

AssayComplete™ Hygromycin B

Materials Provided

Catalog Number:	92-0029
Quantity Shipped:	1 X 20 mL
Stock Concentration:	50 mg/mL
Lot Number:	See product label

Description

Ready-to-use AssayComplete Hygromycin B is a sterile-filtered, selection antibiotic used for the selection and maintenance of eukaryotic cells transfected with the E. coli Hygromycin resistance gene. The effective killing concentration of this antibiotic will vary with the cell type, media, growth conditions, density, and the cell's metabolic rate and position in the cell cycle. For effective concentrations of Hygromycin B to be used in cell culture kits, please refer to cell line specific datasheet.

Product Information

Storage Conditions:	Store at -20°C. Protect from light.
Shelf Life:	See product label for expiration date. If stored properly, this product has data supporting a shelf life as indicated above. For optimal performance, avoid multiple freeze/thaw cycles.
Shipping Conditions:	Frozen on dry ice (-70°C).
Instructions on Use:	Refer to cell line user manual.

Related Products*

AssayComplete Cell Culture Kits (92-31XXG Series)
AssayComplete Puromycin (92-0028)
AssayComplete G418 (92-0030)

*Refer to cell line datasheet to determine catalog numbers for each related product required for the assay.

Notice to Purchaser

AssayComplete Hygromycin B has been optimized for use with DiscoverX cell-based assays.

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