

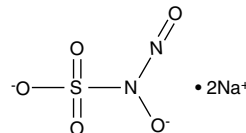
Product Information



Sulpho NONOate

Item No. 83300

Formal Name: disodium (E)-1-sulfonatodiazen-1-ium-1,2-diolate
MF: $\text{N}_2\text{O}_5\text{S} \cdot 2\text{Na}$
FW: 186.1
Purity: $\geq 90\%$
Stability: ≥ 1 year at -80°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 259 nm
Misc: Moisture, light, and oxygen sensitive



Laboratory Procedures

For long term storage, keep Sulpho NONOate sealed under nitrogen at -80°C . It should be stable for at least one year. The crystals are sensitive to moisture and should be stored in an inert atmosphere.

Sulpho NONOate dissociates to sulfate and nitrous oxide in a pH-dependent manner following first order kinetics. The decomposition is also catalyzed by borate anion.¹ Alkaline solutions of NONOates (in 0.01 M NaOH) are very stable and can be stored at 0°C for 24 hours. Sulpho NONOate is highly soluble in water and relatively concentrated solutions can be prepared for further dilution. To initiate the release of nitrous oxide, add a portion of the stock alkaline solution of Sulpho NONOate to excess buffer of pH 7.0-7.4. The half-life of Sulpho NONOate is 7 minutes at 37°C in 0.1 M phosphate buffer (pH 7.4). The decomposition of the NONOate is nearly instantaneous at pH 5.²

The intact Sulpho NONOate has a characteristic UV absorbance with an absorbance maximum at 259 nm ($\epsilon = 8,500 \text{ M}^{-1}\text{cm}^{-1}$), permitting quantitation in aqueous solutions.² The concentration of the basic stock solution of Sulpho NONOate can be measured by UV if there is any uncertainty about the condition under which it was prepared or stored.

Unlike other NONOates, Sulpho NONOate produces nitrous oxide but no nitric oxide at physiological pH. Therefore, Sulpho NONOate may be used as a negative control in experiments using other nitric oxide releasing NONOates.¹⁻³

References

1. Switkes, E.G., Dasch, G.A., and Ackermann, M.N. Kinetics of boric acid catalyzed decomposition of the *N*-nitrosohydroxylamine-*N*-sulfonate ion. *Inorg. Chem.* **12**, 1120-1123 (1973).
2. Maragos, C.M., Morley, D., Wink, D.A., *et al.* Complexes of $\cdot\text{NO}$ with nucleophiles as agents for the controlled biological release of nitric oxide. Vasorelaxant effects. *J. Med. Chem.* **34**, 3242-3247 (1991).
3. Maragos, C.M., Wang, J.M., Hrabie, J.A., *et al.* Nitric oxide/nucleophile complexes inhibit the *in vitro* proliferation of A375 melanoma cells via nitric oxide release. *Cancer Res.* **53**, 564-568 (1993).

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/19/2012