

Seeding Cells

There are three factors to consider when seeding cells:

1: Plates

For most adherent cells that stick to wells tightly (e.g. A431, HeLa, HEK293, CHO), we recommend regular tissue culture microplates with low auto-fluorescence, such as Nunc part no. 167008. For adherent cells that could detach from wells during In-Cell Western™ assay wash steps (e.g. NIH/3T3), we recommend Poly-D-lysine coated 96-well microplates, such as Sigma part no. Z38294-3.

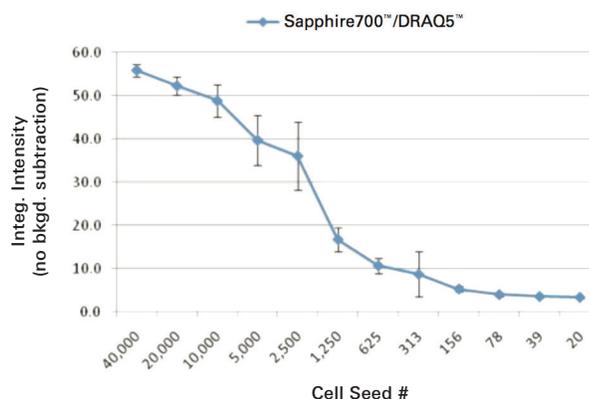
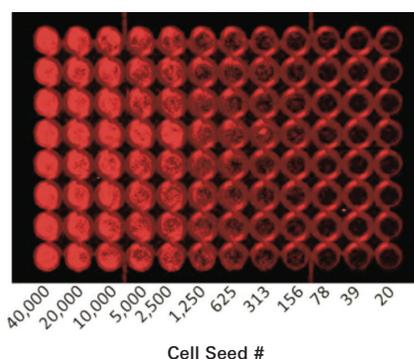
2: Cell seeding density

Typically, 15,000 to 40,000 cells are seeded per well. Two to three days are usually required for cells to reach the appropriate confluency, depending on growth rate. Seeding with low cell numbers is recommended if you plan to culture for several days before use. Plates seeded with higher cell numbers will be ready to use earlier.

3: Confluence

To obtain maximal fluorescent signals, complete or near complete confluency is recommended for cells that stick to wells tightly. For cells that adhere loosely to wells such as NIH/3T3, 70% confluency should be used. Please note that cell type and experimental conditions may affect the acceptable level of growth confluency.

The example below illustrates the importance of cell seeding density for A431 cells. Cell density is more important for some cell lines than others. In particular, cells that depend more on extracellular activity for proliferation (such as epithelial cells) are affected to a greater extent by initial growth conditions. As shown in the corresponding graph, cell growth is greatly inhibited when there are too few neighboring cells.



LI-COR Biosciences

Serving the United States and Canada.

Phone: +1-402-467-0700
Toll free: 800-645-4267
biosales@licor.com

LI-COR GmbH, Germany

Serving Europe, Africa, and the Middle East.

Phone: +49 (0) 6172 17 17 771
bio-eu@licor.com

LI-COR Ltd., United Kingdom

Serving Denmark, Finland, Iceland, Ireland, Norway, Sweden, and UK.

Phone: +44 (0) 1223 422104
bio-eu@licor.com

LI-COR Distributor Network

www.licor.com/bio/distributors

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.

For patent information, visit www.licor.com/patents.

©2017 LI-COR, Inc.

982-10597 08/17 Rev. A