

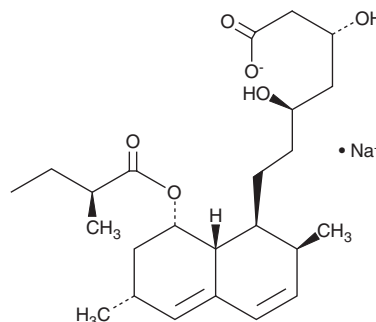
PRODUCT INFORMATION



Lovastatin Hydroxy Acid (sodium salt)

Item No. 10010339

CAS Registry No.: 75225-50-2
Formal Name: 2-methyl-1S,2,3R,7S,8S,8aR-hexahydro-3,7-dimethyl-8-[2-[(2R,4R)-tetrahydro-4-hydroxy-6-oxo-2H-pyran-2-yl]ethyl]-1-naphthalenyl ester, butanoic acid, monosodium salt
MF: C₂₄H₃₇O₆ • Na
FW: 444.5
Purity: ≥95%
UV/Vis.: λ_{max}: 231, 238, 246 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

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

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that 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

Lovastatin hydroxy acid is a potent and competitive inhibitor of HMG-CoA reductase ($K_i = 0.6 \text{ nM}$).¹ It is the more potent, open ring, hydroxy acid form of its prodrug, lovastatin (Item No. 10010338).²

References

1. Alberts, A.W., Chen, J., Kuron, G., *et al.* Mevinolin: A highly potent competitive inhibitor of hydroxymethylglutaryl-coenzyme A reductase and a cholesterol-lowering agent. *Proc. Natl. Acad. Sci. USA* **77**(7), 3957-3961 (1980).
2. Duggan, D.E., Chen, I.W., Bayne, W.F., *et al.* The physiological disposition of lovastatin. *Drug Metab. Dispos.* **17**(2), 166-173 (1989).

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, 09/29/2021

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