

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

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

Product Description

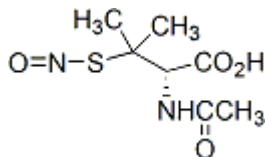
SNAP (S-Nitroso-N-acetylpenicillamine) releases nitric oxide under physiological conditions and also been shown to be a potent vasodilator.

- 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), [DOI: 10.1152/ajplung.1994.267.1.L9](#)
4. Life Sci 54, 1449 (1994), [DOI: 10.1016/0024-3205\(94\)00600-8](#)
5. Neuron 12, 1235 (1994), [DOI: 10.1016/0896-6273\(94\)90440-5](#)

This datasheet was generated on January 17, 2026 at 08:32:27 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/snap-s-nitroso-n-acetylpenicillamine/>