

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: DEA-NONOate** (2-(*N*,*N*-diethylamino)-diazenolate-2-oxide, diethylammonium

salt)

CATALOG #: 00225

MOLECULAR  $C_4H_{10}N_3O_2 \cdot C_4H_{12}N$  Mwt: 206.3

[56329-27-2]

 $H_3C$   $N_2$   $CH_3$   $CH_3$   $CH_3$   $CH_3$   $CH_3$ 

**PROPERTIES:** 

Color & FormOff-white solidPurity $\geq$  98% by HPLC

Solubility Soluble in DMSO (2mg/mL); 100% EtOH (25mg/mL); water; PBS pH 7.2 (2mg/mL)

STORAGE AND HANDLING:

Store desiccated under nitrogen at -20 °C.

**APPLICATION:** Releases nitric oxide spontaneously under physiological conditions. Useful for

reliable NO generation in solution.

Ref: Keefer, et al. Methods in Enzymeology 268, 281(1996); Maragos, et al. J. Med. Chem. 34(11), 3242(1991).

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.