

CellCarrier™ Ultra is now

PhenoPlate™

Superior Image Quality
For You and Your Science



Your trusted CellCarrier™ Ultra microplate has a new name - PhenoPlate™.

While the name may be changing, the manufacturing and ordering process, as well as our unwavering commitment to quality, is not. PhenoPlates will continue to deliver exactly the same high-performance imaging that you are accustomed to.

PhenoPlates are part our growing portfolio of products for high-content imaging, including PhenoVue™ cellular imaging reagents, that have been designed to work together to give you superior results.

All CellCarrier Ultra plates produced on or after September 15, 2021 will feature the name PhenoPlate™. Our website will be updated accordingly. www.perkinelmer.com/PhenoPlates

If you have any questions about the changes, please feel free to contact your local sales representative.

Yours Sincerely,

Fabienne Charrier-Savournin

Product Manager, Imaging Microplates



Part Number	Description
96-well PhenoPlates	
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	
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	
6000031	Lid, clear, for PhenoPlates, sterile, case of 200



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

Q & A

Q Why are you changing the name of CellCarrier Ultra plates?

A Unlike the names of other microplates in our portfolio, the CellCarrier Ultra name does not align with the primary application of the plates, which is cellular imaging and phenotypic screening – hence the new name PhenoPlate. The new name is also aligned with our recently-introduced PhenoVue Cellular Imaging Reagents.

Q Are part numbers and pricing changing?

A No. The changes are to the name only. You will order the plates using the same part numbers as before for easy reference.

Q Have the materials that the plates are made from or the plate coatings changed?

A No. The changes are to the name only. Our PhenoPlate microplates are made to exactly the same standards of manufacture as before, maintaining the same high-quality performance and functionality.

Q Are all CellCarrier plates being renamed?

A All CellCarrier Ultra plates are being renamed as PhenoPlates. CellCarrier Spheroid plates will retain their current names for the time being. Any future changes to this name will be communicated in a similar way.

Q I've received some plates with the new PhenoPlate name, but also some with the old CellCarrier Ultra name. How can I be sure a mistake was not made in preparing my order?

A It may take time for pre-existing stocks of some plate types to be used up. To avoid waste, any plate inventory featuring the CellCarrier Ultra name is not being discarded. There will be a transition period during which plates manufactured before the date of the re-brand will be sold and gradually superseded by plates printed with the new PhenoPlate name. You can verify that your order is correct by checking that you have received plates with the part numbers that you ordered. The part numbers are the same as before.



For a complete listing of our global offices, visit 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.