

# 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 »](#)

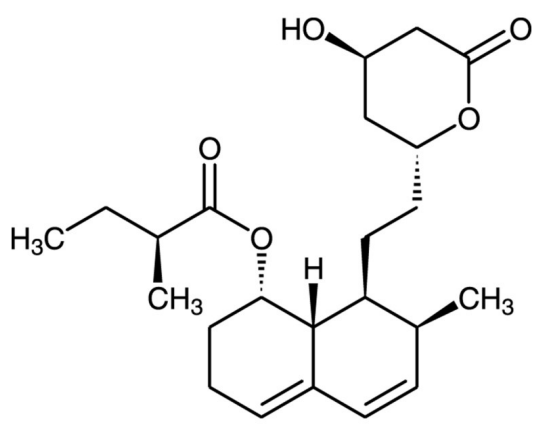
## Ordering Information

[Order Online »](#)

BML-G233-0010	10mg
---------------	------

## 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-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)