

1400W . dihydrochloride

iNOS inhibitor

A slow, tight binding and highly selective inhibitor of inducible nitric oxide synthase (iNOS/NOS II) *in vitro* and *in vivo*. Binds tightly to iNOS with a K_d of 7nM. In contrast, inhibition of nNOS and eNOS is weaker and rapidly reversible with K_i values of 2 and 50 μ M respectively. Inhibits tumor growth *in vivo*. Prevents circulatory failure but not organ injury induced by endotoxin. Displays protective effects in oxygen-deprived brain slices.

Citations: 22

[View Online »](#)

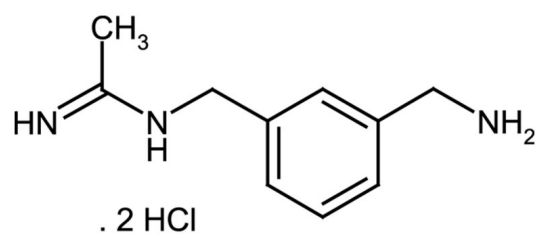
Ordering Information

[Order Online »](#)

ALX-270-073-M005	5mg
ALX-270-073-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	214358-33-5
Couple Target	NOS
Couple Type	Inhibitor
Formula	$C_{10}H_{15}N_3 \cdot 2HCl$
Identity	Determined by NMR.
MW	250.17
Purity	≥95% (TLC)
Solubility	Soluble in water (up to 25mg/ml), methanol or DMSO.
Source	Synthetic.

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)