

Simvastatin

HMG-CoA reductase inhibitor

HMG-CoA reductase inhibitor. Synthetic analog of Lovastatin (Prod. No. BML-G226). Inhibits the biosynthesis of isoprenoids and prevents small G protein activation. Displays lipid lowering and antihypertensive effects. Displays antiinflammatory, immunomodulatory effects as well as protective effects in experimental traumatic brain injury. Activates Rac1-dependent signalling pathway.

Citations: 15

[View Online »](#)

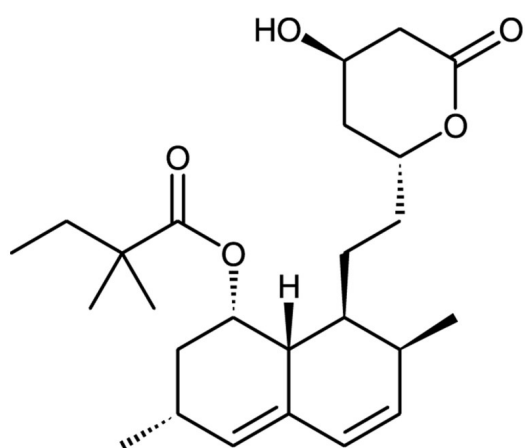
Ordering Information

[Order Online »](#)

BML-G244-0050	50mg
BML-G244-0250	250mg

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

Appearance	White solid.
CAS	79902-63-9
Couple Target	HMG-CoA reductase
Couple Type	Inhibitor
Formula	$C_{25}H_{38}O_5$
MI	14: 8539
MW	418.6
Purity	≥98% (TLC)
RTECS	EK7798000
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml); insoluble in water.

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)