

Research use only. Not for use in diagnostic procedures.

## Technical datasheet: Dimensions of CellCarrier Spheroid ULA 96-well Microplate

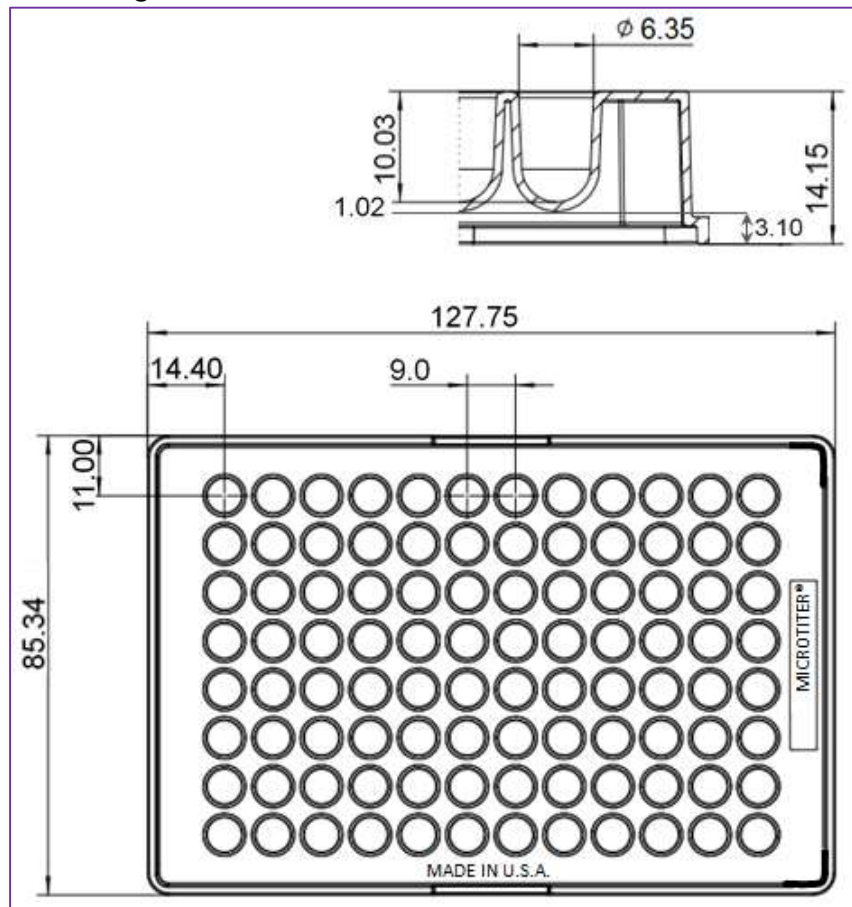
**Part numbers:** PN 6055331 / 6055330 / 6055334 Clear, ULA-treated round-bottom microplates

**Material and treatment:** Round bottom, clear 96-well polystyrene microplate. Plates are coated to turn into Ultra Low Attachment ('ULA') surface for mammalian cells. This ULA coating in round-bottom wells will facilitate formation of round spheroids after incubating seeded cells.

**Packaging:**

Format	Type	Part numbers
Case of 2 plates (Sample Pack)	Individually pouched (with lids)	6055331
Case of 10 plates	Individually pouched (with lids)	6055330
Case of 40 plates	2 bags of 20 plates (with lids)	6055334

**Dimensions -- Customer Drawing:**



MICROTITER is a registered trademark of Thermo Scientific

### 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: (round)

J:  mm

N:  mm

< Back    Next >

