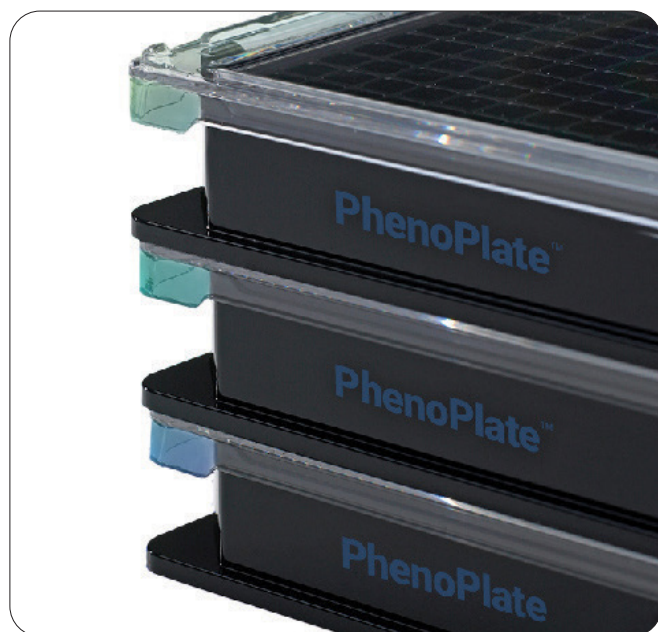


PhenoPlate 96-well microplates

Overview

Revvity PhenoPlates™, formerly known as CellCarrier™ Ultra microplates, have been engineered to deliver superior images and the highest quality data for high-content applications.

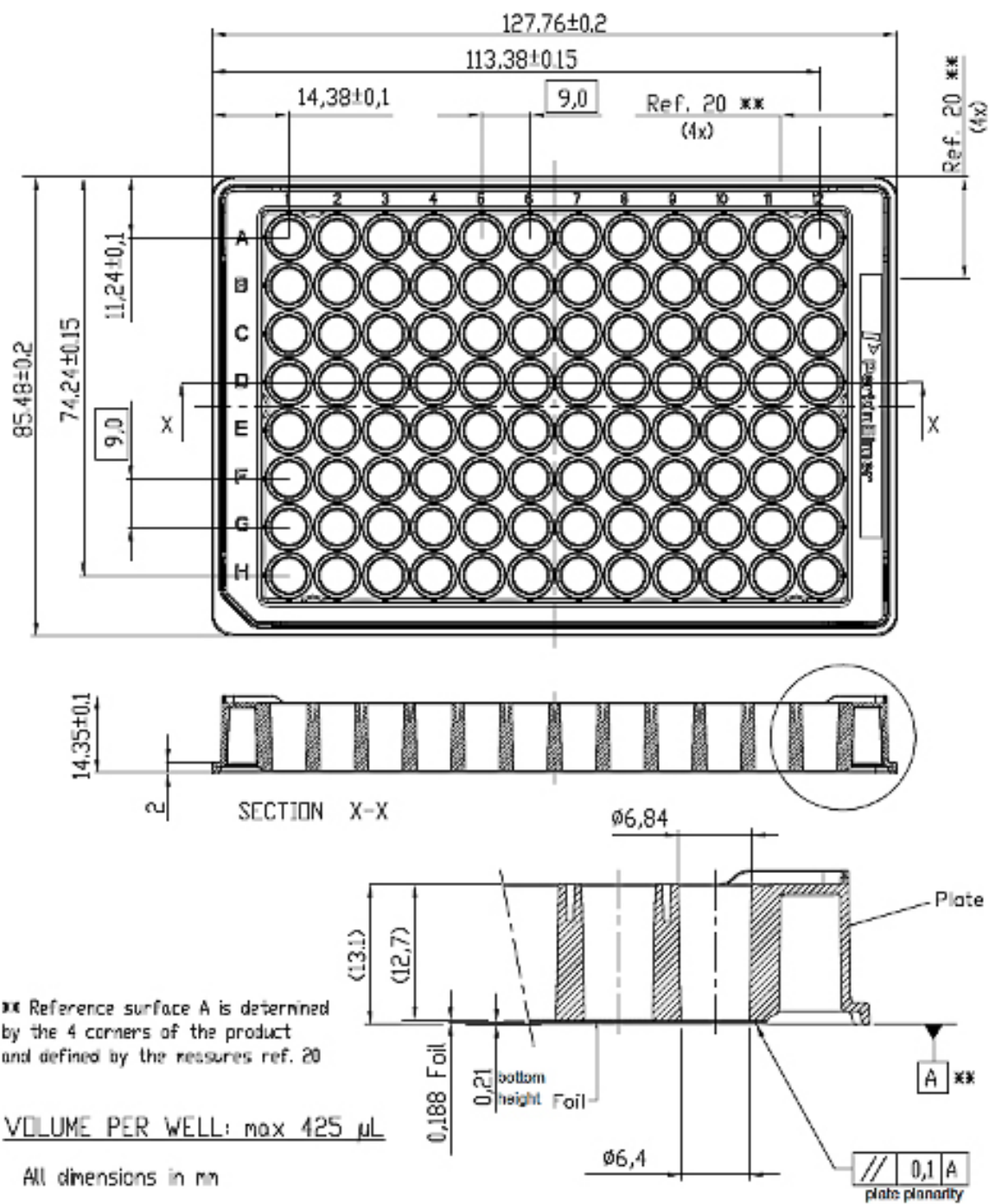
- Optimal clarity and fast autofocusing from excellent flatness of the plate bottom.
- Superior image quality from high optical quality of cyclic olefin imaging surface - black optically clear cyclic olefin microplate (Refractive index of 1.53).
- Better well access when using water immersion and high NA objectives with ultra-low plate bottom ($200 \pm 10 \mu\text{m}$).
- Reduced evaporation from new design, low-profile polystyrene lid.
- Less risk of damaging the imaging surface when stacking with unique corner spacers.
- Available with different coatings to suit your application:
 - Tissue-culture treated.
 - Fibronectin-coated to mimic natural cellular environment and enhance cell adherence.
 - Poly-D-lysine or collagen-coated to facilitate cell adherence and proliferation.
 - ULA-coated to reduce cell adherence and intended for studying 3D spheroids or suspension cells.



Product information

Product name	Part number	Packaging format	Storage temperature
PhenoPlate 96-well, black, optically clear flat-bottom, tissue-culture treated, lids, case of 50	6055300	50 individually pouched plates	RT
PhenoPlate 96-well, black, optically clear flat-bottom, tissue-culture treated, lids, case of 40	6055302	2 bags of 20 plates	RT
PhenoPlate 96-well, black, optically clear flat-bottom, tissue-culture treated, lids, case of 160	6055308	8 bags of 20 plates	RT
PhenoPlate 96-well, black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 40	6055500	2 bags of 20 plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 160	6055508	8 bags of 20 plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, collagen-coated, lids, case of 40	6055700	2 bags of 20 plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, collagen-coated, lids, case of 160	6055708	8 bags of 20 plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, fibronectin-coated, lids, case of 10	6055600	10 individually pouched plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, fibronectin-coated, lids, case of 40	6055602	2 bags of 20 plates	4 °C
PhenoPlate 96-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 10	6055800	10 individually pouched plates	RT
PhenoPlate 96-well, black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 40	6055802	2 bags of 20 plates	RT

Plate dimensions



PhenoPlate dimensions for Harmony® software microplate definition wizard

Plate Type Definition Wizard

Welcome

Define Template

Specify Plate Dimensions

Specify Well Dimensions

Describe Plate

Finish Wizard

Step 4 / 7

Please enter the dimensions of the plate type.

A: 14.38 mm C: 127.76 mm E: 113.38 mm
B: 11.24 mm D: 85.48 mm F: 74.24 mm

< Back Next >

Plate Type Definition Wizard

Welcome

Define Template

Specify Plate Dimensions

Specify Well Dimensions

Describe Plate

Finish Wizard

Step 5 / 7

Side:

G: 14.35 mm
H: 0.21 mm
I: 0.19 mm

Top:

J: 6.84 mm

Bottom:

N: 6.40 mm

< Back Next >

revvity