

# D-NMMA . monoacetate

NOS inhibitor

D-NMMA is inactive as an inhibitor of NOS and may be used to explore non-specific effects of L-NMMA.

Citations: 2

[View Online »](#)

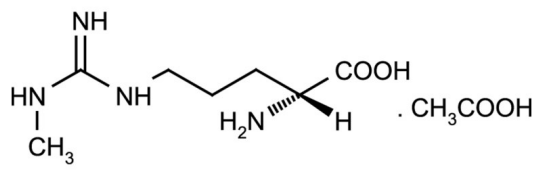
Ordering Information

[Order Online »](#)

ALX-106-002-M025	25mg
------------------	------

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-D-arginine . monoacetate

**Appearance** White solid.

**CAS** 137694-75-8

**Couple Target** NOS

**Couple Type** Inhibitor

**Formula**  $C_7H_{16}N_4O_2 \cdot CH_3COOH$

**MW** 188.2 . 60.1

**Purity** ≥99%

**Solubility** Soluble in water.

**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-  
be@enzolifesciences.com](mailto:info-be@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)