

Glowing products for science

## **SNAP**

Generate nitric oxide and form superoxides spontaneously under physiological conditions and is often used to probe the cell stress response and stimulate calcium-independent synaptic vesicle release.



## Product attributes

Call us: 800-304-5357

CAS number	79032-48-7
Molecular weight	220
Storage Conditions	Store at -10 to -35 °C, Desiccate

## **Product Description**

SNAP (S-Nitroso-N-acetylpenicillamine) releases nitric oxide under physiological conditions and also been shown to be a potent vasodilator.

- Generate nitric oxide and superoxides
- Stimulates cyclic GMP production
- Light green solid soluble in water and DMSO

Fluorescent calcium indicators and SNAP are often used to probe the cellular reactions to cell stress, NO, and trace intracellular calcium. The NO released from SNAP has also been reported to stimulate calcium-independent synaptic vesicle release, which can be detected with SynaptoGreen™ (FM1-43). 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:

FM is a registered trademark of Thermo Fisher Scientific.

## References

- 1. Biochem Biophys Res Comm 221, 163 (1996), DOI: 10.1006/bbrc.1996.0563
- 2. Neuropharmacology 33, 915 (1994), DOI: 10.1016/0028-3908(94)90190-2
- 3. Am J Physiol, L9 (1994), <u>DOI: 10.1152/ajplung.1994.267.1.L9</u> 4. Life Sci 54, 1449 (1994), <u>DOI: 10.1016/0024-3205(94)00600-8</u>
- 5. Neuron 12, 1235 (1994), DOI: 10.1016/0896-6273(94)90440-5

This datasheet was generated on November 27, 2025 at 09:02:11 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/snap-s-nitroso-n-acetylpenicillamine/