

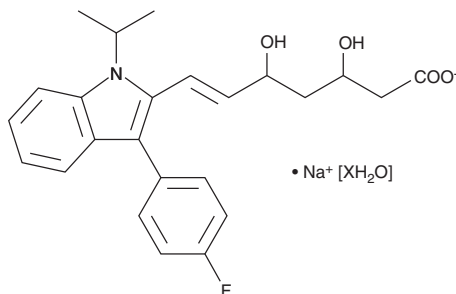
PRODUCT INFORMATION



Fluvastatin (sodium salt hydrate)

Item No. 10010337

Formal Name: 7-[3-(4-fluorophenyl)-1-(1-methylethyl)-1H-indol-2-yl]-3,5-dihydroxy-6-heptenoic acid, sodium salt, hydrate
MF: C₂₄H₂₅FNO₄ • Na [XH₂O]
FW: 410.5
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 305 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Fluvastatin (sodium salt hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the fluvastatin (sodium salt hydrate) in the solvent of choice. Fluvastatin (sodium salt hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of fluvastatin (sodium salt hydrate) 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 hydrate) in PBS, pH 7.2, is approximately 0.2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Fluvastatin is an inhibitor of HMG-CoA reductase (K_i = 0.3 nM for the rat enzyme).^{1,2} It also inhibits the human cytochrome P450 (CYP) isoform CYP2C9 (IC₅₀ = 100 nM).³ Fluvastatin inhibits oxidized LDL-induced ferroptosis and reverses oxidized LDL-induced decreases in glutathione peroxidase 4 (GPX4) and system X_c⁻ cystine-glutamate antiporter levels in human umbilical vein endothelial cells (HUVECs).⁶ *In vivo*, fluvastatin (2 mg/kg per day) decreases serum cholesterol, triglyceride, and phospholipid levels, the formation of thiobarbituric acid-reactive substances (TBARS), and vascular angiotensin-converting enzyme (ACE) activity in rabbits fed a high-cholesterol diet.⁴ It increases survival in a mouse model of myocardial infarction when administered at a dose of 10 mg/kg per day.⁵ Formulations containing fluvastatin have been used in the treatment of hypercholesterolemia and the prevention of cardiovascular disease.

References

1. Istvan, E.S. and Deisenhofer, J. *Science* **292**, 1160-1164 (2001).
2. Corsini, A., Maggi, F.M., and Catapano, A.L. *Pharmacol. Res.* **31(1)**, 9-27 (1995).
3. Transon, C., Leemann, T., and Dayer, P. *Eur. J. Clin. Pharmacol.* **50(3)**, 209-215 (1996).
4. Mitani, H., Bando, T., Ishikawa, J., et al. *Br. J. Pharmacol.* **119(6)**, 1269-1275 (1996).
5. Hayashidani, S., Tsutsui, H., Shiomi, T., et al. *Circulation* **105(7)**, 868-873 (2002).
6. Li, Q., Liu, C., Deng, L., et al. *Exp. Ther. Med.* **22(5)**, 1275 (2021).

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, 10/05/2021

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