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 peropheral sensory nerve endings, specifically unmyelinated afferent neurons (C-fibers). Reversibly inhibits aggregation of platelets.

Citations: 12

View Online »

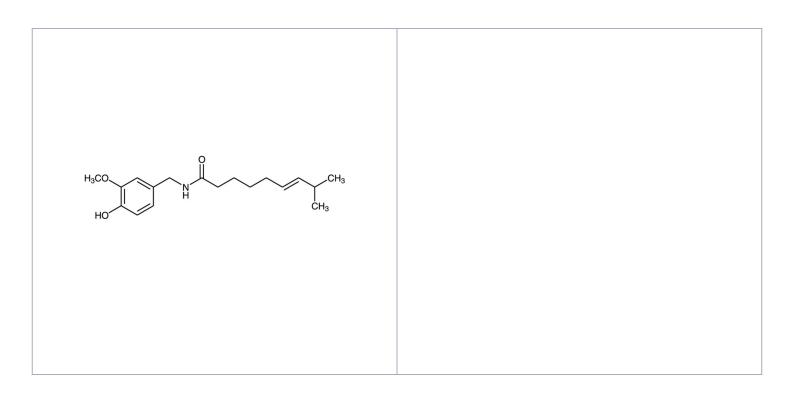
Ordering Information

Order Online »

BML-El125-0200	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 Capsicum fruit.

Last modified: May 29, 2024

