

L-NMMA . monoacetate

NOS inhibitor

Competitively inhibits the generation of NO from arginine and is a useful tool for inhibition of NO mediated effects. *In vitro*, L-NMMA inhibits relaxation of rat aorta rings induced by acetylcholine ($IC_{50}=9.5\mu M$). *In vivo*, it causes a dose dependent increase in mean arterial blood pressure in anesthetized rats ($EC_{50}=16.7mg/kg$).

Citations: 13

[View Online »](#)

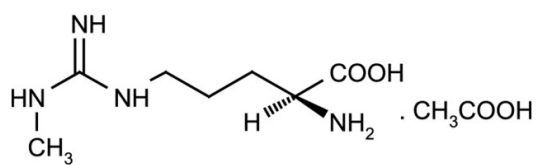
Ordering Information

[Order Online »](#)

ALX-106-001-M025	25mg
ALX-106-001-M100	100mg
ALX-106-001-G001	1g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NG-Monomethyl-L-arginine . monoacetate

Appearance White to off-white solid.

CAS 53308-83-1

Couple Target NOS

Couple Type Inhibitor

Formula $C_7H_{16}N_4O_2 \cdot CH_3COOH$

MW 188.2 . 60.1

Purity ≥99% (TLC)

Solubility Soluble in water, 100% ethanol (1mg/ml) or DMSO (0.25mg/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)