

General Information

It has been recognized that hydrogen sulfide (H_2S) has an important role as a physiological active substance for vasodilation, cytoprotection, and modulation of insulin secretion. H_2S is considered as a gaseous molecule such as NO and CO. However, around 80% of the total sulfide exists as hydrogen sulfide anion (HS^-) under physiological condition, since the pK_a is about 7 (Fig. 1). In addition, H_2S easily converts to various biochemical molecules such as persulfides and polysulfides, which react with sulfhydryl moieties in a living body. The functional mechanism of H_2S has not been well understood. Sodium sulfide (Na_2S) has been widely used as a H_2S donor. Na_2S is readily decomposed and release H_2S when Na_2S is dissolved in H_2O .

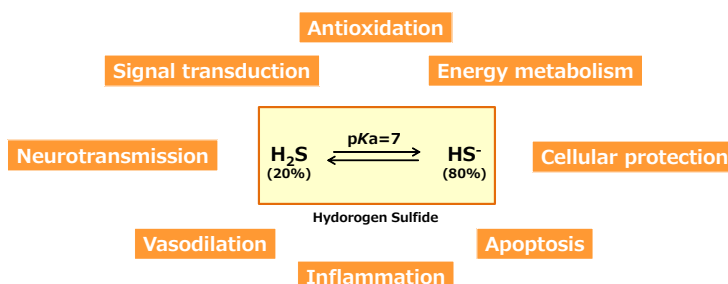


Fig. 1 Physiological functions of hydrogen sulfide

Contents

-SulfoBiotics- Sodium sulfide (Na_2S) : 100 mg x 5

Storage Condition

Store at 0-5 °C

Precaution

- * Use the reagent after bringing to room temperature, because it is moisture sensitive.
- * Use up the reagent soon after opening.

General Protocol

- 1) Dissolve 7.8 mg of Na_2S in 1 ml of deionized H_2O under nitrogen gas (100 mmol/l Na_2S solution).
 - 2) Dilute the 100 mmol/l Na_2S aqueous solution to an appropriate concentration depending on your experiment.
- * Purge deionized H_2O with nitrogen gas for longer than 30 minutes to prevent the oxidation.
 - * Use the aqueous solution as soon as prepared. The solution is not stable enough to store.

Experimental Example

- Detection of hydrogen sulfide by methylene blue method -

- 1) 20 μl of 100 mmol/l Na_2S aqueous solution was added to 980 μl of deionized H_2O to prepare 2 mmol/l Na_2S solution.
- 2) 100 μl of 2 mmol/l Na_2S solution was added to 900 μl of deionized H_2O to prepare 200 $\mu\text{mol/l}$ Na_2S solution.
- 3) 200 $\mu\text{mol/l}$ Na_2S solution was diluted with deionized H_2O to prepare various concentration of Na_2S solution by serial dilution (200, 100, 50, 25, 12.5, 6.3, 3.2, 0 $\mu\text{mol/l}$).
- 4) 300 μl of 1% zinc acetate solution, 50 μl of 20 mmol/l N,N -Dimethyl- p -phenylenediammonium (7.2 mol/l HCl) solution and 50 μl of 30 mmol/l FeCl_3 (1.2 mol/l HCl) solution were added to 250 μl of the Na_2S solutions and mixed using a vortex.
- 5) The solutions were incubated at room temperature for 30 minutes and transferred 200 μl of the solution to each well (96-well plate).
- 6) Measure the absorbance at 650 nm using a microplate reader (Fig. 2).

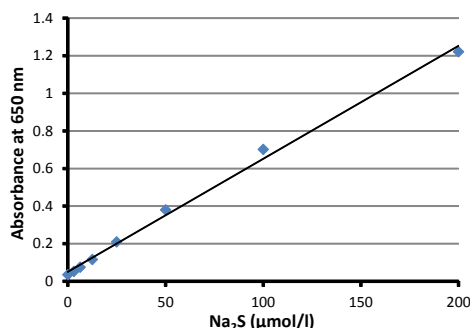


Fig. 2 Absorbance change at 650 nm depending on the concentration of hydrogen sulfide

References

- 1) R. Greiner, Z. Palinkas, K. Basell, D. Becher, H. Antelmann, P. Nagy and T. P. Dick, "Polysulfides link H_2S to protein thiol oxidation", *Antioxid. Redox Signal.*, **2013**, 19, 1749.
- 2) N. S. Lawrence, J. Davis and R. G. Compton, "Analytical strategies for the detection of sulfide: a review", *Talanta*, **2000**, 52, 771.

If you need more information, please contact Dojindo technical service.

Dojindo Laboratories
2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto
861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525
E-mail: info@dojindo.co.jp Web: www.dojindo.co.jp

Dojindo Molecular Technologies, Inc.
Tel: +1-301-987-2667 Web: <http://www.dojindo.com/>
Dojindo EU GmbH
Tel: +49-89-3540-4805 Web: <http://www.dojindo.eu.com/>
Dojindo China Co., Ltd
Tel: +86-21-6427-2302 Web: <http://www.dojindo.cn/>