FTY-720 . HCI

S1P receptor ligand

Novel immune modulator that alters migration and homing of lymphocytes. Acts as a potent agonist of sphingosine 1-phosphate receptors (S1P) after phosphorylation by sphingosine kinase. Also inhibits cytosolic phospholipase A_2 independently of sphingosine 1-phosphate receptors.

Citations: 11

View Online »

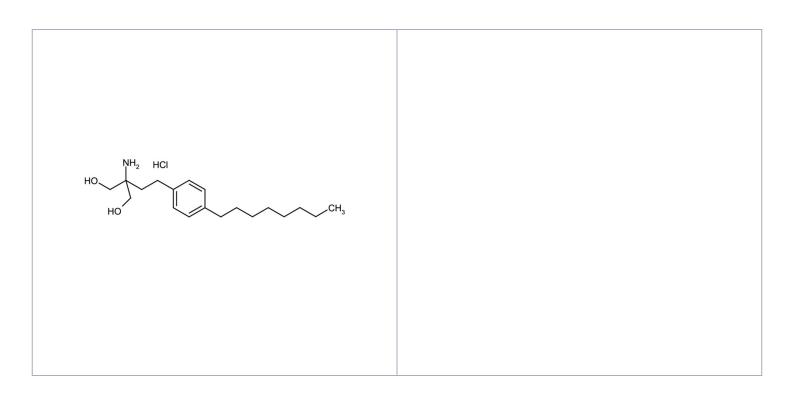
Ordering Information

Order Online »

BML-SL233-0005	5mg
BML-SL233-0025	25mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up

to 3 months at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-Amino-2-[2-(4-octylphenyl)ethyl]-1,3-propanediol . HCl,

Fingolimod . HCl

Appearance White solid.

CAS 162359-56-0

Couple Target Phospholipase, PLA, S1P receptor

Couple Type Ligand

Formula $C_{19}H_{33}NO_2$. HCI

Identity Determined by NMR.

MW 307.4 . 36.5

Purity ≥98% (HPLC)

Soluble in DMSO (10mg/ml), 100% ethanol or dimethyl

formamide.

Last modified: May 29, 2024



usa@enzolifesciences.com