

# Simvastatin . sodium salt

**Competitive inhibitor of HMG-CoA reductase.**  
**Suppresses TNF-induced NF-κB activation.**

Simvastatin is marketed under the trade name Zocor™ and is often prescribed in combination with ezetimibe (Zetia) to treat dyslipidemia. This drug combination is known commercially as Vytorin or Inegy. Simvastatin also suppresses TNF-induced NF-κB activation, which potentiates apoptosis in human myeloid leukemia cells and thus, may be useful in treating cancer.

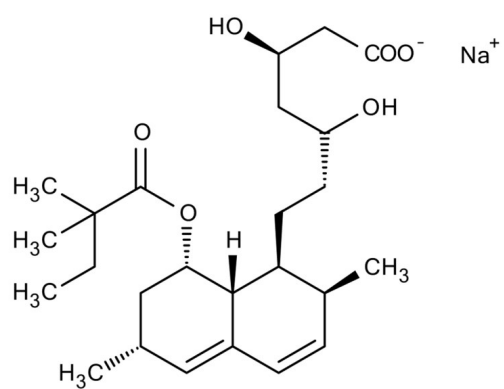
## Ordering Information

[Order Online »](#)

ENZ-CHM142-0005	5mg
ENZ-CHM142-0010	10mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	(βR, αR, 1S, 2S, 6R, 8S, 8aR)-8-[(2, 2-dimethyl-1-oxobutylideneamino)oxy]-1, 2, 6, 7, 8, 8a-hexahydro-β, δ-dihydroxy-2, 6-dimethylnaphthaleneheptanoic acid . Na monosodium salt
Appearance	White powder solid.
CAS	101314-97-0
Couple Target	HMG-CoA reductase
Couple Type	Inhibitor
Formula	C <sub>25</sub> H <sub>39</sub> O <sub>6</sub> . Na
MW	458.6
Purity	≥97% (HPLC)
Solubility	Soluble 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)