

# Capsaicin

TRPV1 agonist

Capsaicin is an agonist at the TRPV1 receptor, a ligand-gated non-selective cation channel. Exhibits neuroprotective effects in a model of transient global cerebral ischemia. Induces release of substance P from afferent nociceptive specific neurons. Constituent of cayenne pepper. Powerful excitant of peripheral sensory nerve endings, specifically unmyelinated afferent neurons (C-fibers). Reversibly inhibits aggregation of platelets.

Citations: 12

[View Online »](#)

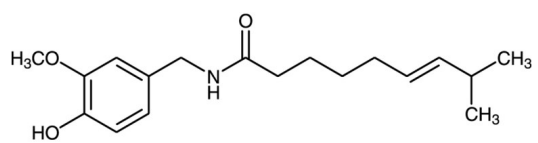
## Ordering Information

[Order Online »](#)

BML-EI125-0200	200mg
BML-EI125-1000	1g

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 2 years. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	8-Methyl-N-vanillyl-6-nonenamide
Appearance	White to pale yellow.
CAS	404-86-4
Couple Target	TRP channel
Couple Type	Activator
Formula	$C_{18}H_{27}NO_3$
MW	305.4
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, methanol or chloroform. Sparingly soluble in water.
Source	Isolated from <i>Capsicum</i> fruit.



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)