SIN-1

Generate nitric oxide and form superoxides spontaneously under physiological conditions and is often used to probe the cell stress response and induce cGMP production.



Product attributes

CAS number	16142-27-1
Molecular weight	206.7
Storage Conditions	Store at -10 to -35 °C or below, Desiccate

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

SIN-1 (Morpholinosydnonimine, hydrochloride) releases nitric oxide and forms superoxide spontaneously under physiological conditions; stimulates cGMP production; and has been shown to be a potent vasodilator.

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

Fluorescent calcium indicators and SIN-1 are often used to probe the cellular reactions to cell stress, NO, and trace intracellular calcium. It has been shown to be a potent vasodilator in vivo and in vitro and to inhibit smooth muscle cell mitogenesis and proliferation. 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:

References

- J Pharmacol Exp Ther 248, 762 (1989), J Pharmacol Exp Ther 248, 762 (1989)
 Nature 364, 626 (1993), DOI: 10.1038/364626a0
 Bioorg Med Chem Lett. 21, 6515 (2011), DOI: 10.1016/j.bmcl.2011.08.071

This datasheet was generated on November 4, 2025 at 10:50:44 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/sin-1-morpholinosydnonimine-hydrochloride/