

FTY720 Phosphate

FTY 720 phosphate, or Fingolimod, is a sphingosine analog and immunomodulating drug currently approved for multiple sclerosis treatment. It is derived from the fungal metabolite myriocin (ISP-1), and acts as a potent modulator for sphingosine-1 phosphate receptors, encouraging recruitment of lymphocytes into the CNS during an immune response.

Materials Provided

Catalog Number:	92-1104
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:	387.5 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:	1 mM
Reconstitution Volume:	2.579 mL
Reconstitution Solvent:	Chloroform

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