

L-NAME . hydrochloride

NOS inhibitor

Inhibitor of nitric oxide synthase (NOS). More soluble than N^G-Nitro-L-arginine (Prod. No. ALX-105-001). Suppresses the release of nitric oxide (NO) from endothelial cells. Has a different effect on the superoxide generation by NOS than L-NMMA (Prod. No. ALX-106-001).

Citations: 18

[View Online »](#)

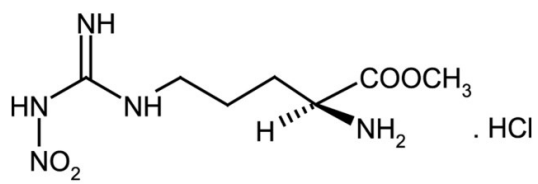
Ordering Information

[Order Online »](#)

ALX-105-003-G005	5g
ALX-105-003-G025	25g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name NG-Nitro-L-arginine-methyl ester . HCl

Appearance White to off-white solid.

CAS 51298-62-5

Couple Target NOS

Couple Type Inhibitor

Formula $C_7H_{15}N_5O_4 \cdot HCl$

MW 233.2 . 36.5

Purity ≥98% (TLC)

Solubility Soluble in water (50mg/ml) or methanol (25g/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)