
Product Manual

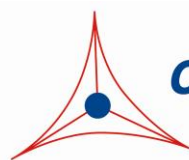
OxiSelect™ Free Hydrogen Sulfide Gas Assay Kit

Catalog Number

XAN-5084

100 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Hydrogen sulfide (H₂S) is a colorless gas containing a foul odor reminiscent of rotten eggs. High levels of H₂S are corrosive, flammable and poisonous. H₂S can be generated from the breakdown of organic matter by microbes when oxygen is not present (typically in swamps and sewers) by a process called anaerobic digestion. H₂S can also be found in natural gas, emissions from volcanoes, and in water from some wells. The human body produces minute levels of H₂S where the molecule functions as a gaseous messenger like carbon monoxide or nitric oxide. H₂S acts as a cerebrovascular dilator and is produced by the enzyme cystathione gamma lyase. Through activation of potassium ATP channels, H₂S can dilate arterioles. Furthermore, H₂S stimulates vascular smooth cell apoptosis and blocks proliferation-based vascular remodeling. Finally, H₂S has been shown to have neuroprotective and cardioprotective properties, lowering oxidative stress by increasing levels of glutathione by enhancing cystine and cysteine transport to neurons.

The OxiSelect™ Free Hydrogen Sulfide Gas Assay Kit is a novel technique for measurement of free H₂S gas emission. Liquid samples in a 96 well microplate release H₂S gas into the air which is captured in a gel spot containing Ag⁺ ions on the underside of the microplate lid just above the sample (Figure 1). The resulting silver sulfide (Ag₂S) produced inside the spot creates a brownish color (Figure 4A) that can be measured colorimetrically on a microplate spectrophotometer. Each kit provides sufficient reagents to perform up to 100 assays, including positive control and unknown samples.

Assay Principle

The OxiSelect™ Free Hydrogen Sulfide Gas Assay Kit is a 96 well microplate based kit for simple detection of H₂S gas emission. First, a solution containing silver nitrate is coated on the underside of a microplate lid to form a thin polymer capture surface (Figure 2). Samples producing H₂S such as those containing an H₂S donor, enzyme, cells, or tissue are pipetted into the wells of the microplate. The coated lid is then placed over the microplate so that the polymer rests just above each sample. H₂S gas escapes from the sample in the well and enters the polymer coating. The Ag⁺ ions in the polymer react with the H₂S gas to form Ag₂S nanoparticles, producing a brown discoloration of the polymer film. After one hour, the coated lid is transferred to another empty microplate and then OD is measured at 405 nm.

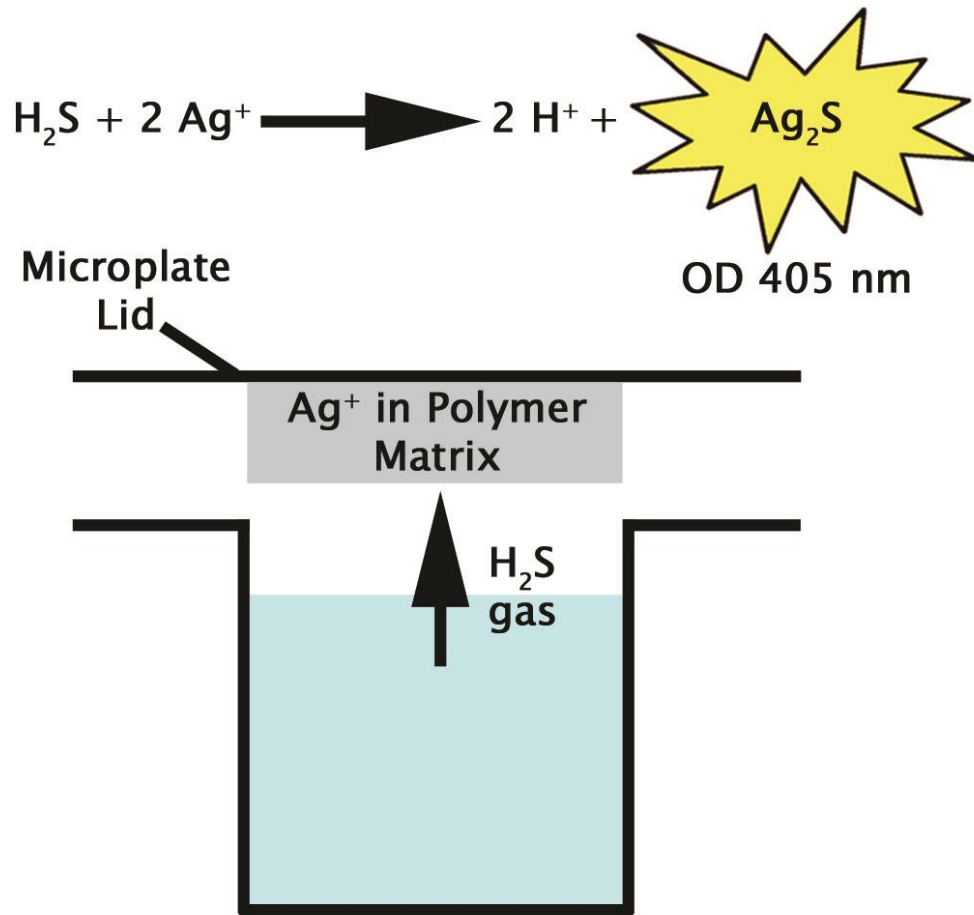
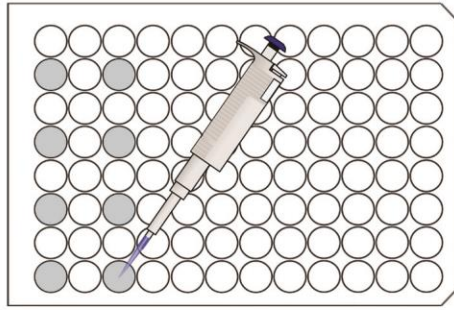
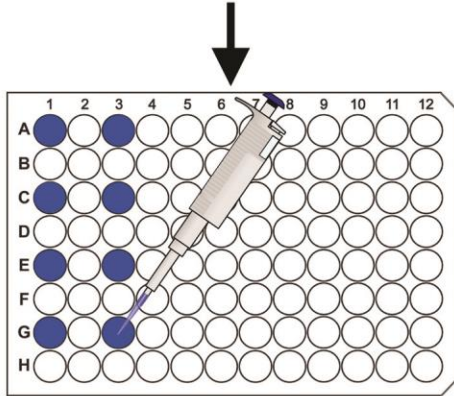


Figure 1: OxiSelect™ Free Hydrogen Sulfide Gas Assay Principle.

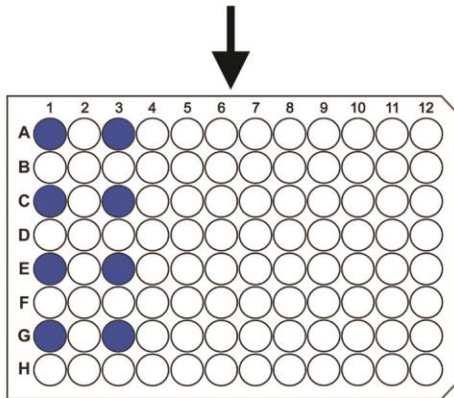
Coat the underside of a 96-well microplate lid with polymer matrix



Pipette sample into the 96-well microplate



Place coated lid on microplate and incubate 1 hour



Transfer coated lid to an empty microplate and read OD 405 nm

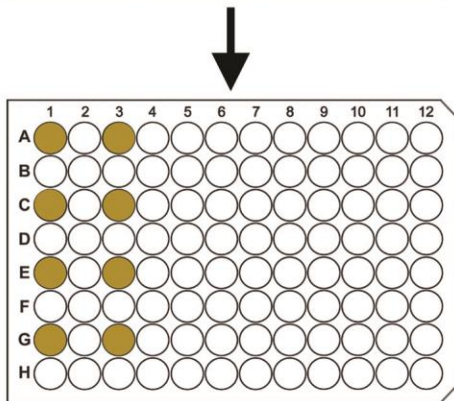


Figure 2: Schematic of Free Hydrogen Sulfide Gas Assay Protocol.

Related Products

1. STA-801: OxiSelect™ In Vitro Nitric Oxide (Nitrite/Nitrate) Assay Kit (Fluorometric)
2. STA-802: OxiSelect™ In Vitro Nitric Oxide (Nitrite/Nitrate) Assay Kit (Colorimetric)
3. XAN-5040: OxiSelect™ Trolox Equivalent Antioxidant Capacity (TEAC) Assay Kit (ABTS)

Kit Components (shipped at room temperature)

1. Sodium Sulfide (Part No. 50841B): One vial containing 100 mg of Na₂S·9H₂O
2. Polymer Matrix (Part No. 50842A): Two 1.5 mL vials.
3. 20X Silver Probe Solution (Part No. 50843A): One 150 µL vial.

Materials Not Supplied

1. 96 well microplate with lid (Costar 96-well plate Cat. No. 3370 recommended)
2. 1 µL to 20 µL adjustable single channel micropipettes with disposable tips
3. 20 µL to 200 µL adjustable single channel micropipettes with disposable tips
4. 200 µL to 1000 µL adjustable single channel micropipettes with disposable tips
5. Spectrophotometric microplate reader capable of reading 405-415 nm
6. Microcentrifuge tubes

Storage

Upon receipt, store the Sodium Sulfide at 4°C. Store all other kit components at room temperature.

Preparation of Reagents

- **Lid Coating Mix:** Transfer the proper amount of Polymer Matrix to a microcentrifuge tube and dilute the 20X Silver Probe Solution 1:20 into the Polymer Matrix. For example add 45 µL of 20X Silver Probe Solution to 855 µL of Polymer Matrix and mix well. This Lid Coating Mix volume is enough for 30 lid spots. The Lid Coating Mix is stable for one day at room temperature in a tightly capped tube.

Note: Prepare only enough for immediate use by scaling the above example proportionally.

- **Sodium Sulfide Positive Control:** Prepare a 10 mM solution of Sodium Sulfide by weighing out 5 milligrams of solid Sodium Sulfide and resuspending the solid at 2.4 mg/mL in deionized water. Then dilute the 10 mM Sodium Sulfide 1:50 into distilled water to a final concentration of 200 µM. For example, for 1 mL of Sodium Sulfide Positive Control, add 20 µL of 10 mM Sodium Sulfide to 980 µL of deionized water. While the 10 mM Sodium Sulfide solution is stable for up to 4 hours in a tightly capped tube, the 200 µM Sodium Sulfide Positive Control should be prepared immediately before use in the assay.

Preparation of Coated Lid

Notes:

- Make sure that the coated circular areas will cover the wells of interest when the lid is properly applied to the microplate.
- To avoid cross contamination of H_2S gas between experimental wells, always arrange the experiment in the 96 well microplate so that a well containing a sample is bordered on every side by a blank well (see Figure 3).

1. Place a 96 well lid upside down on a lab bench, or in a hood.
2. Draw up 30 μ L of Lid Coating Mix (see Preparation of Reagents) with a micropipette and slowly apply about half of the Lid Coating Mix to the center of the circular area that covers one well. Use the pipette tip to evenly spread the Lid Coating Mix throughout the entire circular area. Then apply all but approximately 1-2 μ L of the Lid Coating Mix remaining in the pipette tip to the circular area.

Note: Leaving about 1-2 μ L in the pipette tip ensures that no bubbles will be introduced into the coated circular area. If bubbles are accidentally introduced into the circular area, use the pipette tip to gently aspirate away the bubbles.

3. Allow the inverted lid to dry undisturbed for 1 hour at room temperature. Do not touch or move the lid during this time.

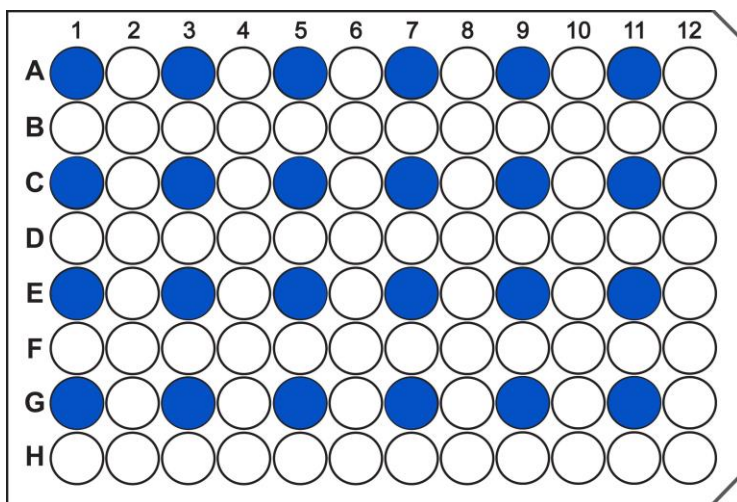


Figure 3. Schematic of a fully used microplate for the Free Hydrogen Sulfide Gas Assay Kit. Wells containing sample are labeled in blue.

Assay Protocol

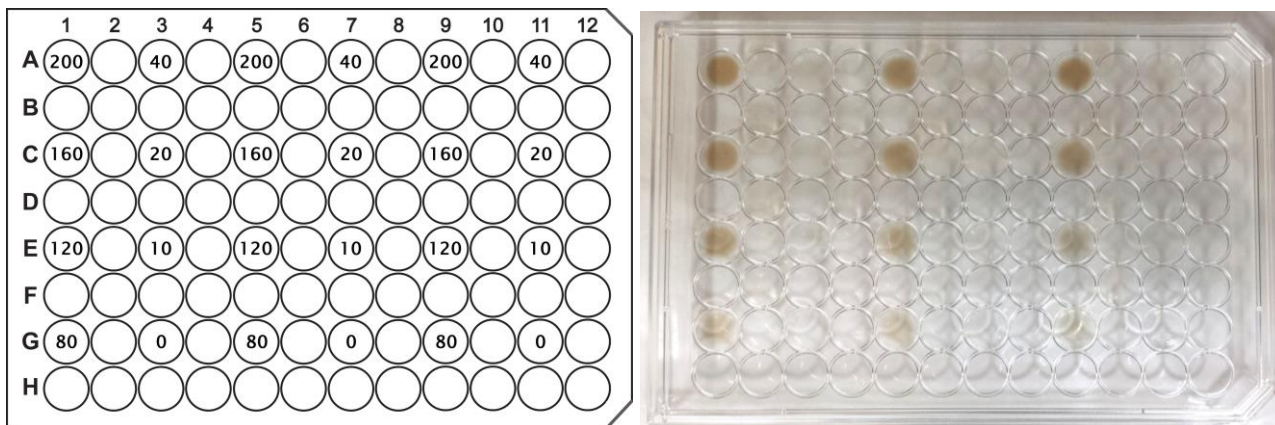
1. Add 300 μ L of unknown sample, Sodium Sulfide positive control, or Negative Control to the wells of a 96 well microplate.
2. Cover the 96 well microplate with the Coated Lid (See Preparation of Coated Lid).
3. Incubate plate at room temperature undisturbed for one hour.
4. Transfer the Coated Lid to a new empty 96 well microplate.

5. Read absorbance of each microwell on a spectrophotometer using 405 nm as the primary wavelength.

Example of Results

The following figures demonstrate typical OxiSelect™ Free Hydrogen Sulfide Gas Assay Kit results. One should use the data below for reference only. This data should not be used to interpret actual results.

A



B

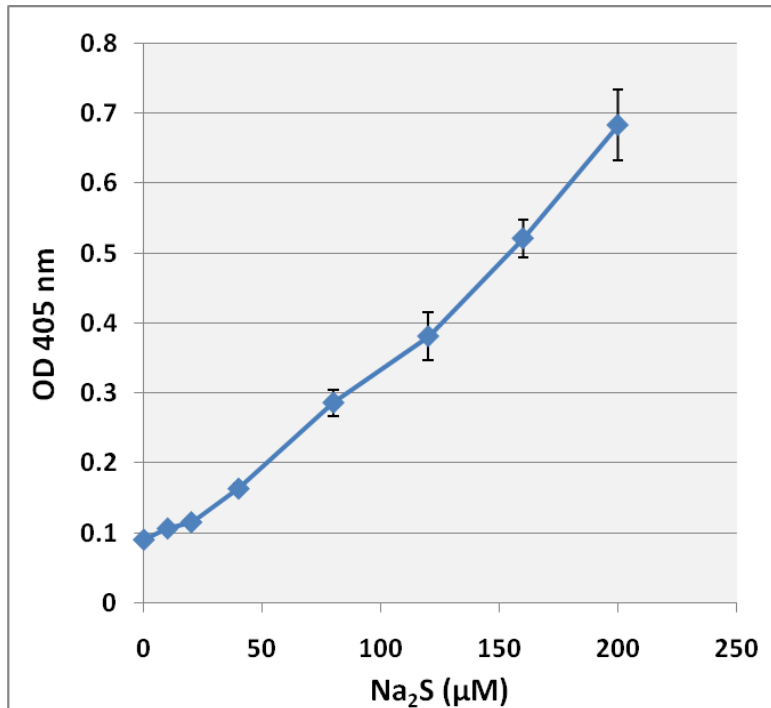


Figure 4. Detection of H₂S from a Hydrogen Sulfide donor. A, Left: Sodium sulfide was prepared at micromolar concentrations indicated and used in the OxiSelect™ Free Hydrogen Sulfide Gas Assay

Kit. **A, Right:** Image of the Coated Lid after one hour of incubation at room temperature according to the Assay Protocol. **B:** Averaged data from (A) plotted with standard deviation.

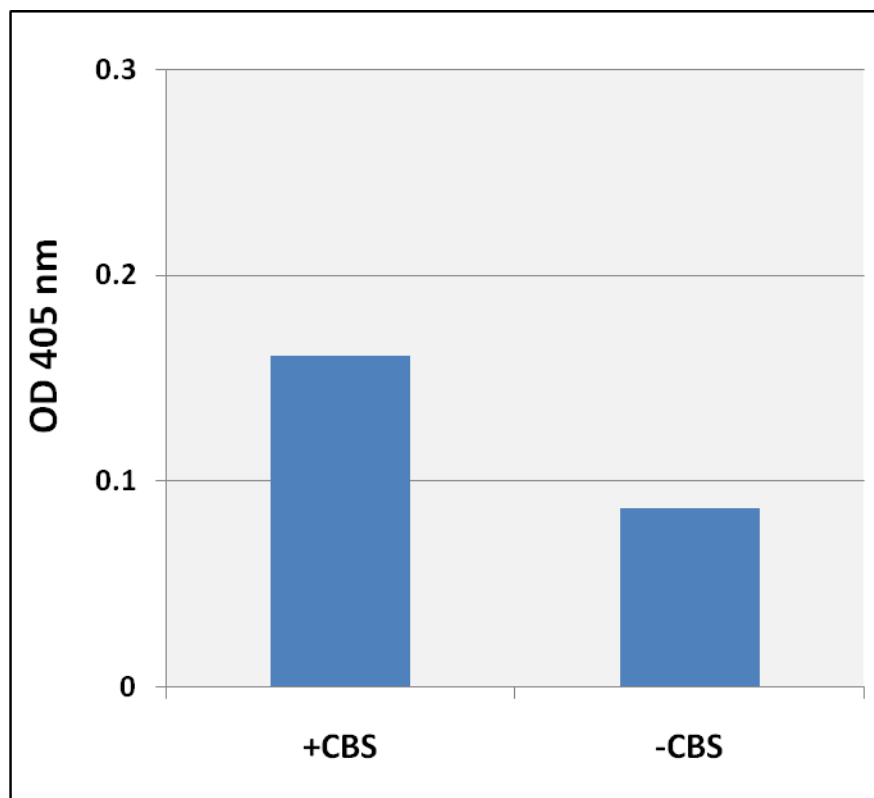


Figure 5. H₂S production by recombinant Cystathionine- β -Synthase (CBS). A Coated Lid was incubated in the absence (-CBS) or presence (+CBS) of 40 μ g of recombinant CBS for 1 hour and then measured with a plate spectrophotometer according to the Assay Protocol.

References

1. Leffler WL, Parfenova H, Jagger JH, and Wang R (2006). Carbon monoxide and hydrogen sulfide: gaseous messengers in cerebrovascular circulation. *J. Appl. Physiol.* **100**, 1065-1076.
2. Greabu M, Totan A, Miricescu D, Radulescu R, Virlan J, and Calenic B. (2016). Hydrogen Sulfide, Oxidative Stress and Periodontal Diseases: A Concise Reivew. *Antioxid. (Basel)*. **5**:3-21.
3. Predmore BL, Lefer DJ, and Gojon G. (2012). Hydrogen sulfide in biochemistry and medicine. *Antioxid. Redox Signal.* **20**: 2118-2120.
4. Kimura H. (2015). Hydrogen sulfide and polysulfides as signaling molecules. *Proc. Jpn. Acad. Ser. B Phys. Biol. Sci.***91**: 131-159.
5. Cheng Y, Ndisang JF, Tang G, Cao K and Wang R. (2004) Sensitive relaxant response of rat mesenteric atery beds to hydrogen sulfide. *Am. J. Physiol.* **287**: H2316-H2323.
6. Yang G, Cao K, and Wang R (2004). Cystathionine gamma lyase overexpression inhibits cell proliferation via a H₂S-dependent modulation of ERK 1/2 phosphorylation and p21^{Cip/WAK-1}. *J. Biol. Chem.* **279**:49199-49205.
7. Kimura H, Shibuya N, and Kimura Y.(2012) Hydrogen sulfide is a signaling molecule and a cytoprotectant. *Antioxid.Redox Signal.***17**:45-57.

Recent Product Citation

Balasubramanian, R. et al. (2022). The Sulfide-Responsive SqrR/BigR Homologous Regulator YgaV of *Escherichia coli* Controls Expression of Anaerobic Respiratory Genes and Antibiotic Tolerance. *Antioxidants (Basel)*. **11**(12):2359. doi: 10.3390/antiox11122359.

Warranty

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS's sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

Contact Information

Cell Biolabs, Inc.
5628 Copley Drive
San Diego, CA 92111
Worldwide: +1 858 271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2017-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.