

## AssayComplete™ Protein Dilution Buffer

### Materials Provided

<b>Catalog Number:</b>	92-0023 Series
<b>Quantity Supplied:</b>	1 X 50 mL (size S)*, 2 X 50 mL (size M) and 2 X 250 mL (size L) * Custom size, available upon request
<b>Lot Number:</b>	See product label

### Description

This ready-to-use AssayComplete Protein Dilution Buffer is formulated and optimized for DiscoverX cell lines and used for protein agonists to help in reconstituting them to the optimal stock concentration. In addition, this buffer is routinely used to dilute proteins and antibodies used in specific DiscoverX assays.

### Product Information

<b>Storage Conditions:</b>	Store at $-20^{\circ}\text{C}$ . Thaw contents at room temperature and mix well by gently inverting the bottle prior to use. Once thawed, store at $4^{\circ}\text{C}$ 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 ( $-70^{\circ}\text{C}$ )
<b>Instructions on Use:</b>	Refer to cell line user manual

### Notice to Purchaser

Each AssayComplete Protein Dilution Buffer has been optimized with DiscoverX cell lines.

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