

Anabesine . hydrochloride

Acetylcholine receptor activator

Tobacco alkaloid. Potent nicotinic acetylcholine receptor agonist.
Insecticide.

Citations: 5

[View Online »](#)

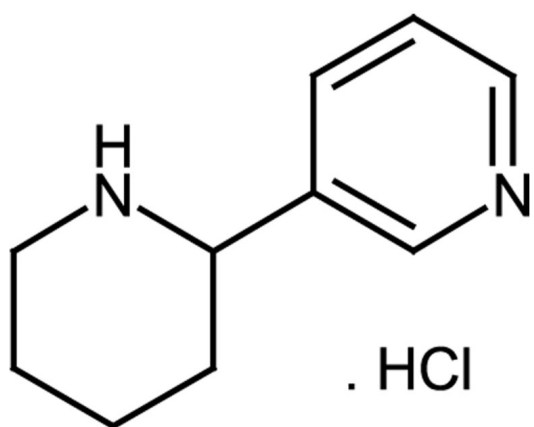
Ordering Information

[Order Online »](#)

ALX-350-112-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-(2-Piperidiny)pyridine, 2-(3-Pyridyl)piperidine, Neonicotine
Appearance	White to off-white solid.
CAS	15251-47-5
Couple Target	Acetylcholine receptor
Couple Type	Activator
Formula	$C_{10}H_{14}N_2 \cdot HCl$
MI	14: 619
MW	162.2 . 36.5
Purity	≥97% (TLC, NMR)
Solubility	Soluble in alcohol or water.
Source	Isolated from <i>Anabasis aphylla</i> .

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com