

## AssayComplete™ Cell Plating Reagent

### Materials Provided

<b>Catalog Number:</b>	93-0563RXXX Series
<b>Quantity Supplied:</b>	100 mL (size A) and 500 mL (size B)
<b>Lot Number:</b>	See product label

### Description

Each DiscoverX cell line has been optimized for use with a specific AssayComplete Cell Plating Reagent to produce the highest possible assay performance. Each cell plating reagent was developed to maintain high cell viability and maximize ligand responsiveness for the duration of the cell-based assay while minimizing high basal activity by eliminating additional hormones and growth components that could impact the assay. These reagents are only to be used when preparing assay plates and are not intended for use during routine growth and passaging of DiscoverX cell lines.

### Product Information

<b>Storage Conditions:</b>	Store at -20°C. Thaw contents at room temperature and mix well by gently inverting the bottle prior to use. Once thawed, store at 4°C for up to 4 weeks. Avoid multiple freeze/thaw cycles.
<b>Shelf Life:</b>	See product label for expiration date
<b>Shipping Conditions:</b>	Frozen on dry ice (-70°C)
<b>Instructions on Use:</b>	Refer to cell line user manual for usage.

### Notice to Purchaser

AssayComplete Cell Plating Reagent has been optimized with DiscoverX cell lines. Occasionally, the reagent may be yellow or pink in color. This indicates a slight change in pH has occurred. It has been determined that using discolored media does not adversely impact cell growth or assay performance. Refer to the FAQ section for a more detailed explanation of this phenomenon.

For research use only. Not intended for use in diagnostic or therapeutic procedures.

### Quality Control Data

Each AssayComplete Cell Plating Reagent has been functionally tested with an appropriate control DiscoverX cell line. The EC<sub>50</sub> and signal-to-background (S:B) ratio was demonstrated to be comparable to the AssayComplete Cell Plating Reagent qualified lot.

This product includes high quality serum that has been tested for optimal performance. Testing includes heat inactivation, sterility, endotoxin ( $\leq 5$  EU/ml), pH ( $\geq 6.9$  to  $\leq 7.8$ ), osmolality ( $\geq 280$  to  $\leq 340$ ), total protein ( $\geq 3.0$  to  $\leq 5.0$ ), hemoglobin ( $\leq 10$  mg/dl), mycoplasma, virus, and bacteriophage levels determination.

### FAQ

#### [After the reagent was thawed, the media was discolored, is this normal?](#)

---

Variations in color of the media reagent can occur. These variations result from CO<sub>2</sub> vapor from the dry ice in the shipping package altering the concentrations of dissolved CO<sub>2</sub> in the media reagents. This can result in slight changes of pH that will induce a yellowing of the phenol red indicator dye present in the media.

DiscoverX has determined that use of media with a yellow color does not affect cell growth or assay performance. Exposure of discolored media to 5% CO<sub>2</sub> in a tissue culture incubator will re-balance the media to proper physiological pH which will normalize the color accordingly.

For additional information or Technical Support, see contact information below.