

Research use only. Not for use in diagnostic procedures.

Dimensions of VisiPlate-24

Part numbers:	PN 1450-602	White, untreated
	PN 1450-603 / 1450-604	White, TC-treated
	PN 1450-605 / 1450-606	Black, TC-treated

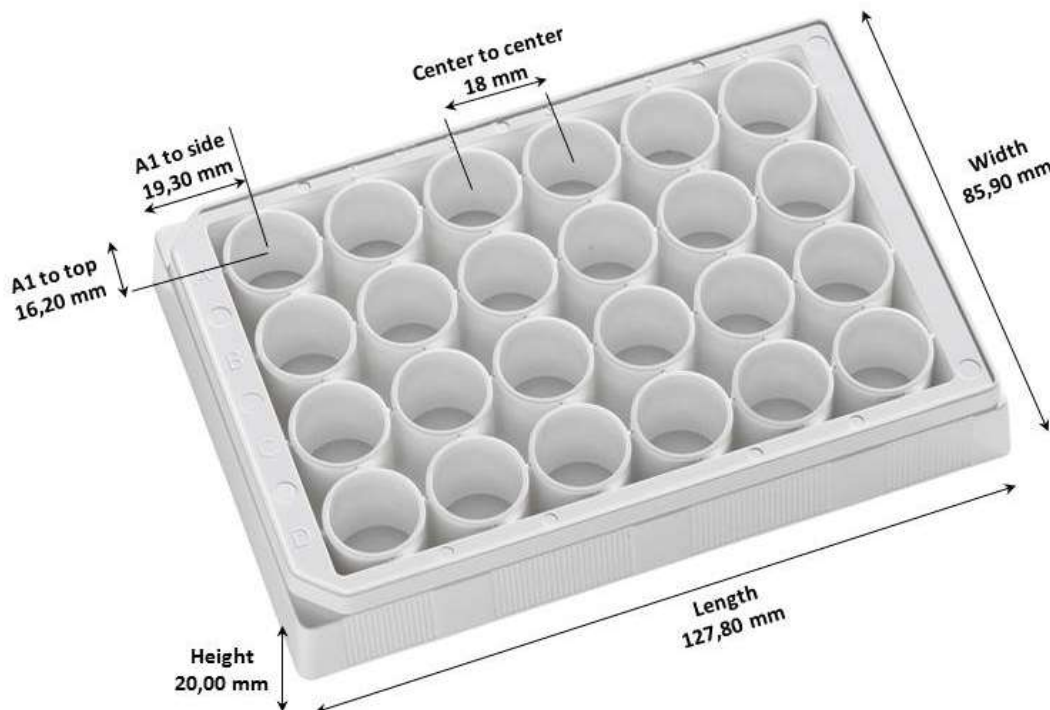
Material and treatment or coating:

Black or white polystyrene plate frame with thicker transparent polystyrene bottom (920 µm). Refractive index of plate bottom is 1.58. Plate is available untreated (white only) or Tissue-Culture treated to facilitate cell adherence and proliferation.

Packaging:

1450-602 is an over-case with 4 boxes of 17 plates each. 1450-603 and 1450-605 is a box with 14 lidded plates that are individually pouched. 1450-604 and 1450-606 is an over-case with 4 boxes containing 14 lidded plates each (individually pouched).

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