

## L-NMMA

A competitive inhibitor of nitric oxide synthase in a wide range of cells.



## **Product attributes**

CAS number	53308-83-1
Molecular weight	248.3
Storage Conditions	Store at 2 to 8 °C, Desiccate

Call us: 800-304-5357 Email: btinfo@biotium.com

## **Product Description**

L-NMMA (NG-Methyl-L-arginine, acetate salt) competitively inhibits nitric oxide synthase in a wide range of cells to study the role of nitric oxide in biochemical and physiological functions.

- Inhibition of the nitric oxide synthase isoforms in a wide range of cells
- White solid soluble in water

If a variety of NOS inhibitors are needed, we offer a NOS Inhibitor Kit which contains four four NOS inhibitors: 7-NI (00240) (50 mg), L-NMMA (00241) (25 mg), L-NIL (00242) (5 mg) and L-NIO (00243) (5 mg). We also have a selection of dyes and probes for detecting nitric oxide or reactive oxygen in cells.

Molecular structure:

## References

- Am J Physiol 269, E709 (1996), <u>DOI: 10.1152/ajpendo.1996.270.4.E709</u>
  Biochem Biophys Res Comm 205, 992 (1994), <u>DOI: 10.1006/bbrc.1994.2764</u>
  Nat Med 17, 1619 (2011). <u>DOI: 10.1038/nm.2544</u>

This datasheet was generated on November 6, 2025 at 11:26:02 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/l-nmma-ng-methyl-l-arginine-acetate-salt/