

AssayComplete™ Puromycin

Materials Provided

Catalog Number:	92-0028
Quantity Shipped:	1 X 1 mL
Stock Concentration:	10 mg/mL
Lot Number:	See product label

Description

Ready-to-use AssayComplete Puromycin is sterile-filtered and belongs to the amino-nucleoside family of antibiotics and is isolated from *S. alboniger*. Puromycin is a broad spectrum antibiotic and antibacterial agent. It is active against Gram-positive microorganisms, less active against acid-fast bacilli, and weakly active against Gram-negative microorganisms. For effective concentrations of Puromycin 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 Hygromycin B (92-0029)
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 Puromycin has been optimized for use with DiscoverX cell-based assays.

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