

## AssayComplete™ Cell Plating 39 Reagent Pack

### Materials Provided

<b>Catalog Number:</b>	93-0563R39 series
<b>Quantity Supplied:</b>	100 mL and 1 mL L-Glutamine (size A) 500 mL and 5 mL L-Glutamine (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 12 weeks or until the expiration date on the product label, whichever comes first. Do not refreeze. Once L-glutamine has been added, the medium is stable for 10 days at 4°C.
<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>	On the day of the assay, add fresh L-glutamine to the cell plating reagent at a final concentration of 2mM (e.g. for 100 mL of cell plating reagent, add 1 mL of 200 mM (100X) L-glutamine). Refer to cell line user manual for additional details.

### Notice to Purchaser

AssayComplete Cell Plating 39 Reagent Pack contents have been optimized for use with DiscoverX cell lines. Occasionally, the reagent may be yellow or pink in color, indicating a slight change in the pH. It has been determined that using discolored media does not adversely impact growth of cultured cell lines or assay performance. Refer to the FAQs section for a 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 are compared to that of the AssayComplete Cell Plating Reagent qualified lot.

### FAQs

#### 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.

We recommend incubating an aliquot of the discolored media in a 5% CO<sub>2</sub> tissue culture incubator until it regains its color, to re-balance the media to its proper pH prior to use in the assay.

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