

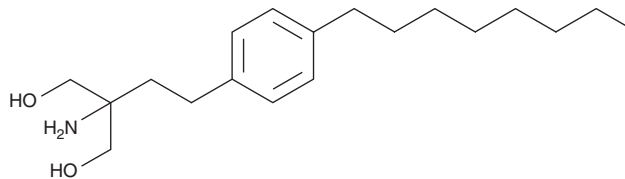
PRODUCT INFORMATION



Fingolimod

Item No. 11975

CAS Registry No.: 162359-55-9
Formal Name: 2-amino-2-[2-(4-octylphenyl)ethyl]-1,3-propanediol
Synonym: FTY720
MF: C₁₉H₃₃NO₂
FW: 307.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Fingolimod is supplied as a crystalline solid. A stock solution may be made by dissolving the fingolimod in the solvent of choice, which should be purged with an inert gas. Fingolimod is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of fingolimod in these solvents is approximately 5, 10, and 20 mg/ml.

Description

Fingolimod is an immune modulator, a prodrug form of the sphingosine-1-phosphate (S1P) receptor agonist FTY720 phosphate (Item No. 10008639), and a derivative of myriocin (Item No. 63150).¹ It selectively inhibits calcium-dependent cytosolic phospholipase A₂ (cPLA₂) over calcium-independent PLA₂ (iPLA₂) and secretory PLA₂ (sPLA₂) (Type V) when used at a concentration of 800 pmol and inhibits S1P lyase when used at concentrations ranging from 0.3 to 30 μM.^{2,3} Fingolimod (1 mg/kg) decreases the number of circulating lymphocytes in mice.³ It increases skin allograft survival in rats when administered at doses ranging from 0.3 to 3 mg/kg.⁴ Fingolimod (0.3 mg/kg) prevents disease development in a rat model of experimental autoimmune encephalomyelitis (EAE) induced by immunization with an emulsion of bovine spinal cord in complete Freund's adjuvant.¹ Formulations containing fingolimod have been used in the treatment of multiple sclerosis.

References

1. 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).
2. Payne, S.G., Oskeritizian, 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).
3. Bandhuvula, P., Tam, Y.Y., Oskouian, B., *et al.* The immune modulator FTY720 inhibits sphingosine-1-phosphate lyase activity. *J. Biol. Chem.* **280**(40), 33697-33700 (2005).
4. Shimizu, H., Takahashi, M., Kaneko, T., *et al.* KRP-203, a novel synthetic immunosuppressant, prolongs graft survival and attenuates chronic rejection in rat skin and heart allografts. *Circulation* **111**(2), 222-229 (2005).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/17/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM