

LPI, L- α -Lysophosphatidylinositol

L- α -Lysophosphatidylinositol (LPI) is an endogenous phospholipid and endocannabinoid neurotransmitter. Novel ligand for the G protein-coupled receptor 55 (GPR55), promoter of RhoA-dependent calcium signaling, and NFAT activator in embryonic kidney 293 cells.

Materials Provided

Catalog Number:	92-1120
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:	621.4 g/mol
Source:	Soybean
Purity:	$\geq 98\%$
Endotoxin Level:	N/A
Formulation:	N/A
Storage:	$\leq -20^{\circ}\text{C}$ Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	1 mM
Reconstitution Volume:	1.609 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