



CellCarrier Ultra Microplates for High-Content Screening

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 has developed and validated the CellCarrier™ Ultra line of microplates. 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

CellCarrier Ultra plates 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® High Content Imaging and Opera Phenix™ High Content Screening Systems, software, and automation.

Description	Part No.
CellCarrier Ultra 96 / Case of 50	6055300
CellCarrier Ultra 96 / Case of 40 (2x20)	6055302
CellCarrier-96 ultra / 8x20	6055308
PDL-Coated CellCarrier-96 ultra / 2x20	6055500
PDL-Coated CellCarrier-96 ultra / 8x20	6055508
COL-Coated CellCarrier-96 ultra / 2x20	6055700
COL-Coated CellCarrier-96 ultra / 8x20	6055708
CellCarrier-384 ultra / Case of 50 (50x1)	6057300
CellCarrier-384 ultra / Case of 160 (8x20)	6057308
PDL-Coated CellCarrier-384 ultra / Case of 40 (2x20)	6057500
PDL-Coated CellCarrier-384 ultra / Case of 160 (8x20)	6057508
COL-Coated CellCarrier-384 ultra / Case of 40 (2x20)	6057700
COL-Coated CellCarrier-384 ultra	6057708

Please ask a representative for more information or for a free sample pack. Alternatively, visit www.perkinelmer.com/microplates

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


PerkinElmer[®]
For the Better

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2014-2016, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.

011732D_01

PKI