

Gabapentin

Ca²⁺ channel blocker

Gabapentin is a GABA analog which does not act at GABA receptors but which inhibits L-type calcium channels via binding to the α_2 - δ subunit and neuropathic pain relieving actions. Ligands acting at the α_2 - δ subunit lead to anticonvulsant, anxiolytic and neuropathic pain-relieving actions.

Citations: 3

[View Online »](#)

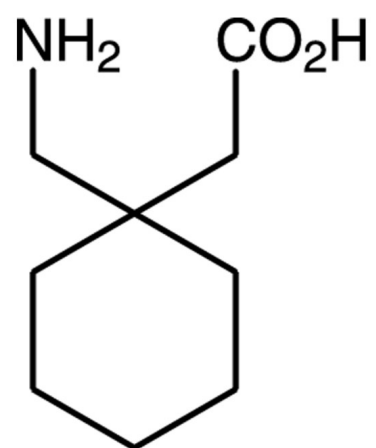
Ordering Information

[Order Online »](#)

BML-CA234-0100	100mg
----------------	-------

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	White solid.
CAS	60142-96-3
Couple Target	Calcium channel
Couple Type	Blocker
Formula	$C_9H_{17}NO_2$
MW	171.2
Purity	≥98% (HPLC)
Solubility	Soluble in water.

Last modified: May 29, 2024



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)