

Mevastatin

HMG-CoA reductase inhibitor

Mevastatin inhibits isoprenoid biosynthesis by inhibition of HMG-CoA reductase (K_i for acid form is 1 nM) and therefore blocks protein isoprenylation and reduces plasma cholesterol levels in humans. It causes cells to arrest early in the G1 phase. Mevastatin is a close structural analog of lovastatin and both agents have the same biochemical and pharmacological activities. Mevastatin is inactive in cell-free assays. In cells however, it is hydrolyzed to the active free acid form by intracellular esterases.

Citations: 8

[View Online »](#)

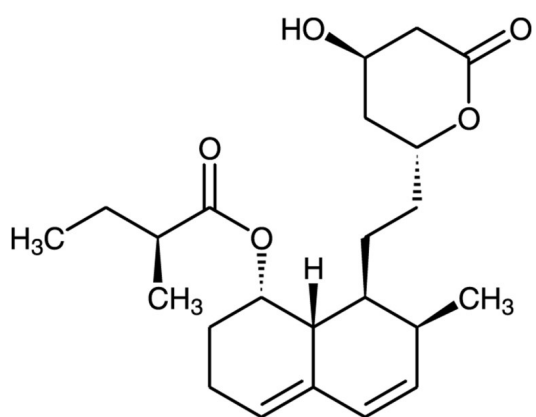
Ordering Information

[Order Online »](#)

BML-G233-0010	10mg
BML-G233-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Compactin

Appearance White to off-white solid.

CAS 73573-88-3

Couple Target HMG-CoA reductase

Couple Type Inhibitor

Formula $C_{23}H_{34}O_5$

Identity Determined by ¹H-NMR.

MW 390.5

Purity ≥95% (HPLC)

Solubility Soluble in DMSO (20mg/ml) or 100% ethanol (25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)