

# PRODUCT INFORMATION



## Lovastatin-d<sub>9</sub> Item No. 26461

**Formal Name:** 2S-(methyl-d<sub>3</sub>)-butanoic acid-2,3,3,4,4,4-d<sub>6</sub>, 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

**MF:** C<sub>24</sub>H<sub>27</sub>D<sub>9</sub>O<sub>5</sub>

**FW:** 413.6

**Chemical Purity:** ≥95% (Lovastatin; mixture of diastereomers)

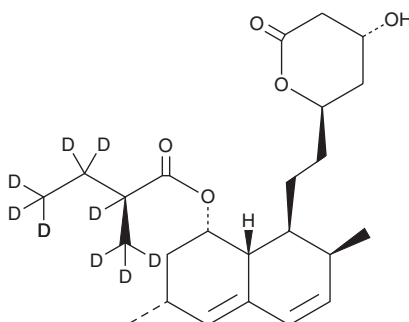
**Deuterium Incorporation:** ≥99% deuterated forms (d<sub>1</sub>-d<sub>9</sub>); ≤1% d<sub>0</sub>

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Synthetic



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

### Laboratory Procedures

Lovastatin-d<sub>9</sub> is intended for use as an internal standard for the quantification of lovastatin (Item No. 10010338) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Lovastatin-d<sub>9</sub> is supplied as a solid. A stock solution may be made by dissolving the lovastatin-d<sub>9</sub> in the solvent of choice, which should be purged with an inert gas. Lovastatin-d<sub>9</sub> is slightly soluble in chloroform and methanol.

### Description

Lovastatin-d<sub>9</sub> is intended for use as an internal standard for the quantification of lovastatin (Item No. 10010338) by GC- or LC-MS. Lovastatin is a fungal metabolite that has been found in *A. terreus* and an inhibitor of HMG-CoA reductase ( $K_i = 1.4$  nM).<sup>1,2</sup> It is also a prodrug form of the HMG-CoA reductase inhibitor lovastatin hydroxy acid (Item No. 10010339).<sup>2</sup> Lovastatin (8 mg/kg per day) reduces plasma cholesterol levels in dogs. It suppresses TNF-induced NF- $\kappa$ B activation ( $IC_{50} = \sim 15$   $\mu$ M) and potentiates apoptosis in human myeloid leukemia cells.<sup>3</sup> Lovastatin also increases cellular lipid peroxidation and decreases glutathione peroxidase 4 (GPX4) levels in cancer cells.<sup>4</sup> Formulations containing lovastatin have been used in the treatment of hypercholesterolemia.

### References

1. Endo, A. The discovery and development of HMG-CoA reductase inhibitors. *J. Lipid Res.* **33(11)**, 1569-1582 (1992).
2. 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).
3. Ahn, K.S., Sethi, G., and Aggarwal, B.B. Reversal of chemoresistance and enhancement of apoptosis by statins through down-regulation of the NF- $\kappa$ B pathway. *Biochem. Pharmacol.* **75(4)**, 907-913 (2008).
4. Viswanathan, V.S., Ryan, M.J., Dhruv, H.D., et al. Dependency of a therapy-resistant state of cancer cells on a lipid peroxidase pathway. *Nature* **547(7664)**, 453-457 (2017).

#### 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, 12/07/2022

#### 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