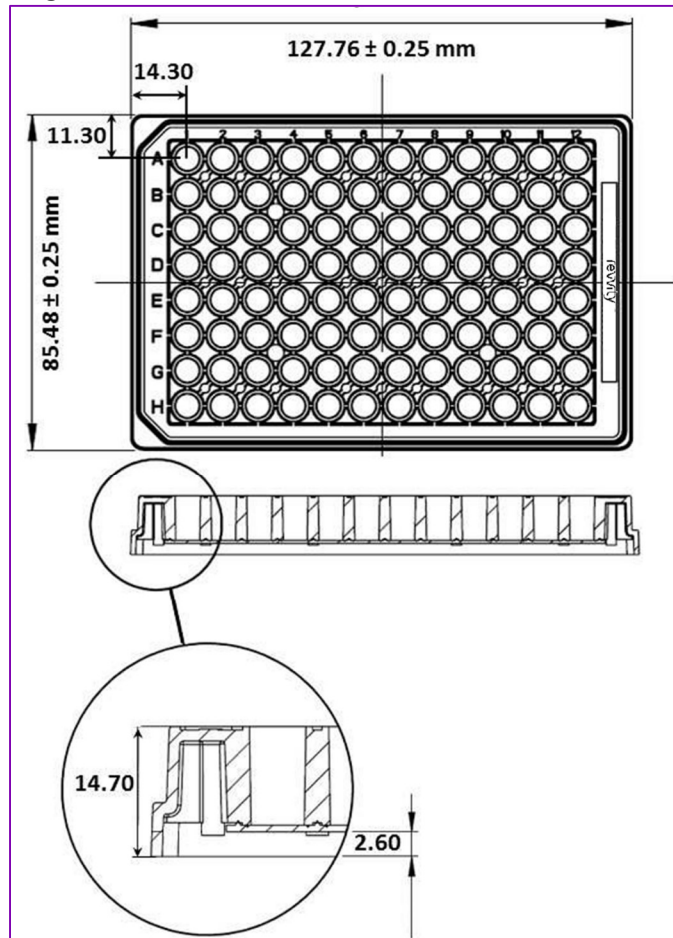
Part numbers: PN 6005182 / 6005225 black, TC-treated
 PN 6005181 white, TC-treated
 PN 6005810 / 6005818 black, Collagen-coated
 PN 6005710 black, PDL-coated

Material and treatment: Black or white polystyrene plate frame with thicker transparent polystyrene bottom (980 µm). Refractive index of plate bottom is 1.58. Plate is available Tissue-Culture treated and PDL or Collagen (COL)-coated to facilitate cell adherence and proliferation.

Packaging:

Format	Type	Part numbers
Case of 50 plates	Individually pouched (with lids)	6005181 / 6005182
Case of 50 plates	10 bags of 5 plates (with lids)	6005225
Case of 40 coated plates (PDL/COL)	2 bags of 20 plates (with lids)	6005810 / 6005710
Case of 160 coated plates (COL)	8 bags of 20 plates (with lids)	6005818

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="14.30"/> mm	C:	<input type="text" value="127.76"/> mm	E:	<input type="text" value="113.30"/> mm
B:	<input type="text" value="11.30"/> mm	D:	<input type="text" value="85.48"/> mm	F:	<input type="text" value="74.30"/> mm

Plate Type Definition Wizard

Step 5 / 7

Side:

G:	<input type="text" value="14.70"/> mm
H:	<input type="text" value="2.60"/> mm
I:	<input type="text" value="0.98"/> mm

Top:

Bottom:

J:	<input type="text" value="6.50"/> mm	N:	<input type="text" value="6.10"/> mm
----	--------------------------------------	----	--------------------------------------

revvity