

Galanthamine . hydrobromide

Acetylcholine esterase inhibitor

Specific, competitive and reversible acetylcholinesterase inhibitor.
Alzheimer's disease therapeutic.

Citations: 7

[View Online »](#)

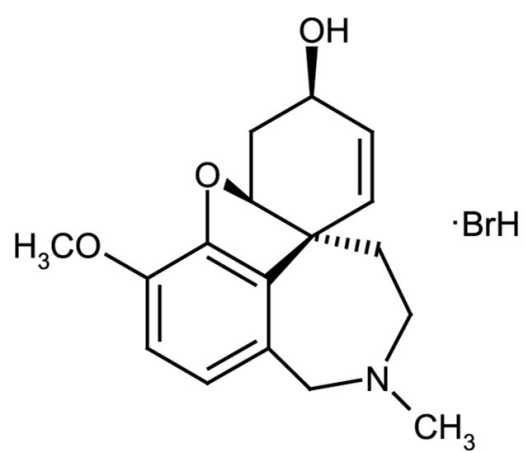
Ordering Information

[Order Online »](#)

ALX-550-336-M050	50mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from moisture.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Galantamine, Nivalin
Appearance	White to off-white crystalline solid.
CAS	1953-04-4
Couple Target	Acetylcholine esterase
Couple Type	Inhibitor
Formula	$C_{17}H_{21}NO_3 \cdot HBr$
MI	14: 4340
MW	287.4 . 80.9
Purity	≥98% (HPLC)
RTECS	DF8075000
Solubility	Soluble in DMSO (10mg/ml) or 100% ethanol (4mg/ml); sparingly 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-
fr@enzolifesciences.com](mailto:info-fr@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)