

(±)Epibatidine

Nicotinic acetylcholine receptor ligand

Epibatidine is an analgesic alkaloid isolated from the skin of the poison frog *Epipedobates tricolor*. It is the most potent nicotinic acetylcholine receptor agonist known. The natural (+) enantiomer is only 2-fold more potent than the (-) enantiomer with respect to analgesic activity in mice. Both enantiomers bind with high affinity ($K_i=0.045$ and 0.058 nM respectively) to rat brain membrane nicotinic receptors.

Citations: 2

[View Online »](#)

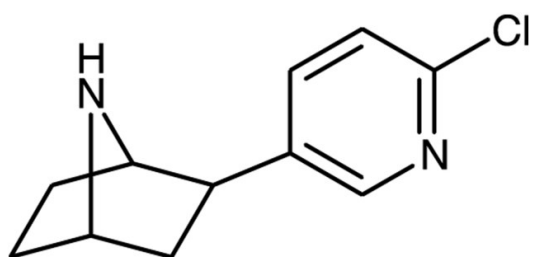
Ordering Information

[Order Online »](#)

BML-C114-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White solid.

CAS 140111-52-0

Couple Target Acetylcholine receptor

Couple Type Ligand

Formula $C_{11}H_{13}ClN_2$

MW 208.7

Purity ≥99% (HPLC)

Solubility Soluble in water (5 mg/ml) or DMSO (20 mg/ml).

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)