

Sphingosine 1-phosphate

Sphingosine 1-phosphate is a signaling sphingolipid and bioactive lipid mediator that is responsible for mobilizing intracellular calcium stores, cell proliferation and motility, and decreasing cellular cAMP. Agonist for sphingosine 1-phosphate receptors, and activator of GPR3, GPR6 and GPR12. Also regulates effectors MAPK, PLC, adenylyl cyclase, myosin light chain phosphatase and FAK.

Materials Provided

Catalog Number:	92-1118
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:	379.5 g/mol
Source:	Synthetic
Purity:	≥ 98%
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:	2.5 mM
Reconstitution Volume:	1.054 mL
Reconstitution Solvent:	0.2 N NaOH

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