

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: NOS Inhibitor Kit

**CATALOG #:** 00249

MOLECULAR INFORMATION:

See Individual Components

**PROPERTIES:** 

Color & FormSee Individual ComponentsPuritySee Individual ComponentsSolubilitySee Individual Components

STORAGE AND HANDLING:

See Individual Components

APPLICATION: This kit 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).

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.