# **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 »

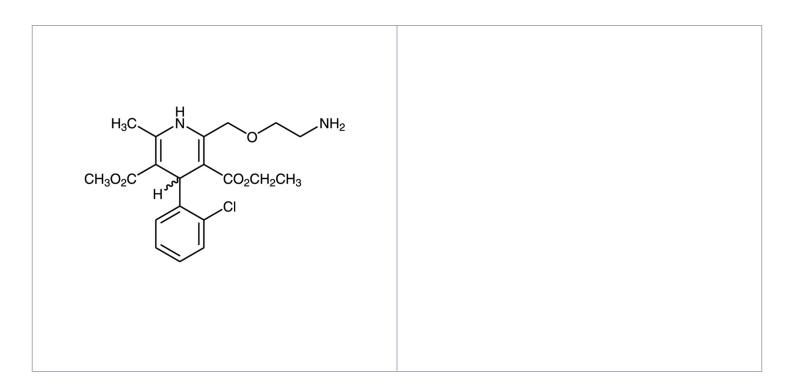
## **Ordering Information**

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

Blocker **Couple Type** 

**Formula**  $C_{20}H_{25}CIN_2O_5$ 

MW 408.9

**Purity** ≥98% (Assay)

Solubility Soluble in methanol, DMSO (25mg/ml), 100% ethanol

(25mg/ml), dimethyl formamide (25mg/ml) or chloroform.

Last modified: May 29, 2024

info-

eu@enzolifesciences.com