

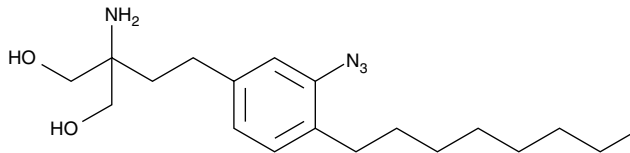
Product Information



azido-FTY720

Item No. 10008612

CAS Registry No.: 881914-35-8
Formal Name: 2-amino-2-[2-(4-(1-azido-2-octylphenyl))ethyl]-1,3-propanediol
MF: C₁₉H₃₂N₄O₂
FW: 348.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A solution in ethanol
UV/Vis.: λ_{max}: 213, 252, 284 nm



Laboratory Procedures

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

azido-FTY720 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of azido-FTY720 in these solvents is approximately 20 mg/ml.

azido-FTY720 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of azido-FTY720 should be diluted with the aqueous buffer of choice. azido-FTY720 has a solubility of approximately 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 potent immunosuppressive agent 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 azido-FTY720 is a highly photoreactive analog of FTY720 that can be used to identify receptor binding sites for this ligand.⁴

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. Sun, C. and Bittman, R. A photoreactive analogue of the immunosuppressant FTY720. *J. Org. Chem.* **71**, 2200-2202 (2006).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008612

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

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 Safety Data Sheet, which has been sent *via* email to 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, 12/12/2013

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