

Technical Data Sheet: PhenoPlate™ 96-wells

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.

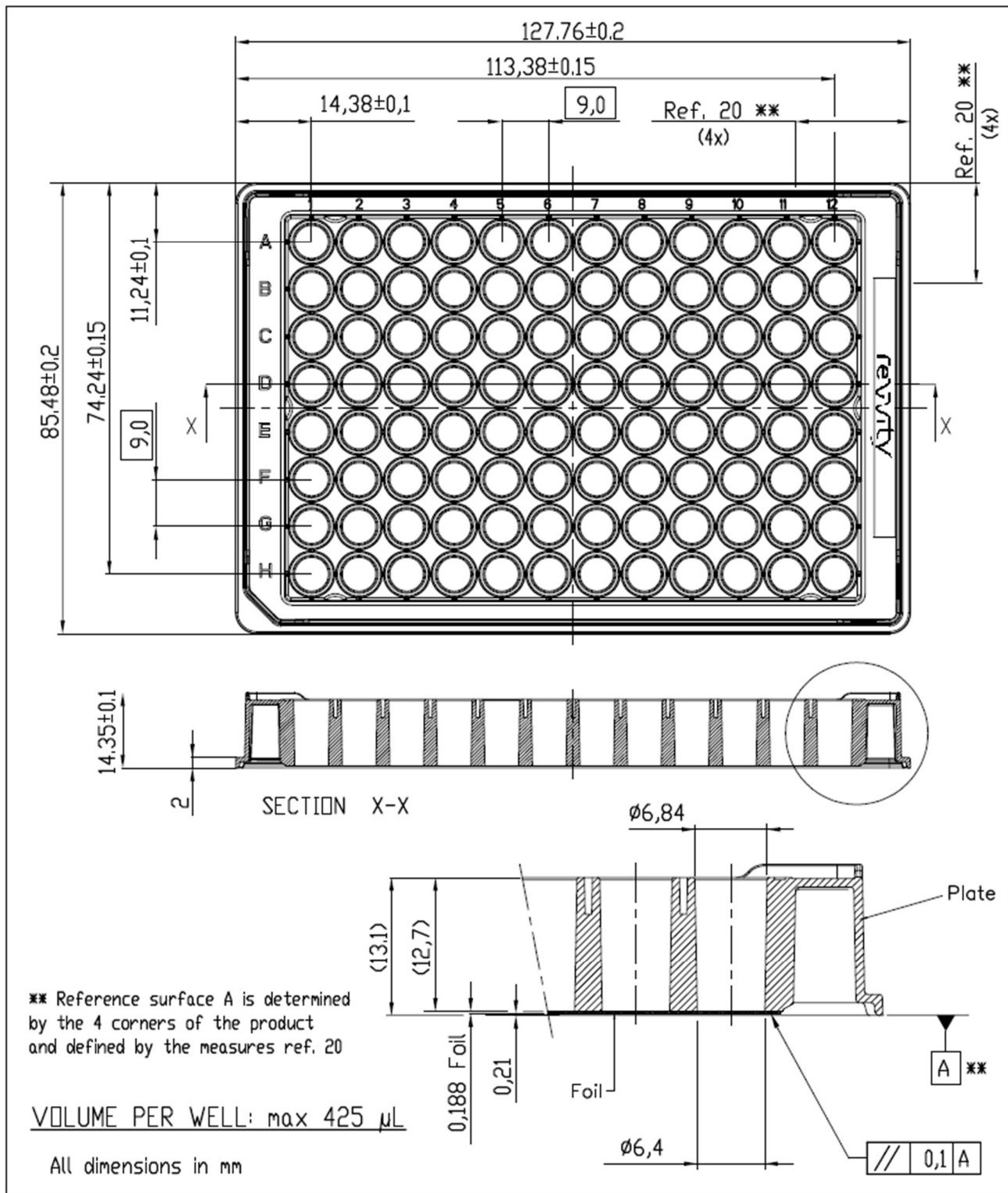
- Optimal clarity and fast autofocusing from excellent flatness of the plate bottom.
- Superior image quality from high optical quality of cyclic olefin imaging surface - black optically clear cyclic olefin microplate (Refractive index of 1.53).
- Better well access when using water immersion and high 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 (TC) treated.
 - Poly-D-lysine (PDL), collagen (COL) or Fibronectin (FN)-coated to facilitate cell adherence and proliferation.
 - Ultra Low Attachment (ULA)-coated to reduce cell adherence and intended for studying 3D spheroids or suspension cells.



Product Information:

Product Name	Packaging Format	Storage Temperature	Part Number
PhenoPlate 96-well, black, optically clear flat-bottom, TC-treated, lids, case of 50	50 individually pouched plates	RT	6055300
PhenoPlate 96-well, black, optically clear flat-bottom, TC-treated, lids, case of 40	2 bags of 20 plates	RT	6055302
PhenoPlate 96-well, black, optically clear flat-bottom, TC-treated, lids, case of 160	8 bags of 20 plates	RT	6055308
PhenoPlate 96-well, black, optically clear flat-bottom, PDL-coated, lids, case of 40	2 bags of 20 plates	4 °C	6055500
PhenoPlate 96-well, black, optically clear flat-bottom, PDL-coated, lids, case of 160	8 bags of 20 plates	4 °C	6055508
PhenoPlate 96-well, black, optically clear flat-bottom, COL-coated, lids, case of 40	2 bags of 20 plates	4 °C	6055700
PhenoPlate 96-well, black, optically clear flat-bottom, COL-coated, lids, case of 160	8 bags of 20 plates	4 °C	6055708
PhenoPlate 96-well, black, optically clear flat-bottom, ULA-coated, lids, case of 10	10 individually pouched plates	RT	6055800
PhenoPlate 96-well, black, optically clear flat-bottom, ULA-coated, lids, case of 40	2 bags of 20 plates	RT	6055802
PhenoPlate 96-well, black, optically clear flat-bottom, FN-coated, lids, case of 10	10 individually pouched plates	4 °C	6055600
PhenoPlate 96-well, black, optically clear flat-bottom, FN-coated, lids, case of 40	2 bags of 20 plates	4 °C	6055602

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: mm C: mm E: mm
 B: mm D: mm F: 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: mm
 H: mm
 I: mm

Top: Bottom:

J: mm N: mm

< Back Next >

