

General Information

It has been recognized that hydrogen sulfide (H₂S) has an important role as a physiological active substance for vasodilation, cytoprotection, modulation of insulin secretion. Although H₂S is considered as a gaseous molecule such as NO and CO, about 80% of the total sulfide exists as hydrogen sulfide anion (HS⁻) under physiological condition. In addition, H₂S easily converts to various biochemical molecules such as persulfides and polysulfides, which react with sulfhydryl moieties in a living body. Therefore, the precise action mechanism of H₂S has not been cleared.

Hydrogen sulfide donor 5a, 8l, and 8o developed by M. Xian *et al.*, release H₂S upon reaction with reducing agents such as cysteine and glutathione *in vivo*¹⁻³⁾. The donors are stable in aqueous solutions, whereas they decompose and release H₂S in blood, cells and tissues containing sulfhydryl molecules. These hydrogen sulfide donors are expected as useful materials for clarification of H₂S roles *in vivo*.

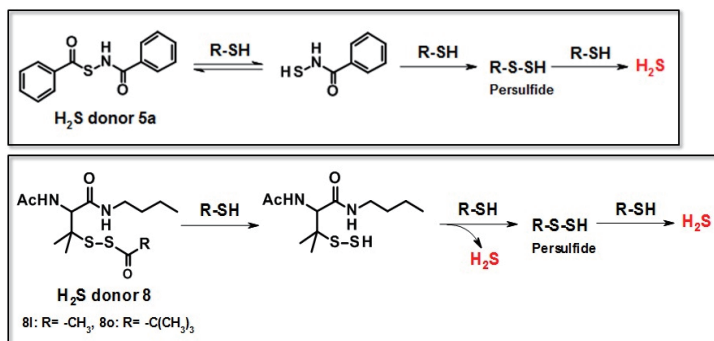


Fig. 1 H₂S release from H₂S donor 5a, 8l, 8o by reaction with thiol molecules

Contents	SB07 -SulfoBiotics- H ₂ S donor 5a 10 mg x 1
	SB08 -SulfoBiotics- H ₂ S donor 8l 10 mg x 1
	SB09 -SulfoBiotics- H ₂ S donor 8o 10 mg x 1

Storage Condition Store at 0-5 °C

Precaution **These reagents release a toxic hydrogen sulfide gas by reaction with thiol molecules. Refer to the material safety data sheet (MSDS) before using the reagent.**

General Protocol 1) Dissolve H₂S donor 10 mg in DMSO (5a: 1.95 ml, 8l: 1.56 ml, 8o: 1.38 ml) by pipetting to prepare 20 mmol/l H₂S donor Stock Solution.

*Store at -20 °C and use within two months. Aliquot the solution for longer storage.

2) Dilute the 20 mmol/l H₂S donor Stock Solution with a buffer or a medium to an appropriate concentration for your experimental condition.

Experimental Example - H₂S release profile of H₂S donor 5a, 8l, 8o by reaction with glutathione-

1) Ten microliter (10 μl) of 20 mmol/l H₂S donor Stock Solution was added to 2 ml of PBS to prepare 100 μmol/l H₂S donor (PBS) solution.

2) Glutathione was added to the 100 μmol/l H₂S donor (PBS) solution to be a final concentration of 5 mmol/l in a screw top vial. Then, the solution was incubated at room temperature.

2) The amount of yielded hydrogen sulfide anion in the solution was quantified by methylene blue method.

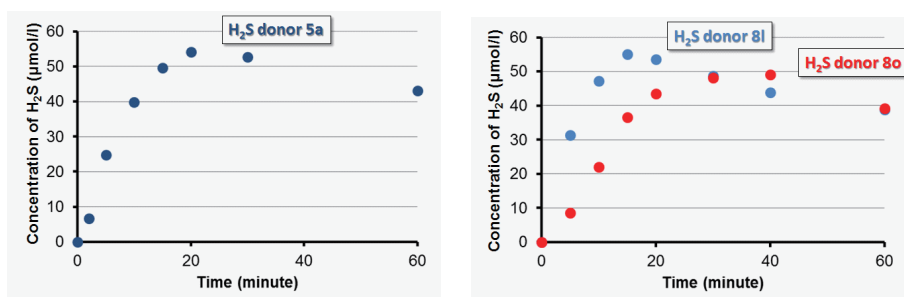


Fig. 2 H₂S release profile of 100 μmol/l H₂S donor 5a, 8l, 8o by reaction with 5 mmol/l glutathione

References

- 1) Y. Zhao, H. Wang and M. Xian, "Cysteine-Activated Hydrogen Sulfide (H₂S) Donors", *J. Am. Chem. Soc.*, **2011**, 133, 15.
- 2) C-T. Yang, Y. Zhao, M. Xian, J-H. Li, Q. Dong, H-B. Bai, J-D. Xu, M-F. Zhang, "A Novel Controllable Hydrogen Sulfide-Releasing Molecule Protects Human Skin Keratinocytes Against Methylglyoxal-Induced Injury and Dysfunction", *Cell. Physiol. Biochem.*, **2014**, 34, 1304.
- 3) Y. Zhao, S. Bhushan, C. Yang, H. Otsuka, J. D. Stein, A. Pacheco, B. Peng, N. O. Devarie-Baez, H. C. Aguilar, D. J. Lefer and M. Xian, "Controllable Hydrogen Sulfide Donors and Their Activity against Myocardial Ischemia-Reperfusion Injury", *Chem. Biol.*, **2013**, 8, 1283.

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

Dojindo Molecular Technologies, Inc.
 30 West Gude Dr., Suite 260, Rockville, MD 20850, USA
 Toll free: 1-877-987-2667 Phone: 301-987-2667 Fax: 301-987-2687
 E-mail: info@dojindo.com Web: www.dojindo.com