

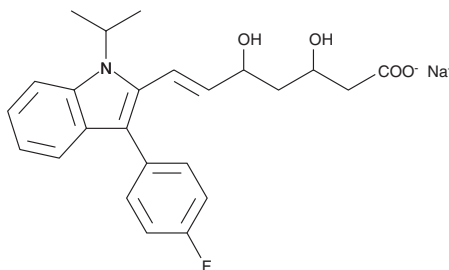
Product Information



Fluvastatin (sodium salt)

Catalog No. 10010337

CAS Registry No.: 93957-55-2
Formal Name: 7-[3-(4-fluorophenyl)-1-(1-methylethyl)-1H-indol-2-yl]-3,5-dihydroxy-6-heptenoic acid, sodium salt
MF: C₂₄H₂₅FNO₄ • Na
FW: 433.5
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that fluvastatin (sodium salt) be stored as supplied at -20°C. It should be stable for at least two years.

Fluvastatin (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the fluvastatin (sodium salt) in an organic solvent purged with an inert gas. Fluvastatin (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fluvastatin (sodium salt) in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. The solubility of fluvastatin (sodium salt) in PBS, pH 7.2, is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Hydroxymethylglutaryl-coenzyme A (HMG-CoA) reductase is the rate-limiting enzyme in the cholesterol biosynthetic pathway and the target of the 'statin' class of cholesterol-lowering drugs.¹ Fluvastatin acts as a competitive inhibitor of HMG-CoA reductase with respect to binding of the substrate HMG-CoA (K_i = 0.3 nM) but not with respect to binding of NADPH.^{2,3} It is marketed under the trade name Lescol™ for the treatment of hypercholesterolemia in the prevention of cardiovascular disease. When fluvastatin was administered in a clinical trial to patients undergoing percutaneous coronary intervention, LDL cholesterol levels were reduced by 27% after six weeks of treatment with a dose of 40 mg/kg twice a day compared to placebo.⁴

References

1. Tobert, J.A. Lovastatin and beyond: The history of the HMG-CoA reductase inhibitors. *Nature Reviews Drug Discovery* **2**, 517-526 (2003).
2. Istvan, E.S. and Deisenhofer, J. Structural mechanism for statin inhibition of HMG-CoA reductase. *Science* **292**, 1160-1164 (2001).
3. Corsini, A., Maggi, F.M., and Catapano, A.L. Pharmacology of competitive inhibitors of HMG-CoA reductase. *Pharmacol. Res.* **31(1)**, 9-27 (1995).
4. Serruys, P.W.J.C., de Feyter, P., Macaya, C., et al. Fluvastatin for prevention of cardiac events following successful first percutaneous coronary intervention: A randomized controlled trial. *JAMA* **287(24)**, 3215-3222 (2002).

Related Products

Berberine - Cat. No. 10006427 • PCSK9 (human) Polyclonal Antibody - Cat. No. 10007185 • LDL Receptor Polyclonal Antibody - Cat. No. 10007665 • PCSK9 (murine) Polyclonal Antibody - Cat. No. 10008811 • SREBP-2 Cell-Based Translocation Assay Kit - Cat. No. 10009239 • PCSK9 Western Ready Control - Cat. No. 10009567 • Lovastatin - Cat. No. 10010338 • Lovastatin (sodium salt) - Cat. No. 10010339 • Mevastatin - Cat. No. 10010340 • Mevastatin (sodium salt) - Cat. No. 10010341 • Pravastatin - Cat. No. 10010342 • Pravastatin (sodium salt) - Cat. No. 10010343 • Simvastatin - Cat. No. 10010344 • Simvastatin (sodium salt) - Cat. No. 10010345

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

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 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, 05/20/2010

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