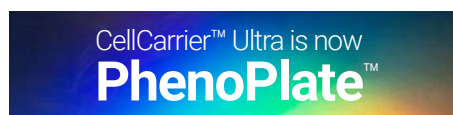


## PhenoPlates for High-Content Screening

### Better Microplates Means Better Results

To get better results from your high-content imaging system, you need better imaging microplates.



Drawing on many years of experience in high-content screening, PerkinElmer's expert team have developed and validated the PhenoPlate line of microplates (formerly known as CellCarrier™ Ultra plates). These microplates have been engineered to deliver both performance and superior images 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
- Better well access when using water immersion and high NA objectives with ultra-low plate bottom
- Reduced evaporation from new design, low-profile polystyrene lid
- Avoid damaging the imaging surface when stacking with unique corner spacers
- Available with different coatings to suit your application

PhenoPlates are part of a range of microplates for high content applications, and belong to our complete solution for high content imaging that includes the Operetta®CLS™ high-content analysis and Opera Phenix®Plus high-content screening systems, software, and automation.

96-well PhenoPlates	
Part Number	Description
6055300	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 50
6055302	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 40
6055308	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 160
6055500	Black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 40
6055508	Black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 160
6055700	Black, optically clear flat-bottom, collagen-coated, lids, case of 40
6055708	Black, optically clear flat-bottom, collagen-coated, lids, case of 160
6055800	Black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 10
6055802	Black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 40

384-well PhenoPlates	
Part Number	Description
6057300	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 50
6057302	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 40
6057308	Black, optically clear flat-bottom, tissue-culture treated, lids, case of 160
6057328	Black, optically clear flat-bottom, tissue-culture treated, non-irradiated, lids, case of 160
6057500	Black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 40
6057508	Black, optically clear flat-bottom, poly-D-lysine-coated, lids, case of 160
6057700	Black, optically clear flat-bottom, collagen-coated, lids, case of 40
6057708	Black, optically clear flat-bottom, collagen-coated, lids, case of 160
6057800	Black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 10
6057802	Black, optically clear flat-bottom, ultra-low-attachment-coated, lids, case of 40

PhenoPlate Lids	
Part Number	Description
6000031	Lid, clear, for PhenoPlates, sterile, case of 200

Please ask a representative for more information or for a free sample pack.  
Alternatively, visit [www.perkinelmer.com/phenoplates](http://www.perkinelmer.com/phenoplates)

PerkinElmer, Inc.  
940 Winter Street  
Waltham, MA 02451 USA  
P: (800) 762-4000 or  
(+1) 203-925-4602  
[www.perkinelmer.com](http://www.perkinelmer.com)



For a complete listing of our global offices, visit [www.perkinelmer.com/ContactUs](http://www.perkinelmer.com/ContactUs)

Copyright ©2021 PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.