

PRODUCT INFORMATION

3 α -hydroxy Pravastatin (sodium salt)

Item No. 35748

CAS Registry No.: 81093-43-8
Formal Name: (β R, δ R,1S,2R,3S,8S,8aR)-1,2,3,7,8,8a-hexahydro- β , δ ,3-trihydroxy-2-methyl-8-[(2S)-2-methyl-1-oxobutoxy]-1-naphthaleneheptanoic acid, monosodium salt

Synonyms: 3 α -Isopravastatin, R-416

MF: C₂₃H₃₆O₇ • Na

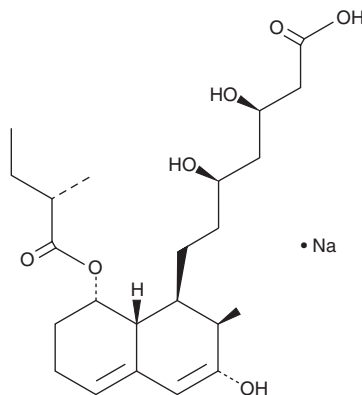
FW: 447.5

Purity: $\geq$ 95%

Supplied as: A solid

Storage: -20°C

Stability: $\geq$ 4 years



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

Laboratory Procedures

3 α -hydroxy Pravastatin (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the 3 α -hydroxy pravastatin (sodium salt) in the solvent of choice, which should be purged with an inert gas. 3 α -hydroxy Pravastatin (sodium salt) is slightly soluble in methanol.

3 α -hydroxy Pravastatin (sodium salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

3 α -hydroxy Pravastatin is a metabolite of the HMG-CoA reductase inhibitor pravastatin (Item No. 10010343).¹

Reference

1. Kawabata, K., Samata, N., and Urasaki, Y. Quantitative determination of pravastatin and R-416, its main metabolite in human plasma, by liquid chromatography-tandem mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **816**(1-2), 73-79 (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, 05/04/2023

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