

## 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

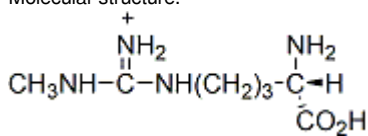
## Product Description

L-NMMA (N<sup>G</sup>-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 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

1. Am J Physiol 269, E709 (1996), [DOI: 10.1152/ajpendo.1996.270.4.E709](#)
2. Biochem Biophys Res Comm 205, 992 (1994), [DOI: 10.1006/bbrc.1994.2764](#)
3. Nat Med 17, 1619 (2011). [DOI: 10.1038/nm.2544](#)

This datasheet was generated on January 7, 2026 at 04:32:35 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/>