

PRODUCT INFORMATION



L-NMMA (citrate)

Item No. 80200

Formal Name: N⁵-[imino(methylamino)methyl]-L-ornithine, citrate

Synonym: L-NG-monomethyl Arginine citrate

MF: 3[C₇H₁₆N₄O₂] • C₆H₈O₇

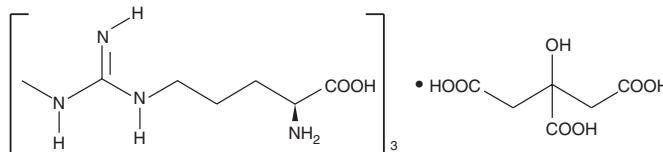
FW: 756.8

Purity: ≥99%

Supplied as: A crystalline 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

L-NMMA (citrate) is supplied as a crystalline solid. Aqueous solutions of L-NMMA (citrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of L-NMMA (citrate) in PBS (pH 7.2) is approximately L-NMMA (citrate) mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

L-NMMA is a relatively non-selective inhibitor of all NOS isoforms. It is the archetypal NOS inhibitor to which other inhibitors are most often compared.¹ The K_i values for nNOS (rat), eNOS (human), and iNOS (mouse) are approximately 0.18, 0.4, and 6 μM, respectively.^{2,3}

References

1. Griffith, O.W. and Kilbourn, R.G. Nitric oxide synthase inhibitors: Amino acids. *Methods Enzymol.* **268**, 375-392 (1996).
2. Frey, C., Narayanan, K., McMillan, K., *et al.* L-Thiocitrulline. A stereospecific, heme-binding inhibitor of nitric-oxide synthases. *J. Biol. Chem.* **269**(42), 26083-26091 (1994).
3. Garvey, E.P., Tuttle, J.V., Covington, K., *et al.* Purification and characterization of the constitutive nitric oxide synthase from human placenta. *Arch. Biochem. Biophys.* **311**(2), 235-241 (1994).

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, 04/08/2024

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