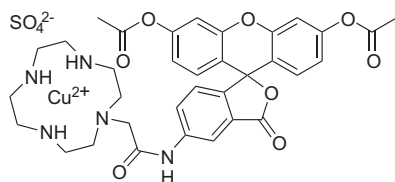


Technical Manual (Japanese version) is available at <http://www.dojindo.co.jp/manual/sb22.pdf>

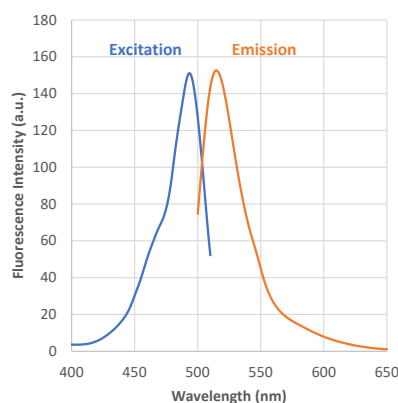
General Information

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 nitric oxide and carbon monoxide. However, around 80% of the total sulfide exists as hydrogen sulfide anion (HS^-) under physiological condition, since the pK_a is about 7. In addition, HS^- easily converts to various biochemical molecules such as persulfides and polysulfides, which react with sulfhydryl moieties in a living body. -SulfoBiotics- HSip-1 DA is cell membrane permeable and it enables fluorescent imaging of intracellular H_2S .



HSip-1 DA
(MW: 803.29)

Fig. 1 Chemical structures of HSip-1 DA



λ_{ex} : 491 nm
 λ_{em} : 516 nm

< Recommended filter >
Ex : 470 ~ 500 nm
Em : 500 ~ 550 nm

Fig. 2 Excitation and emission spectra of HSip-1 reacted with H_2S

Contents 50 μg x 1

Storage Conditions Store at $-20^\circ C$.

Required Equipment and Materials

- Dimethyl sulfoxide (DMSO)
- Serum-free medium
- HBSS
- Micropipettes

Preparation of Solutions **Preparation of 1 mmol/l HSip-1 DA stock solution**
Add 62 μl of DMSO to a tube containing 50 μg of HSip-1 DA and dissolve it by pipetting.
***Store at $-20^\circ C$. The reconstituted solution is stable at $-20^\circ C$ for 1 month.**

Experimental Example

Fluorescence imaging of hydrogen sulfide with HSip-1 DA

- 1) HeLa cells were seeded on μ -slide 8 well (Ibidi) and cultured at $37^\circ C$ overnight in a 5% CO_2 incubator.
- 2) The culture medium was discarded and the cells were washed with a serum-free medium (MEM) twice.
- 3) HSip-1 DA stock solution (1 mmol/l) was diluted with a serum-free medium (MEM) to prepare 5 $\mu mol/l$ HSip-1 DA working solution.
***Please optimize the final concentration of HSip-1 DA depending on the cell lines.**
- 4) HSip-1 DA working solution (5 $\mu mol/l$, 200 μl) was added to the cells, and the cells were cultured at $37^\circ C$ for 30 minutes in a 5% CO_2 incubator.
- 5) The supernatant was discarded, and the cells were washed with HBSS twice.
- 6) Na_2S solution (200 $\mu mol/l$, 200 μl) was added to the each well, and the cells were cultured at $37^\circ C$ for 30 minutes in a 5% CO_2 incubator.
- 7) The supernatant was discarded and the cells were washed with HBSS twice.
- 8) HBSS (200 μl) were added, and the cells were observed by confocal fluorescence microscopy.

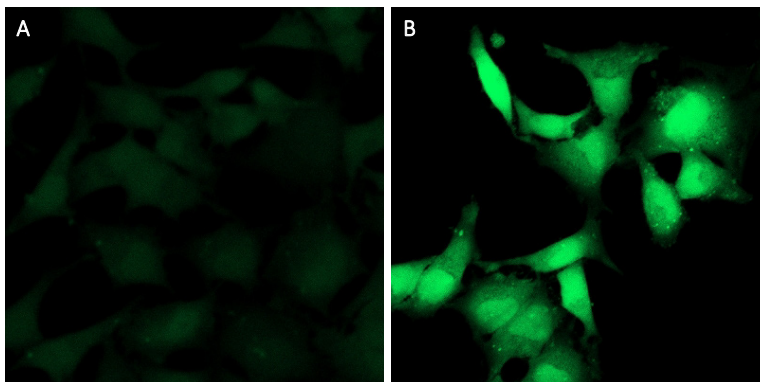


Fig.3 Detection of hydrogen sulfide using HSip-1 DA in HeLa cells treated with Na₂S.
(A: Control, B: 200 μmol/l Na₂S treated)

These products were commercialized under the advisory of Dr. Tetsuo Nagano and Dr. Kenjiro Hanaoka (The University of Tokyo).

Reference

- 1) K. Sasakura, K. Hanaoka, N. Shibuya, Y. Mikami, Y. Kimura, T. Komatsu, T. Ueno, T. Terai, H. Kimura, and T. Nagano, "Development of a Highly Selective Fluorescence Probe for Hydrogen Sulfide", *J. Am. Chem. Soc.*, **2011**, *133*, 18003.

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/>