

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

CATALOG #: 00223

**MOLECULAR**  $C_{10}H_{16}N_4O_7S$  **INFORMATION:** Mwt: 336

[57564-91-7]

**PROPERTIES:** 

Color & Form Hygroscopic pink solid

**Purity**  $\geq 98\%$  by TLC

**Soluble** in water or DMSO

Absorption 336nm Extinction 900

STORAGE AND HANDLING:

Spontaneously decomposes in solution. Store desiccated at -20 °C.

APPLICATION:

Similar to SNAP, S-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.

**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.