

## 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	336
Storage Conditions	Store at -10 to -35 °C, Desiccate

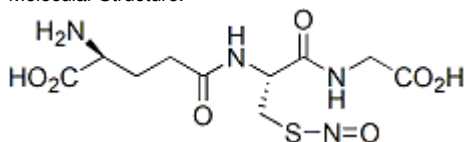
## Product Description

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](https://doi.org/10.1021/bi00021a032)
2. Br J Pharmacol 107, 745 (1992), DOI: [10.1111/j.1476-5381.1992.tb14517.x](https://doi.org/10.1111/j.1476-5381.1992.tb14517.x)
3. Chem. Res. Toxicol. 23, 2 (2010), DOI: [10.1021/tx900387k](https://doi.org/10.1021/tx900387k)

This datasheet was generated on January 17, 2026 at 10:01:57 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/s-nitroso-l-glutathione/>