

(+)-Himbacine

Muscarinic acetylcholine receptor antagonist

Potent muscarinic antagonist that displays selectivity for the M2 or M4 receptors.

Citations: 8

[View Online »](#)

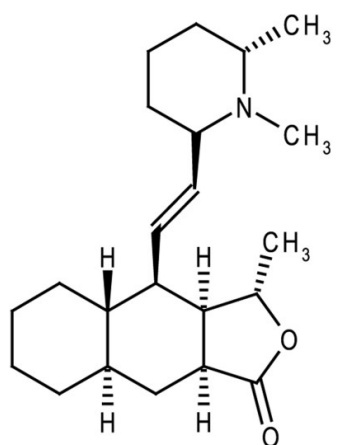
Ordering Information

[Order Online »](#)

ALX-550-061-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Handling Hygroscopic.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (3S,3aR,4R,4aS,8aR,9aS)-4-[(1E)-2-[(2R,6S)-1,6-Dimethyl-2-piperdiny]ethenyl]decahydro-3-methylnaphtho[2,3-c]furan-1(3H)-one

Appearance White to off-white powder.

CAS 6879-74-9

Couple Target Acetylcholine receptor

Couple Type Inhibitor

Formula $C_{22}H_{35}NO_2$

MW 345.5

Purity $\geq 98\%$

RTECS QL0843500

Solubility Soluble in methanol, 100% ethanol or methylene chloride; insoluble in water.

Source Isolated from the Australian pine *Galbulimima baccata*.



ENZO LIFE SCIENCES,
Last modified: **May 29, 2024**
Phone: 800.942.0430
info-usa@enzolifesciences.com

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