

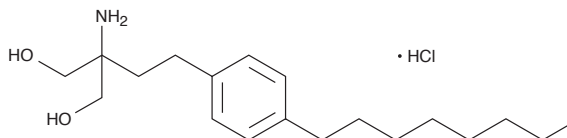
Product Information



FTY720

Catalog No. 10006292

CAS Registry No.: 162359-56-0
Formal Name: 2-amino-2-[2-(4-octylphenyl)ethyl]-1,3-propanediol, hydrochloride
MF: C₁₉H₃₃NO₂ • HCl
FW: 343.9
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that FTY720 be stored as supplied at -20°C. It will be stable for at least two years.

FTY720 is supplied as a crystalline solid. A stock solution may be made by dissolving the FTY720 in an organic solvent purged with an inert gas. FTY720 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of FTY720 in these solvents is at least 20 mg/ml.

FTY720 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FTY720 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. FTY720 has a solubility of 0.2 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

FTY720 is a derivative of ISP-1 (myriocin), a fungal metabolite of the Chinese herb *Iscaria sinclairii* as well as a structural analog of sphingosine. It is a novel immune modulator that prolongs allograft transplant survival in numerous models by inhibiting lymphocyte emigration from lymphoid organs.¹ FTY720 is phosphorylated by sphingosine kinase, which then acts as a potent agonist at four of the sphingosine-1-phosphate (S1P) receptors (S1P₁, S1P₃, S1P₄, and S1P₅).² Down-regulation of S1P₁ receptors on T and B lymphocytes by FTY720 results in defective egress of these cells from spleen, lymph nodes, and Peyer's patch.³ FTY720 also enhances the activity of the sphingosine transporter Abcb1 and the leukotriene C₄ transporter Abcc1 and inhibits cytosolic phospholipase A₂ activity.^{4,5}

References

1. Brinkmann, V., Pinschewer, D.D., Feng, L., *et al.* FTY720: Altered lymphocyte traffic results in allograft protection. *Transplantation* **72**, 764-769 (2001).
2. Brinkmann, V., Davis, M.D., Heise, C.E., *et al.* The immune modulator FTY720 targets sphingosine 1-phosphate receptors. *J. Biol. Chem.* **277**(24), 21453-21457 (2002).
3. Matloubian, M., Lo, C.G., Cinamon, G., *et al.* Lymphocyte egress from thymus and peripheral lymphoid organs is dependent on S1P receptor 1. *Nature* **427**, 355-360 (2004).
4. Honig, S.M., Fu, S., Mao, X., *et al.* FTY720 stimulates multidrug transporter- and cysteinyl leukotriene-dependent T cell chemotaxis to lymph nodes. *J. Clin. Invest.* **111**(5), 627-637 (2003).
5. Payne, S.G., Oskeritzian, C.A., Griffiths, R., *et al.* The immunosuppressant drug FTY720 inhibits cytosolic phospholipase A₂ independently of sphingosine-1-phosphate receptors. *Blood* **109**(3), 1077-1085 (2007).

Related Products

Sphingosine-1-phosphate - Cat. No. 62570 • Fumonisin B₁ - Cat. No. 62580 • S1P₁ Polyclonal Antibody - Cat. No. 10005228 • S1P₃ Polyclonal Antibody - Cat. No. 10006373 • FTY720 Phosphoric Acid Ester - Cat. No. 10008639

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 04/26/2007