

Lovastatin . sodium salt

Active in whole cells as well as cell-free assays

Active carboxylate form of Lovastatin (Prod. No. BML-G226). Active in whole cells as well as cell-free assays. Inhibits cholesterol/ isoprenoid biosynthesis by inhibition of HMG-CoA reductase (K_i for acid form is 1 nM) and blocks protein isoprenylation and reduces plasma cholesterol levels in humans. Causes cells to arrest early in the G1 phase.

Citations: 4

[View Online »](#)

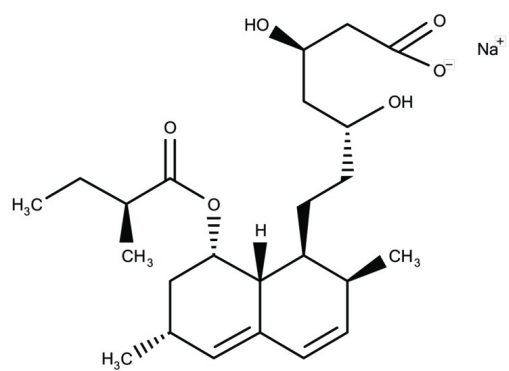
Ordering Information

[Order Online »](#)

BML-G245-0005	5mg
---------------	-----

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 . sodium salt, 6- α -Methylcompactin . sodium salt
Appearance	White solid.
CAS	75225-50-2
Couple Target	HMG-CoA reductase
Couple Type	Inhibitor
Formula	$C_{24}H_{37}O_6 \cdot Na$
MI	14: 5586
MW	444.5
Purity	$\geq 97\%$ (HPLC)
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol (10mg/ml).



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-
be@enzolifesciences.com](mailto:info-be@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)