

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:	Nitric Oxide Generation Kit
CATALOG #:	00239
MOLECULAR INFORMATION:	See information for individual components on the following pages.
PROPERTIES: Color & Form Purity Solubility	See Individual Components See Individual Components See Individual Components
STORAGE AND HANDLING:	See Individual Components
APPLICATION:	This kit contains 10 mg each of the following nitric oxide donors: SIN-1 (00221), SNAP (00222), S-nitrosoglutathione (00223), spermine NONOate (00224), and DEA-NONOate (00225).
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.



www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:

IE: SIN-1 (morpholinosydnonimine, hydrochloride)

CATALOG #: 00221 (00221-1)

MOLECULAR INFORMATION: $\begin{array}{l} C_{6}H_{10}N_{4}O_{2}.\\ Mwt: 206.7.\\ [16142-27-1]. \end{array}$



PROPERTIES:

Color & Form	White solid
Purity	\geq 95% by TLC
Solubility	Soluble in water, methanol or DMSO.

STORAGE AND Store desiccated at -20 °C. **HANDLING:**

APPLICATION: SIN-1 releases nitric oxide and forms superoxide spontaneously under physiological conditions and stimulates cyclic GMP production. It has been shown to be a potent vasodilator *in vivo* and *in vitro* and to inhibit smooth muscle cell mitogenesis and proliferation.

The relationship between NO generated from SIN-1 and intracellular calcium has been studied using fluorescent calcium indictors, which are also available from Biotium.

Also see **SNAP** (00222).

Ref: 1) Nature **364**, 626(1993); 2) J. Pharmacol. Exp. Ther. **248**, 762(1989); 3) Am J Physiol, L9 (1994); 4) Life Sci **54**, 1449 (1994); 5) Neuron **12**, 1235 (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.



www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:

ME: SNAP (*S*-nitroso-*N*-acetylpenicillamine)

CATALOG NO. 00222

MOLECULAR INFORMATION: $\begin{array}{l} C_7 H_{12} N_2 O_4 S \\ Mwt: \ 220 \\ [79032-48-7] \end{array}$



PROPERTIES: Color & Form Purity Solubility

Light green solid \geq 98% by TLC Soluble in water and DMSO

Store desiccated at -20 °C.

STORAGE AND HANDLING:

APPLICATION:

SNAP releases nitric oxide and forms superoxide spontaneously under physiological conditions and stimulates cyclic GMP production. It has been shown to be a potent vasodilator *in vivo* and *in vitro* and to inhibit smooth muscle cell mitogenesis and proliferation.

The relationship between NO generated from SNAP and intracellular calcium has been studied using fluorescent calcium indictors, which are also available from Biotium. The NO released from SNAP has also been reported to stimulate calcium-independent synaptic vesicle release, which can be detected with FM1-43 (also available from Biotium)

Also see SIN-1 (00221)

Ref: 1) Nishio, et al. Biochem. Biophys. Res. Comm. 221(1), 163(1996); 2) Gross, et al. Neuropharcology 33(7), 915(1994). 3) Am J Physiol, L9 (1994); 4) Life Sci 54, 1449 (1994); 5) Neuron 12, 1235 (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.



www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:

S-Nitrosoglutathione (N-(N-L-γ-glutamyl-S-nitroso-L-cysteinyl)-glycine)

00223 CATALOG #:

MOLECULAR INFORMATION:

 $C_{10}H_{16}N_4O_7S$ Mwt: 336 [57564-91-7]



PROPERTIES:

Color & Form Purity Solubility Absorption Extinction Coefficient	Hygroscopic pink solid ≥ 98% by TLC Soluble in water or DMSO 336nm 900
STORAGE AND HANDLING:	Spontaneously decomposes in solution. Store desiccated at -20 °C.
APPLICATION:	Similar to SNAP, <i>S</i> -nitrosogluthathione releases nitric oxide spontaneously under physiological conditions. Also see SNAP (00222)
	Ref: Biochemistry 34, 7177(1995); Br. J. Pharmacol. 107, 745(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.



www.biotium.com

PRODUCT AND SAFETY DATA SHEET

Sperimine NONOale
00224
$\begin{array}{c} C_{10}H_{26}N_{6}O_{2} \\ Mwt: 262.4 & O \\ [136587-13-8] & & // \\ N \end{array}$
H ₂ N N N N N OH
Crystalline solid ≥ 98% Soluble in water 252nm 8,500
Hygroscopic. Store desiccated at -20 °C. For long term storage, keep sealed at -80 °C, and it will be stable for at least one year
Spermine NONOate dissociates to the free amine and nitric oxide in a pH-dependent manner following first order kinetics. To initiate the release of nitric oxide, add a portion of the stock alkaline solution of Spermine NONOate to excess buffer of pH 7.0-7.4. The half-life of Spermine NONOate is 39 minutes at 37°C in 0.1 M phosphate buffer (pH 7.4). The decomposition of the NONOate is nearly instantaneous at pH 5. ¹
Ref: 1) J. Med. Chem 34 , 3242-3247 (1991); 2) Cancer Res. 53 , 564-567 (1993); 3) Proc. Natl. Acad. Sci. USA 90 , 9813-9817 (1993)
Unknown
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.



www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME:	DEA-NONOate (2-(<i>N</i> , <i>N</i> -diethylamino)-diazenolate-2-oxide, diethylammonium salt)
CATALOG #:	00225
MOLECULAR INFORMATION:	$\begin{array}{c} C_{4}H_{10}N_{3}O_{2} \cdot C_{4}H_{12}N \\ Mwt: 206.3 \\ [56329-27-2] \end{array} \\ H_{3}C \underbrace{ N_{2} }_{H_{2}}CH_{3} \underbrace{ O_{1} }_{H_{3}} \underbrace{ N_{2} }_{H_{3}}CH_{3} \underbrace{ O_{1} }_{H_{3}} \underbrace{ N_{2} }_{H_{3}}CH_{3} \\ \end{array}$
PROPERTIES: Color & Form Purity Solubility	Off-white solid ≥ 98% by HPLC 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.