

Vinyl-L-NIO

nNOS inhibitor

Very selective and potent neuronal nitric oxide synthase (nNOS/NOS I) inhibitor.

Citations: 10

[View Online »](#)

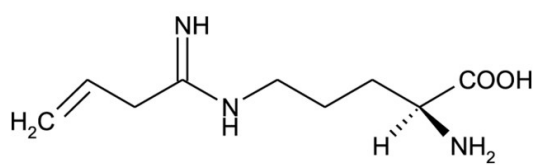
Ordering Information

[Order Online »](#)

ALX-270-216-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. Open only in anhydrous atmosphere.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N5-(1-Imino-3-butenyl)-L-ornithine, L-VNIO
Appearance	Off-white solid to clear film.
CAS	728944-69-2
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_9H_{17}N_3O_2 \cdot HCl$
MW	235.7
Purity	≥95% (TLC)
Solubility	Soluble in water, 100% ethanol (30mg/ml), DMSO (50mg/ml) or dimethyl formamide (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)