

1-Oleoyl-LPA

1-Oleoyl lysophosphatidic acid is a lysophosphatidic acid (LPA) analog that potently induces calcium mobilization in A431 cells and growth stimulation in various cell lines. It is an agonist of LPA1 and LPA2, and inhibits neural stem cell (NSC) differentiation into neurons.

Materials Provided

Catalog Number:	92-1041
Quantity Supplied:	1 vial(s), 1 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:	436.9 g/mol
Source:	Synthetic
Purity:	≥ 95%
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:	22.9 mM
Reconstitution Volume:	Follow Handling Instructions
Reconstitution Solvent:	N/A

Handling Instructions:

Starting from the 22.9 mM ethanol stock, make a 2 mM intermediate stock in 100% ethanol. From the 2 mM intermediate stock, make a 500 μ M solution using PBS/0.1% BSA as the diluent (one part 2 mM ethanol stock : 3 parts PBS/0.1% BSA). Generate 3-fold serial dilutions from the 500 μ M solution using a diluent consisting of 75% PBS/0.1% BSA and 25% ethanol.

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