

(±)-Huperzine A

Acetylcholine esterase inhibitor

Racemic modification of (-)-Huperzine A (Prod. No. ALX-550-065) with acetylcholinesterase inhibitor activity. Has a weaker biological activity than (-)-Huperzine A. Neurotoxic.

Citations: 11

[View Online »](#)

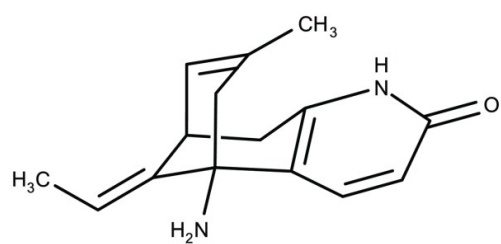
Ordering Information

[Order Online »](#)

ALX-550-064-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (±)-Selagine

Appearance Off-white solid.

CAS 120786-18-7

Couple Target Acetylcholine esterase

Couple Type Inhibitor

Formula $C_{15}H_{18}N_2O$

MI **14:** 4755

MW 242.3

Purity ≥95%

Solubility Soluble in dilute aqueous acids or chloroform.

Source Synthetic.

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)