

nor-NOHA . acetate

Arginase inhibitor

Potent and reversible arginase inhibitor. Specifically interacts with the manganese-cluster of the arginase active site. Does not act as a substrate or as an inhibitor of NOS.

Citations: 2

[View Online »](#)

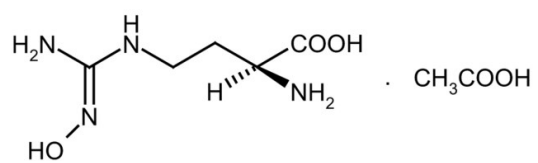
Ordering Information

[Order Online »](#)

ALX-270-467-M001	1mg
ALX-270-467-M005	5mg
ALX-270-467-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Very hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N ω -Hydroxy-nor-L-arginine . acetate
Appearance	White to slightly yellowish powder.
CAS	189302-40-7
Couple Target	Arginase
Couple Type	Inhibitor
Formula	C ₅ H ₁₂ N ₄ O ₃ . CH ₃ COOH
MW	176.2 . 60.1
Purity	≥97% (TLC)
Solubility	Soluble in water (50mg/ml).

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)