

U-18666A

U 18666A is a small molecule inhibitor of cholesterol transport from endosomes/lysosomes to the endoplasmic reticulum in a variety of cell types, including macrophages and sympathetic neurons. Also weakly inhibits hedgehog signaling. Implicated in multiple diseases including Neimann-Pick disease type C (NPC), alzheimer's, cataracts, and atherosclerosis.

Materials Provided

Catalog Number:	92-1094
Quantity Supplied:	1 vial(s), 5 mg per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	424.06 g/mol
Source:	Synthetic
Purity:	≥ 98% by HPLC
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	≤ -20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	10 mM
Reconstitution Volume:	1.17 mL
Reconstitution Solvent:	DMSO

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com