

FTY-720 . HCl

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 A₂ independently of sphingosine 1-phosphate receptors.

Citations: 11

[View Online »](#)

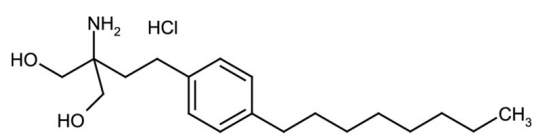
Ordering Information

[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	$C_{19}H_{33}NO_2 \cdot HCl$
Identity	Determined by NMR.
MW	307.4 . 36.5
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/ml), 100% ethanol or dimethyl formamide.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)