

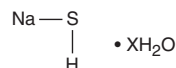
# PRODUCT INFORMATION



## Sodium Hydrogen Sulfide (hydrate)

Item No. 10012555

**CAS Registry No.:** 207683-19-0  
**Formal Name:** sodium sulfide, hydrate  
**Synonyms:** NaHS, NSC 158264, Sodium Hydrosulfide  
**MF:** HNaS • XH<sub>2</sub>O  
**FW:** 56.1  
**Purity:** ≥65%  
**UV/Vis.:** λ<sub>max</sub>: 224 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Sodium hydrogen sulfide (hydrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the sodium hydrogen sulfide (hydrate) in the solvent of choice, which should be purged with an inert gas. Sodium hydrogen sulfide (hydrate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of sodium hydrogen sulfide (hydrate) in these solvents is approximately 3 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of sodium hydrogen sulfide (hydrate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of sodium hydrogen sulfide (hydrate) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Hydrogen sulfide (H<sub>2</sub>S) is, like nitric oxide, an important gaseous mediator that has significant effects on the immunological, neurological, cardiovascular and pulmonary systems of mammals. Sodium hydrogen sulfide is an H<sub>2</sub>S donor commonly used in cellular and whole animal experimental systems. For example, it has been used to suggest that H<sub>2</sub>S promotes neutrophil migration, reduces airway inflammation, and protects neurites, heart, and intestine from chemical or ischemic-reperfusion damage.<sup>1-5</sup>

### References

1. Dal-Secco, D., Cunha, T.M., Freitas, A., *et al.* Hydrogen sulfide augments neutrophil migration through enhancement of adhesion molecule expression and prevention of CXCR2 internalization: Role of ATP-sensitive potassium channels. *J. Immunol.* **181(6)**, 4287-4298 (2008).
2. Chen, Y.H., Wu, R., Geng, B., *et al.* Endogenous hydrogen sulfide reduces airway inflammation and remodeling in a rat model of asthma. *Cytokine* **45(2)**, 117-123 (2009).
3. Hu, L.F., Lu, M., Wu, Z.Y., *et al.* Hydrogen sulfide inhibits rotenone-induced apoptosis via preservation of mitochondrial function. *Mol. Pharmacol.* **75(1)**, 27-34 (2009).
4. Yong, Q.C., Lee, S.W., Foo, C.S., *et al.* Endogenous hydrogen sulphide mediates the cardioprotection induced by ischemic postconditioning. *Am. J. Physiol. Heart Circ. Physiol.* **295(3)**, H1330-H1340 (2008).
5. Liu, H., Bai, X.B., Shi, S., *et al.* Hydrogen sulfide protects from intestinal ischaemia-reperfusion injury in rats. *J. Pharm. Pharmacol.* **61(2)**, 207-212 (2009).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/27/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD

ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM