

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

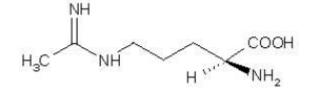
PRODUCT NAME:

L-NIO (L-*N*⁵-(1-iminoethyl)ornithine, dihydrochloride)

CATALOG #: 00243

MOLECULAR INFORMATION:

C₇H₁₇Cl₂N₃O₂ Mwt: 246.1 [36889-13-1]



PROPERTIES: Color & Form Purity Solubility	Off-white solid. Very hygroscopic. ≥ 98% by TLC (single spot) Soluble in water or MeOH (100mg/ml)
STORAGE AND HANDLING:	Stored desiccated at -20°C upon receipt and protect from moisture at all times.
APPLICATION:	A potent, irreversible inhibitor of neutrophil nitric oxide synthase.
	Ref: McCall, et al. Br. J. Pharmacol. 102(1), 234(1991); Mulligan, et al. Br. J. Pharmacol. 107(4), 1159(1992)
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.*