

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

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

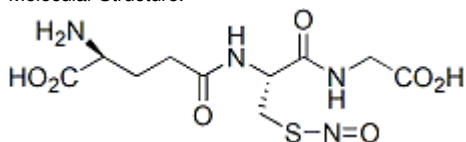
Product Description

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-nitrosogluthathione ([00223](#)), spermine NONOate ([00224](#)), and DEA-NONOate ([00225](#)).

Molecular Structure:



References

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

This datasheet was generated on June 19, 2026 at 06:03:58 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/s-nitroso-l-glutathione/>