

# Amlodipine

Ca<sup>2+</sup> channel blocker

Amlodipine is a third generation L-type calcium channel blocker with a unique spectrum of actions, over and above its antihypertensive activity which include stimulation of NO synthase, inhibition of monocyte-endothelial adhesion, inhibition of oxidative stress and sustained duration of action.

Citations: 6

[View Online »](#)

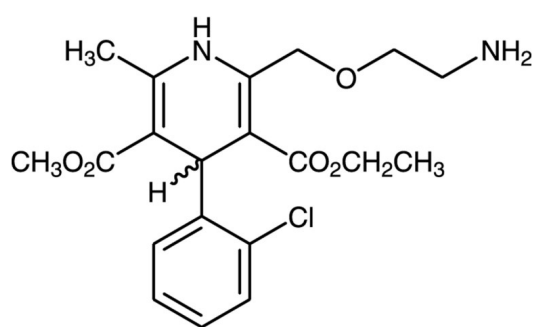
## Ordering Information

[Order Online »](#)

BML-CA202-0100	100mg
BML-CA202-1000	1g

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** Ambient

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** Almost white crystal powder

**CAS** 88150-42-9

**Couple Target** Calcium channel

**Couple Type** Blocker

**Formula**  $C_{20}H_{25}ClN_2O_5$

**MW** 408.9

**Purity** ≥98% (Assay)

**Solubility** Soluble in methanol, DMSO (25mg/ml), 100% ethanol (25mg/ml), dimethyl formamide (25mg/ml) or chloroform.



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)