

L-NIL . dihydrochloride

NOS inhibitor

Selective inhibitor of inducible nitric oxide synthase (iNOS; NOS II). Shows anti-inflammatory activity.

Citations: 6

[View Online »](#)

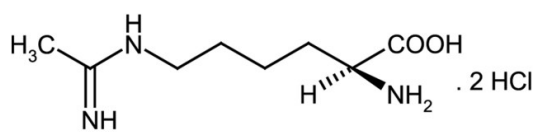
Ordering Information

[Order Online »](#)

ALX-270-010-M010	10mg
ALX-270-010-M050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	L-N6-(1-Iminoethyl)-lysine . 2HCl
Appearance	White to off-white solid.
CAS	159190-45-1
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_8H_{17}N_3O_2 \cdot 2HCl$
MW	187.2 . 73.0
Purity	≥97% (TLC)
Solubility	100mg/ml soluble in water or methanol.

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)