

# Technical Data Sheet:

## PhenoPlate™ 384-wells

### Introduction:

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	Part Number
PhenoPlate 384-well, black, optically clear flat-bottom, TC-treated, lids, case of 50	50 individually pouched plates	6057300
PhenoPlate 384-well, black, optically clear flat-bottom, TC-treated, lids, case of 40	2 bags of 20 plates	6057302
PhenoPlate 384-well, black, optically clear flat-bottom, TC-treated, lids, case of 160	8 bags of 20 plates	6057308
PhenoPlate 384-well, black, optically clear flat-bottom, TC-treated, non-irradiated, lids, case of 160	8 bags of 20 plates	6057328
PhenoPlate 384-well, black, optically clear flat-bottom, PDL-coated, lids, case of 40	2 bags of 20 plates	6057500
PhenoPlate 384-well, black, optically clear flat-bottom, PDL-coated, lids, case of 160	8 bags of 20 plates	6057508
PhenoPlate 384-well, black, optically clear flat-bottom, COL-coated, lids, case of 40	2 bags of 20 plates	6057700
PhenoPlate 384-well, black, optically clear flat-bottom, COL-coated, lids, case of 160	8 bags of 20 plates	6057708
PhenoPlate 384-well, black, optically clear flat-bottom, ULA-coated, lids, case of 10	10 individually pouched plates	6057800
PhenoPlate 384-well, black, optically clear flat-bottom, ULA-coated, lids, case of 40	2 bags of 20 plates	6057802
PhenoPlate 384-well, black, optically clear flat-bottom, FN-coated, lids, case of 10	10 individually pouched plates	6057600
PhenoPlate 384-well, black, optically clear flat-bottom, FN-coated, lids, case of 40	2 bags of 20 plates	6057602



# PhenoPlate 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 >

