

N- ω -Propyl-L-arginine

nNOS inhibitor

Potent and selective inhibitor of neuronal nitric oxide synthase (nNOS; NOS I) relative to inducible nitric oxide synthase (iNOS; NOS II) (3158-fold) and endothelial nitric oxide synthase (eNOS; NOS III) (149-fold).

Citations: 9

[View Online »](#)

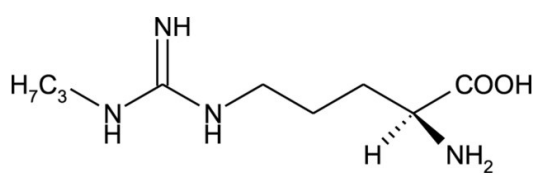
Ordering Information

[Order Online »](#)

ALX-270-203-M025	25mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year when stored at -20°C. We do not recommend storing aqueous solutions for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N5-[Imino(propylamino)methyl]-L-ornithine
Appearance	White to off-white solid.
CAS	137361-05-8
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_9H_{20}N_4O_2$
MW	216.3
Purity	≥98% (TLC)
Solubility	Soluble in PBS, pH 7.2 (25mg/ml), 100% ethanol (10mg/ml), DMSO (50µg/ml) or dimethyl formamide (50µg/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)