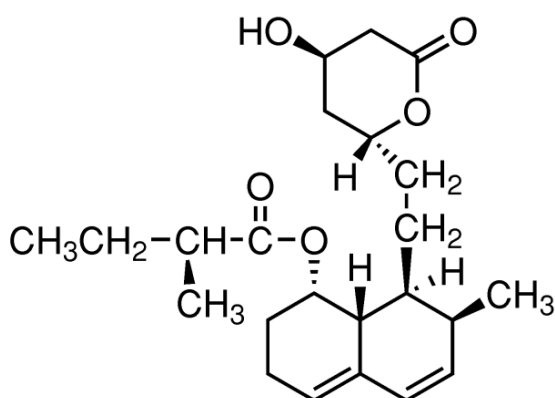


## Mevastatin

### statin

### HMG-CoA reductase inhibitor

Cat. No.	Amount
INH-014	1 mg



For *in vitro* use only  
 Quality guaranteed for 12 months  
 Store at 4°C

#### Synonym

Compactin

#### Form

White solid

#### Molecular weight

390.51 g/mol

#### Molecular Formula

C<sub>23</sub>H<sub>34</sub>O<sub>5</sub>

#### Purity

>96% by HPLC

#### CAS Number

73573-88-3

#### Description

Potent inhibitor of 3-hydroxy-3-methylglutaryl-CoA reductase, HMG-CoA reductase.

#### Selected References:

Nguyen *et al.* (2009) Analysis of the eukaryotic prenylome by isoprenoid affinity tagging. *Nature Chemical Biology*. **5** (4):227

Hasumi *et al.* (1987) Overaccumulation of 3-hydroxy-3-methylglutaryl-coenzyme-A reductase in a compactin(ML-236B)-resistant mouse cell line with defects in the regulation of its activity. *Eur J Biochem*. **164** (3):547