

AssayComplete™ Cell Detachment Reagent

Materials Provided

Catalog Number:	92-0009
Quantity Supplied:	50 mL
Lot Number:	See product label

Description

Ready-to-use AssayComplete Cell Detachment Reagent is formulated and optimized to be used with DiscoverX cell lines to sustain intact cell membrane morphology upon detachment of cells from tissue culture plastic ware. It is used for routine cell passaging and preparation of assay plates.

Product Information

Storage Conditions:	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.
Shelf Life:	See product label for expiration date
Shipping Conditions:	Frozen on dry ice (-70°C)
Instructions on Use:	Refer to cell line user manual

Notice to Purchaser

AssayComplete Cell Detachment Reagent has been optimized with DiscoverX cell lines.

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

Quality Control Data

AssayComplete Cell Detachment Reagent is tested for sterility, pH, enzyme activity, and the ability to successfully dissociate cells from tissue culture plastic ware.

DiscoverX® name and the DiscoverX® logo, are a registered trademark of DiscoverX Corporation.
© 2016 DiscoverX Corporation. All rights reserved.