

# 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:** C<sub>24</sub>H<sub>37</sub>O<sub>6</sub> • Na

**FW:** 444.5

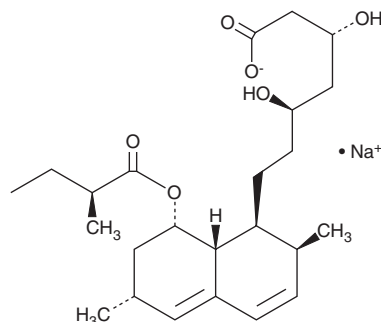
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 231, 238, 246 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

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 the solvent of choice, which should be 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.

### Description

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

### 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, 12/14/2022

#### 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