

PRODUCT INFORMATION

D-NMMA (acetate)

Item No. 14186

CAS Registry No.: 137694-75-8

Formal Name: N⁵-[imino(methylamino)methyl]-D-ornithine, monoacetate

Synonyms: D-N^G-monomethyl Arginine Acetate

MF: C₇H₁₆N₄O₂ • C₂H₄O₂

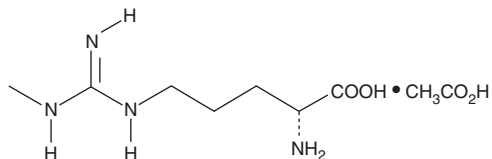
FW: 248.3

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

D-NMMA (acetate) is supplied as a solid. A stock solution may be made by dissolving the D-NMMA (acetate) in the solvent of choice, which should be purged with an inert gas. D-NMMA (acetate) is soluble in methanol and water. We do not recommend storing the aqueous solution for more than one day.

Description

D-NMMA is the inactive enantiomer of the competitive NOS inhibitor, L-NMMA (Item No. 10005031). It fails to inhibit NOS at concentrations as high as 100 μM and thus, may be used to explore non-specific effects of L-NMMA.^{1,2}

References

1. Rees, D.D., Palmer, R.M.J., Hodson, H.F., *et al.* A specific inhibitor of nitric oxide formation from L-arginine attenuates endothelium-dependent relaxation. *Br. J. Pharmacol.* **96**(2), 418-424 (1989).
2. Bogle, R.G., Moncada, S., Pearson, J.D., *et al.* Identification of inhibitors of nitric oxide synthase that do not interact with the endothelial cell L-arginine transporter. *Br. J. Pharmacol.* **105**(4), 768-770 (1992).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/21/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM