

PNU-282987

Nicotinic acetylcholine receptor activator

A highly selective $\alpha 7$ nicotinic acetylcholine receptor agonist and prototypical novel therapeutic for schizophrenia.

Citations: 2

[View Online »](#)

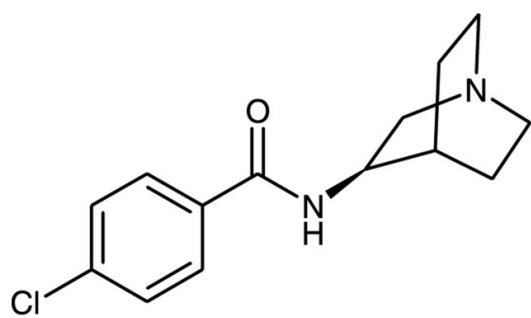
Ordering Information

[Order Online »](#)

BML-C120-0010	10mg
---------------	------

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 one year. Store solutions at -20°C for up to three months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white solid
CAS	123464-89-1
Couple Target	Acetylcholine receptor
Couple Type	Activator
Formula	$C_{14}H_{17}ClN_2O$
MW	264.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (20 mg/ml), water (6 mg/ml) or ethanol (6 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)