

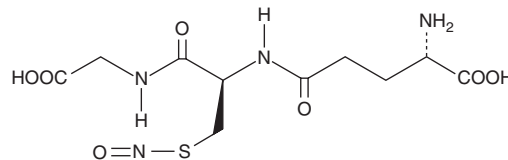
# PRODUCT INFORMATION



## S-Nitroso-L-glutathione

Item No. 82240

CAS Registry No.: 57564-91-7  
Formal Name: L-γ-glutamyl-S-nitroso-L-cysteinyl-glycine  
Synonym: GSNO  
MF:  $C_{10}H_{16}N_4O_7S$   
FW: 336.3  
Purity: ≥98%  
Supplied as: A crystalline solid  
Storage: -80°C  
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

S-Nitroso-L-glutathione is supplied as a crystalline solid. Aqueous solutions of S-nitroso-L-glutathione can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of S-nitroso-L-glutathione in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

S-Nitroso-L-glutathione is an S-nitrosothiol NO donor based on naturally-occurring glutathione. It is a potent smooth muscle relaxant and inhibitor of platelet aggregation. Its stability in solution varies from seconds to hours depending on temperature, buffer composition, and metal content.<sup>1-4</sup>

### References

1. Radomski, M.W., Rees, D.D., Dutra, A., *et al.* S-Nitroso-glutathione inhibits platelet activation *in vitro* and *in vivo*. *Br. J. Pharmacol.* **107**(3), 745-749 (1992).
2. Butler, A.R. and Rhodes, P. Chemistry, analysis, and biological roles of S-nitrosothiols. *Anal. Biochem.* **249**(1), 1-9 (1997).
3. Gasco, A., Fruttero, R., and Sorba, G. NO-donors: An emerging class of compounds in medicinal chemistry. *Farmaco* **51**(10), 617-635 (1996).
4. Cook, J.A., Kim, S.Y., Teague, D., *et al.* Convenient colorimetric and fluorometric assays for S-nitrosothiols. *Anal. Biochem.* **238**(2), 150-158 (1996).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 04/25/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM