

S-Nitroso-L-glutathione

Releases nitric oxide spontaneously under physiological conditions and plays an essential role in the biochemical and physiological functions of nitric oxide pathways.



Product attributes

Call us: 800-304-5357

CAS number	57564-91-7	
Molecular weight	336	
Storage Conditions	Store at -10 to -35 °C, Desiccate	

Product Description

Similar to SNAP, S-nitrosogluthathione (SNOG or GSNO) releases nitric oxide (NO) spontaneously under physiological conditions and plays an essential role in the biochemical and physiological functions of NO pathways.

- Generate nitric oxide and superoxides
- Pinkish solid soluble in water and DMSO

NO concentration is often assessed by measuring nitrite level using the Griess Reagent. If a variety of nitric oxide generators are needed, we offer a Nitric Oxide Generation Kit which has 10mg each of SIN-1 (00221), SNAP (00222), S-nitrosoglutathione (00223), spermine NONOate (00224), and DEA-NONOate (00225).

Molecular Structure:

$$H_2N$$
 HO_2C
 HO_2C
 HO_2C
 HO_2C

References

- Biochemistry 34, 7177 (1995), <u>DOI: 10.1021/bi00021a032</u>
 Br J Pharmacol 107, 745 (1992), <u>DOI: 10.1111/j.1476-5381.1992.tb14517.x</u>
 Chem. Res. Toxicol. 23, 2 (2010), <u>DOI: 10.1021/tx900387k</u>

This datasheet was generated on November 27, 2025 at 09:02:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/s-nitroso-l-glutathione/