

DIMENSIONS OF CELLCARRIER SPHEROID ULA 96 (PN 6055331 / 6055330 / 6055334)

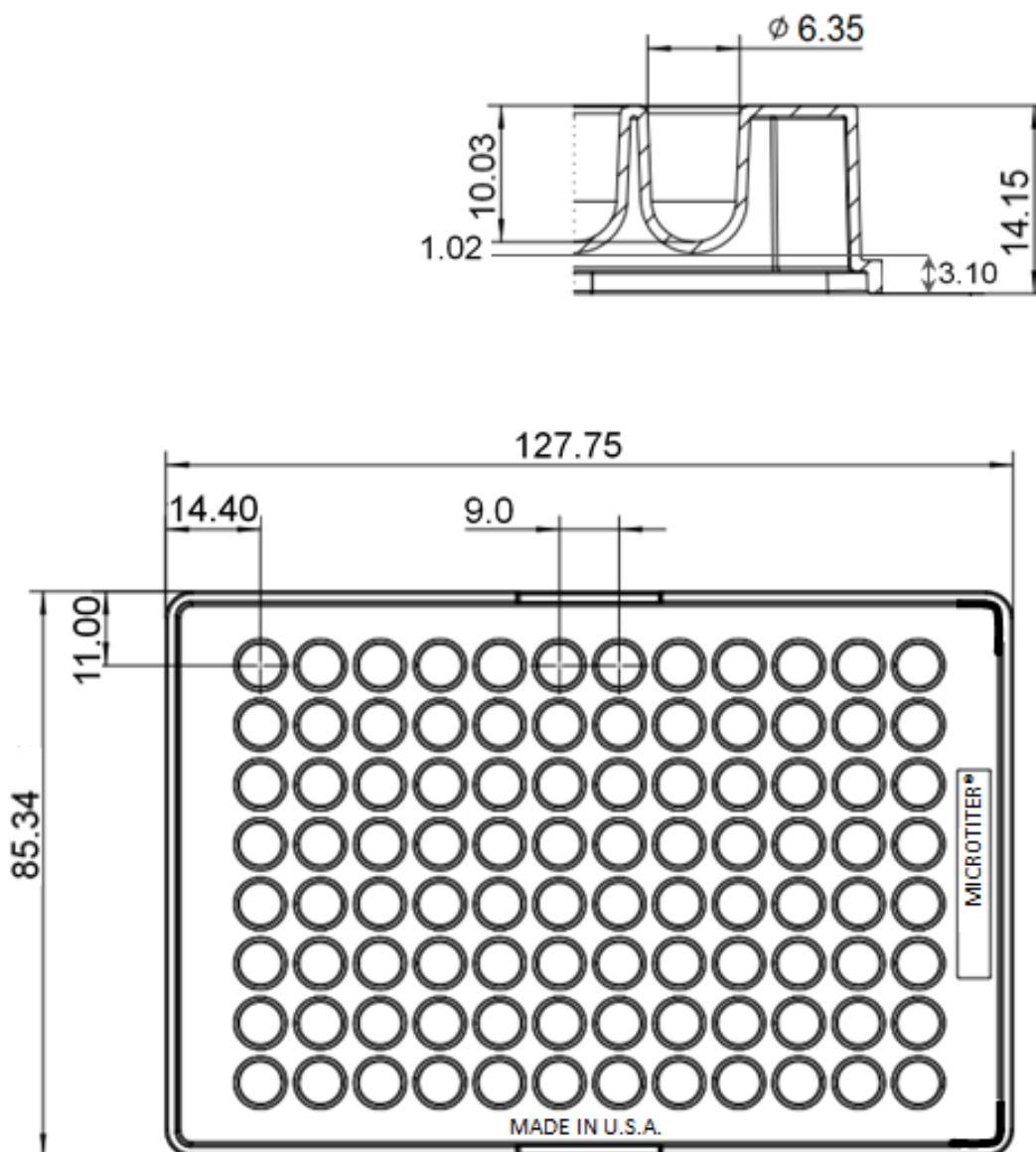
Material and treatment

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

6055331 is a sample pack with 2 individually pouched plates with lids. 6055330 is a pack of 10 individually pouched plates and 6055334 is a case with 40 lidded plates packed as 2 bags with 20 plates each.

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: **14.40** mm C: **127.75** mm E: **113.50** mm
 B: **11.00** mm D: **85.34** mm F: **74.15** mm

< Back Next >

Plate Type Definition Wizard

Step 5 / 7

Side:

G: **14.15** mm
 H: **3.10** mm
 I: **1.02** mm

Top: Bottom: (round)

J: **6.35** mm

N: **0.60** mm

< Back Next >

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