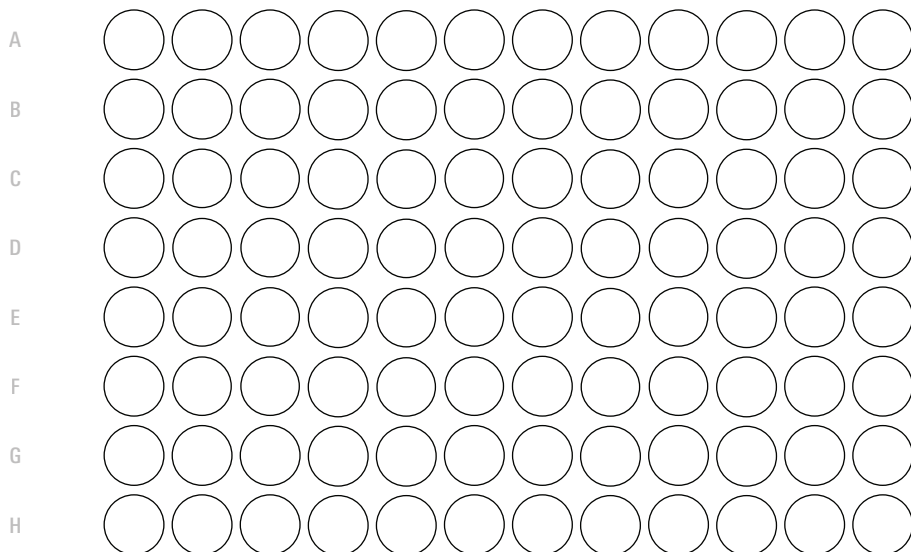
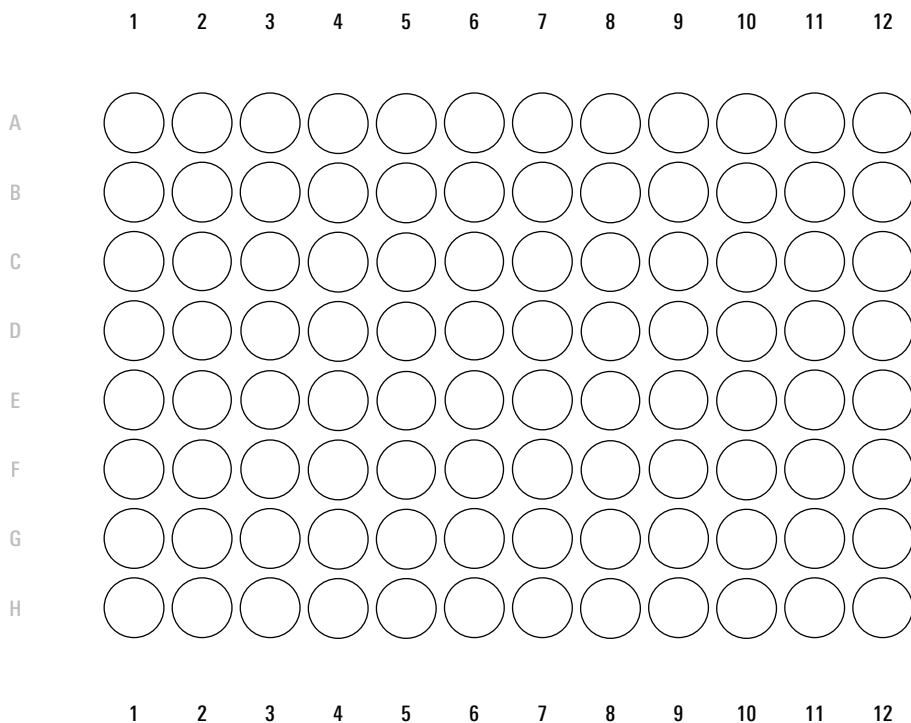


Blank Plate Loading Guides

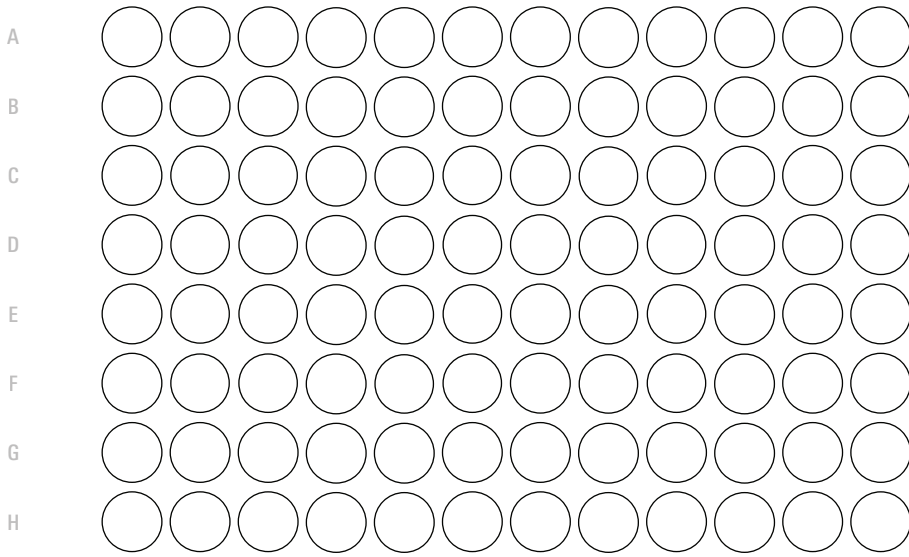
These blank plate layouts can be placed under a 96-well multiwell plate to make it easier to pipette the correct solutions into the correct wells. The loading guides below are intended to fit under Greiner Bio-One 96-Well μ CLEAR[®] Plates (LI-COR PN 926-19156, 8 pack; 926-19157, 32 pack). Label the rows and columns or color in the wells as appropriate for your experiment.

To easily create a custom plate layout, use the Experiment Designer workflow in Empiria Studio[®] Software. The Experiment Designer allows you to create a Plate Template that can be printed for use as a loading guide and used in Empiria Studio for quick multiwell plate analysis. See licor.com/empiria for more information.

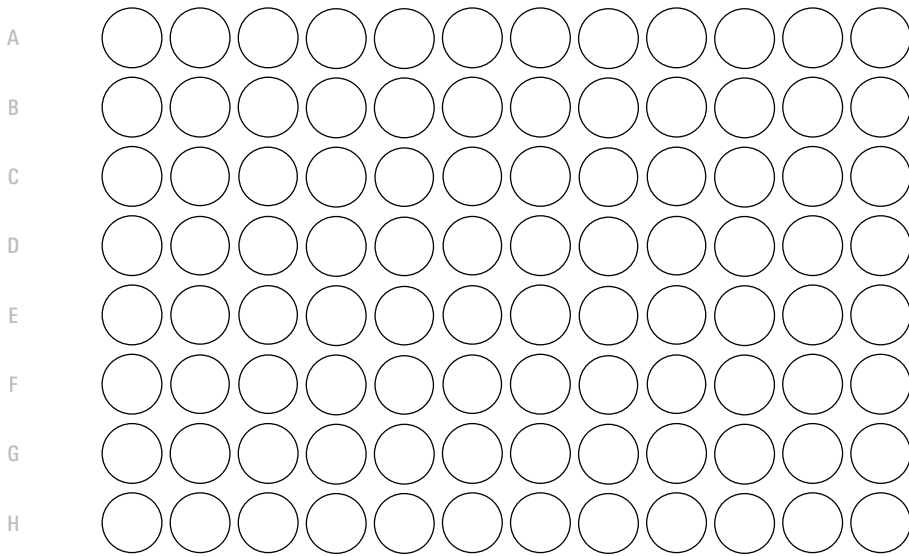
🖨 Printing this document? The correct print settings may vary depending on your printer. Look for an option called "Actual Size" for the best likelihood of printing the plates at a size that will fit under Greiner Bio-One 96-Well μ CLEAR[®] Plates.



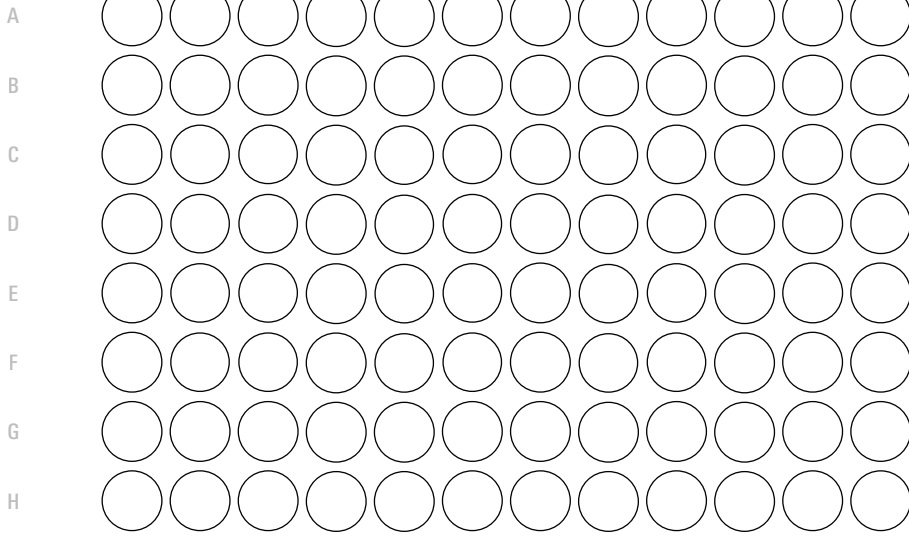
1 2 3 4 5 6 7 8 9 10 11 12



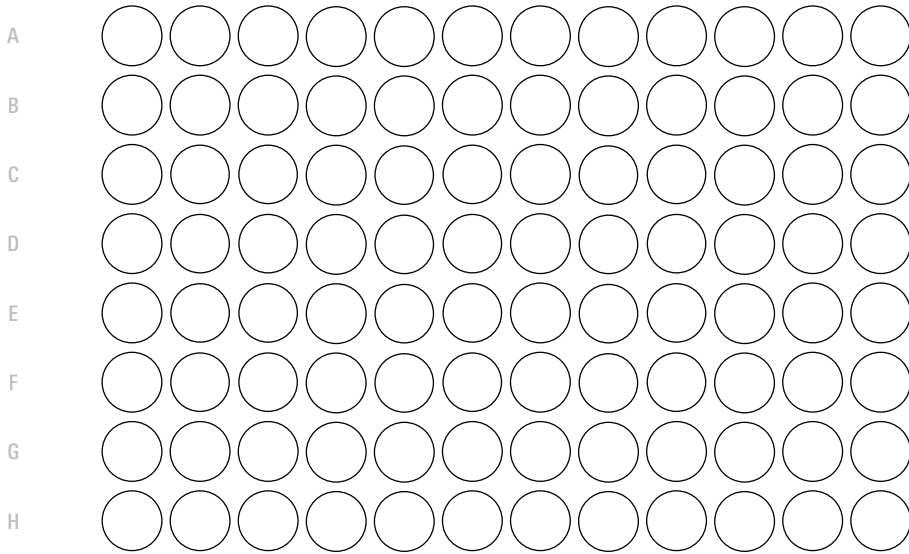
1 2 3 4 5 6 7 8 9 10 11 12



1 2 3 4 5 6 7 8 9 10 11 12



1 2 3 4 5 6 7 8 9 10 11 12





© 2021 LI-COR, Inc. LI-COR is a registered trademark of LI-COR, Inc. in the United States and other countries. All other trademarks belong to their respective owners.

LI-COR Biosciences

4647 Superior Street
Lincoln, NE 68504
Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com
licor.com/bio

Regional Offices

LI-COR Biosciences GmbH
Siemensstraße 25A
61352 Bad Homburg
Germany
Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Biosciences UK Ltd.
St. John's Innovation Centre
Cowley Road • Cambridge
CB4 0WS • United Kingdom
Phone: +44 (0) 1223 422104
bio-eu@licor.com