

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** L-NMMA ( $N^G$ -methyl-L-arginine, acetate salt)

**CATALOG #:** 00241

MOLECULAR $C_9H_{20}N_4O_4$ INFORMATION:Mwt: 248.3

[53308-83-1]

CH<sub>3</sub>NH - C - NH(CH<sub>2</sub>)<sub>3</sub> - C - OH

**PROPERTIES:** 

Color & FormWhite solid, hygroscopicPurity $\geq 98\%$  by elemental analysis.

**Soluble** in water

STORAGE AND HANDLING:

Stored desiccated at 4°C upon receipt. Protect from moisture.

**APPLICATION:** Competitively inhibits nitric oxide synthase in a wide range of cells.

Ref: Baron, et al. Am. J. Physiol. 269, E709(1995); Oddis, et al. Biochem. Biophys. Res. Comm. 205, 992(1994)

TOXICITY: Unknown

FIRST AID: Potentially harmful; avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas

with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.