

## DIMENSIONS OF CellCarrier-96 Ultra

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

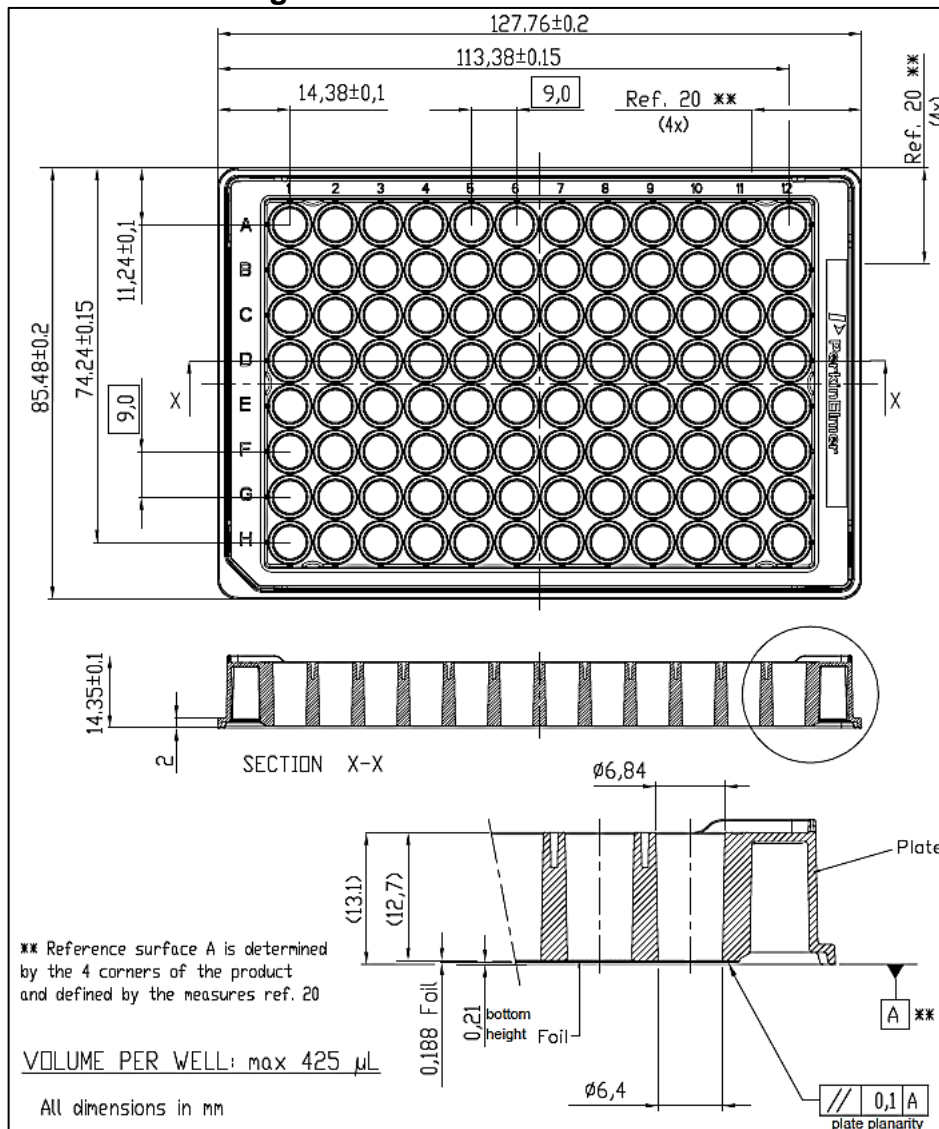
### Material and treatment

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 box with 50 individually pouched plates with lids. 6055302, 6055700 and 6055500 are boxes with 40 lidded plates, packed as 2 bags of 20 each. 6055308, 6055708 and 6055508 holds 160 lidded plates as 8 bags of 20. 6055800 holds 10 single-pouched, lidded 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