

Naloxone . hydrochloride

Opioid receptor antagonist

Opioid antagonist.

Citations: 1

[View Online »](#)

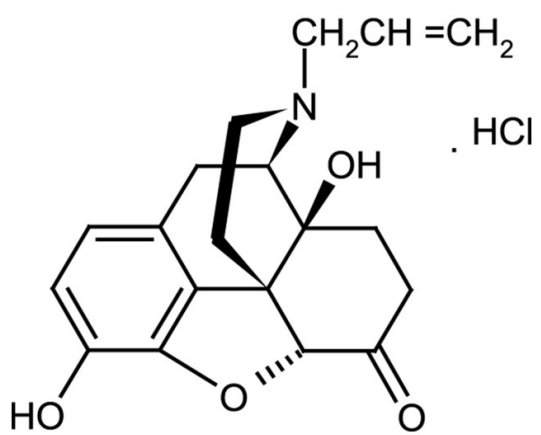
Ordering Information

[Order Online »](#)

ALX-550-055-M050	50mg
------------------	------

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	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(5α)-17-allyl-4,5-epoxy-3,14-dihydroxymorphinan-6-one hydrochloride
Appearance	White to off-white solid.
CAS	357-08-4
Couple Target	Opioid receptor
Couple Type	Inhibitor
Formula	$C_{19}H_{21}NO_4 \cdot HCl$
MI	14: 6362
MW	363.8
Purity	≥98% (HPLC)
Solubility	Soluble in water or 100% ethanol.

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)