

Lovastatin

HMG-CoA reductase inhibitor

Inhibits cholesterol/isoprenoid biosynthesis by inhibition of HMG-CoA reductase (K_i for acid form is 1nM) and blocks protein isoprenylation and reduces plasma cholesterol levels in humans. Causes cells to arrest early in the G1 phase. Induces apoptosis in cancer cells.

Citations: 13

[View Online »](#)

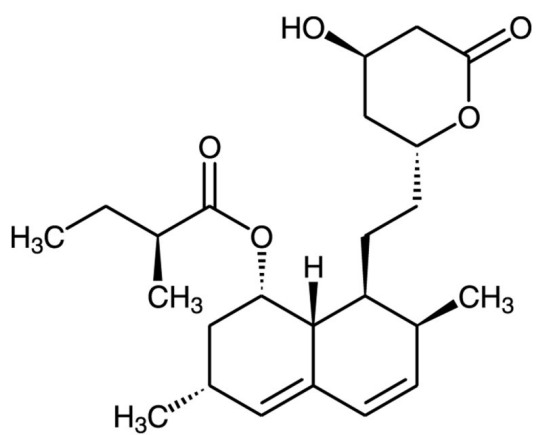
Ordering Information

[Order Online »](#)

BML-G226-0010	10mg
BML-G226-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Mevinolin, 6- α -Methylcompactin
Appearance	White solid.
CAS	75330-75-5
Couple Target	HMG-CoA reductase
Couple Type	Inhibitor
Formula	$C_{24}H_{36}O_5$
MI	14: 5586
MW	404.6
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>35mg/ml); insoluble in water.



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)