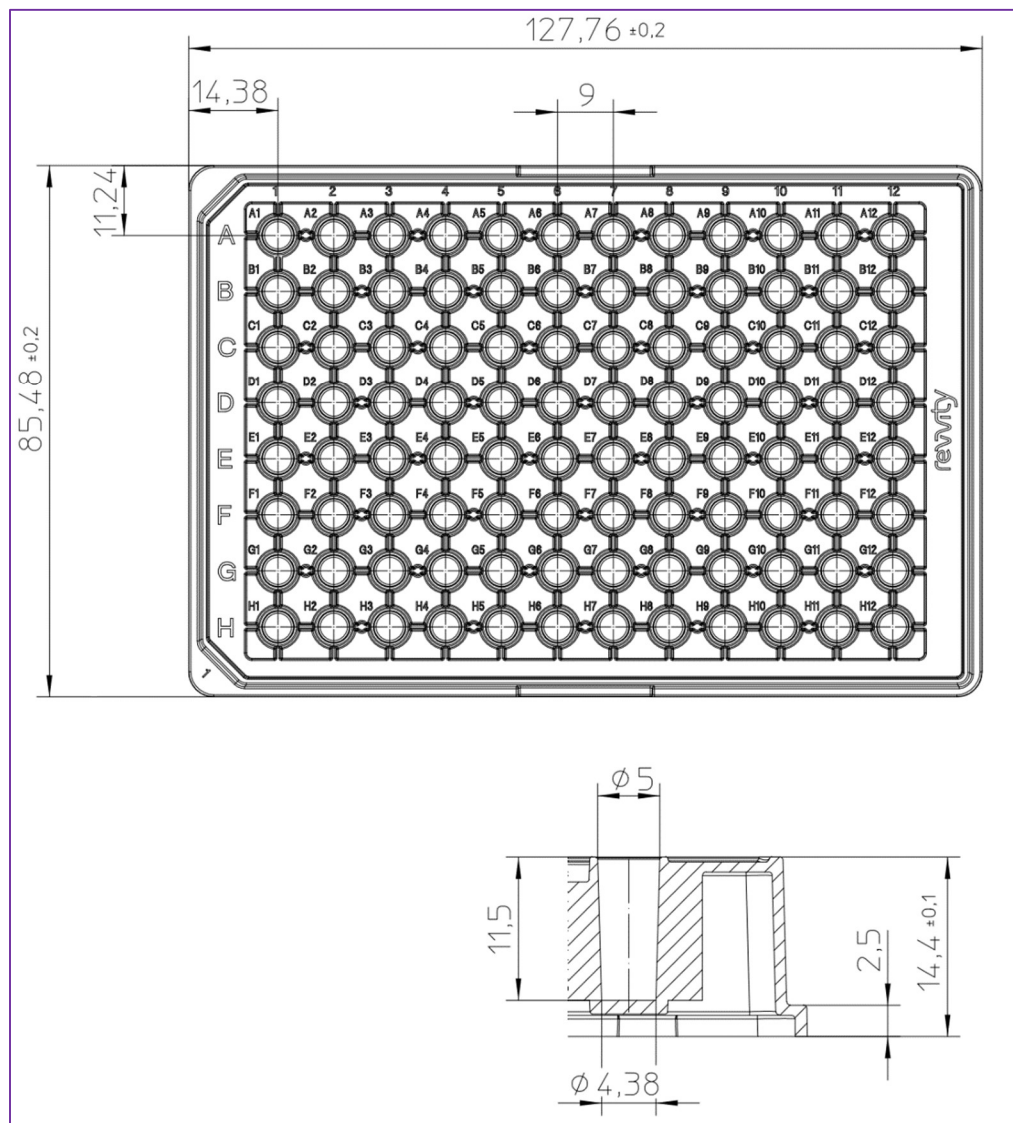


Dimensions of 1/2Area ViewPlate-96

- Part numbers:** PN 6005760 / 6005768 White, TC-treated
- Material and treatment:** White polystyrene plate frame with transparent polystyrene film bottom (190 µm). Refractive index of plate bottom is 1.58. Plate is available Tissue-Culture (TC) treated.
- Packaging:** 6005760 is a case with 40 TC-treated, lidded plates packed as 2 bags of 20 plates each. 6005768 holds 160 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="14.38"/>	mm	C:	<input type="text" value="127.76"/>	mm	E:	<input type="text" value="113.38"/>	mm
B:	<input type="text" value="11.24"/>	mm	D:	<input type="text" value="85.48"/>	mm	F:	<input type="text" value="74.24"/>	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="5.00"/>	mm	N:	<input type="text" value="4.38"/>	mm
----	-----------------------------------	----	----	-----------------------------------	----

< Back Next >

