

DIMENSIONS OF CellCarrier-96 Ultra

PN 6055300 / 6055302 / 6055308 --- TC-treated
PN 6055700 / 6055708 --- Collagen-coated
PN 6055500 / 6055508 --- PDL-coated
PN 6055800 / 6055802 --- Ultra Low Attachment (ULA)-coated

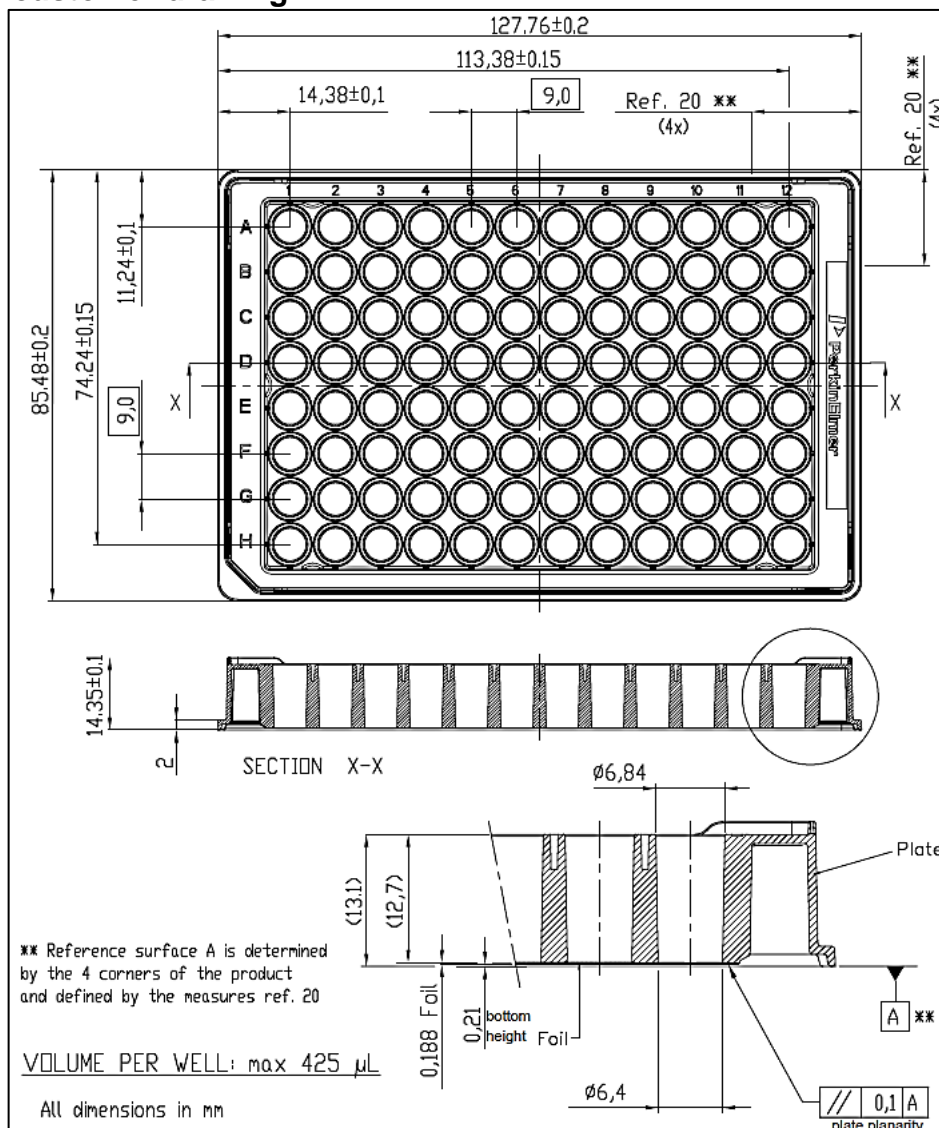
Material and treatment

Low bottom height ($200 \pm 10 \mu\text{m}$), black cyclic olefin microplate with optically clear cyclic olefin foil bottom ($188 \mu\text{m}$). Refractive index of plate bottom is 1.53. Plates are Tissue-Culture treated, PDL- or COL-coated to facilitate cell adherence and proliferation. The ULA-coated plates prevent all cell adherence and are intended for studying 3D spheroids or suspension cells.

Packaging

6055300 is a case with 50 single pouched plates with lids. 6055302, 6055802, 6055700 and 6055500 are boxes with 40 lidded plates, packed as 2 bags of 20 each. 6055308, 6055708 and 6055508 hold 160 lidded plates as 8 bags of 20. 6055800 holds 10 single pouched plates.

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 >

Product manufactured and distributed by:
 PerkinElmer Health Sciences B.V.
 The Netherlands