
Product Manual

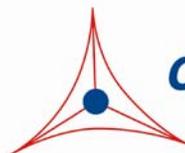
OxiSelect™ Hydrogen Peroxide Assay Kit (Colorimetric)

Catalog Number

STA-343

500 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Oxidative stress is a physiological condition where there is an imbalance between concentrations of reactive oxygen species (ROS) and antioxidants. Research has shown that excessive ROS accumulation will lead to cellular injury, such as damage to DNA, proteins, and lipid membranes. Peroxides, such as hydrogen peroxide (H₂O₂), is one of the most well documented ROS produced under oxidative stress conditions. Hydrogen peroxide is an ROS that is a toxic product of normal aerobic metabolism and pathogenic ROS production involving oxidase and superoxide dismutase reactions. Hydrogen peroxide is poisonous to eukaryotic cells and in high doses can initiate oxidation of DNA, lipids, and proteins, which can lead to mutagenesis and cell death. The cellular damage caused by peroxides have been implicated in the development of many pathological conditions, such as ageing, asthma, arthritis, diabetes, cardiovascular disease, atherosclerosis, Down's Syndrome, and neurodegenerative diseases.

Cell Biolabs' OxiSelect™ Hydrogen Peroxide Assay Kit is a quantitative assay for measuring hydrogen peroxides from both aqueous and lipid solutions. The kit employs a simple HTS-compatible assay for measuring hydrogen peroxide concentrations in biological samples without any need for pretreatment. It can measure aqueous hydrogen peroxides and lipid hydrogen peroxides, and no extraction step is needed for lipid assay format. Absorbance values are proportional to the hydrogen peroxide levels within the samples. The kit has a detection sensitivity limit of 1 μM. Each kit provides sufficient reagents to perform up to 500 assays, including standard curve and unknown samples.

Assay Principle

The OxiSelect™ Hydrogen Peroxide Assay Kit is a quantitative assay for measuring aqueous peroxides and lipid peroxides. For aqueous samples, sorbitol first converts peroxide to a peroxy radical, which oxidizes Fe⁺² into Fe⁺³. For lipid samples, peroxide converts Fe⁺² into Fe⁺³ directly. Then Fe⁺³ reacts with an equal molar amount of xylenol orange in the presence of acid to create a purple product that absorbs maximally between 540-600 nm. The antioxidant BHT is provided to prevent further undesirable chain peroxidation. The peroxide content in unknown samples is determined by comparison with the predetermined H₂O₂ standard curve.

Related Products

1. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
2. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
3. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
4. STA-332: OxiSelect™ MDA ELISA Kit
5. STA-334: OxiSelect™ HNE Adduct ELISA Kit
6. STA-340: OxiSelect™ Superoxide Dismutase Activity Assay
7. STA-341: OxiSelect™ Catalase Activity Assay Kit
8. STA-342: OxiSelect™ ROS Assay Kit
9. STA-345: OxiSelect™ ORAC Activity Assay

Kit Components

1. Xylenol Orange Dye (Part No. 234301): One 1.5 mL amber vial of a 12.5 mM solution.
2. AFS Reagent (Part No. 234302): One 1.5 mL vial of AFS in H₂SO₄.
3. Sorbitol Solution (Part No. 234303): Three 1.5 mL vials of a 4 M solution.
4. Hydrogen Peroxide (Part No. 234102): One 100 µL amber vial of an 8.8 M solution.
5. 1000X BHT Solution (Part No. 234305): One 1.0 mL vial of 5% Butylated hydroxytoluene (BHT) in methanol.
6. TCEP Solution (Part No. 234304): Two 1.5 mL vials of a 10 mM Tris(2-carboxyethyl) phosphine solution in methanol.

Materials Not Supplied

1. Standard 96-well microtiter plates for use in microplate reader.
2. Distilled or deionized water
3. 1X PBS for sample dilutions
4. Methanol for preparing lipid-based assays
5. 10 µL to 1000 µL adjustable single channel micropipettes with disposable tips
6. 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
7. 96-well microtiter plate
8. Multichannel micropipette reservoir
9. Microplate reader capable of reading 560 nm

Storage

Upon receipt, store the BHT and Hydrogen Peroxide at -20°C. Avoid multiple freeze/thaw cycles. Store the remaining kit components at 4°C.

Preparation of Reagents

- **Aqueous Working Reagent:** Prepare the Working Reagent for Aqueous assays by diluting the Xylenol Orange 1:100, the Sorbitol 1:40, and the AFS Reagent 1:100 together with deionized water (eg. For a 25 mL Working Reagent solution: add 0.250 mL of Xylenol Orange, 0.625 mL Sorbitol, and 0.250 mL AFS Reagent together and QS with deionized water to 25 mL. This is enough for 100 assays. Stir or vortex to homogeneity. The Working Reagent is stable for 1 day. Prepare only enough for immediate applications.

Note: AFS Reagent contains acid. Use caution when handling.

- **Lipid Working Reagent:** Prepare the Working Reagent for Lipid assays by diluting the Xylenol Orange 1:100, AFS Reagent 1:100 and BHT 1:1000 together with methanol (90% minimum)(eg. For a 25 mL Working Reagent solution: add 0.250 mL of Xylenol Orange, 0.250 mL AFS Reagent and 50 µL BHT together and QS with methanol to 25 mL). Stir or vortex to

homogeneity. The Working Reagent is stable for 1 day. Prepare only enough for immediate applications.

Sample Preparations

All samples should be assayed immediately or store at -80°C for up to 1-2 months. The assay can be used on cell culture supernatants, serum, plasma, urine, as well as other biological fluids. High levels of interfering substances may cause variations in results. Run proper controls as necessary. Always run a standard curve with samples. Media containing ferrous salts should be avoided, as they will interfere with the assay. Use PBS for dilution and preparation of samples.

Cells: Resuspend cells at $1-2 \times 10^7$ cells/mL in PBS containing 1X BHT. Homogenize or sonicate the cells on ice. Use the whole homogenate in the assay. Cell supernatants can be assayed undiluted.

Cell Culture Supernatants: To remove insoluble particles, spin at 10,000 g for 5 min. The supernatant can be assayed directly or stored at -80°C as necessary.

Serum, Plasma or Urine: To remove insoluble particles, spin at 10,000 g for 5 min. The supernatant can be assayed directly or stored at -80°C as necessary.

Notes:

- *High Hydrogen peroxide ($>200 \mu\text{M}$) can bleach color from the dye, resulting in a lower absorbance value. The working range for this assay is $200 \mu\text{M}$ to $1 \mu\text{M}$. If sample concentrations are unknown, perform several 1:10 serial dilutions of the sample in order to achieve a value within the assay's working range. If absorbance values of diluted samples are the same or higher than the original values, this indicates excessive hydrogen peroxide. Repeat dilutions until the sample values are within range.*
- *If endogenous iron or other transition metals are suspected to be present in the sample, they may cause an increase in the signal. Compensate for the high signal by preparing a Working Reagent Blank (Working Reagent without AFS Reagent). The result will then be obtained by subtracting the sample value obtained with the Working Reagent from the sample value obtained with the Working Reagent Blank.*
- *Several chemicals are known to interfere and should be avoided in sample preparation. These include fructose, sorbitol, sucrose, glucose, formic acid and detergents such as SDS, Tween-20, NP-40 and Triton X-100.*

Preparation of Standard Curve

1. To prepare the Hydrogen Peroxide standards, first perform a 1:1000 dilution of the stock Hydrogen Peroxide in **deionized water (aqueous assay) or methanol (lipid assay)**. Use only enough for immediate applications (eg. Add $5 \mu\text{L}$ of Hydrogen Peroxide to 4.995 mL deionized water). This solution has a concentration of 8.8 mM .
2. Use the $8.8 \text{ mM H}_2\text{O}_2$ solution to prepare standards in the concentration range of $0 \mu\text{M} - 100 \mu\text{M}$ by further diluting in **water (aqueous assay) or methanol (lipid assay)** (see Table 1). H_2O_2 diluted solutions and standards should be prepared fresh. Use the table below as a reference only. The volumes and concentrations of the standard may be adjusted by the user.

Standard Tubes	8.8 mM H ₂ O ₂ Standard (μL)	Deionized Water or MeOH (μL)	H ₂ O ₂ (μM)
1	23	977	200
2	500 of Tube #1	500	100
3	500 of Tube #2	500	50
4	500 of Tube #3	500	25
5	500 of Tube #4	500	12.5
6	500 of Tube #5	500	6.25
7	500 of Tube #6	500	3.125
8	500 of Tube #7	500	1.56
9	500 of Tube #8	500	0.78
10	0	500	0

Table 1. Preparation of H₂O₂ Standards

Assay Protocol

I. For Plasma or Serum Samples

Plasma or serum samples contain both hydrogen peroxide and lipid hydroperoxide. TCEP pretreatment will virtually reduce all the lipid hydroperoxide present.

1. Prepare and mix all reagents thoroughly before use. Prepare the hydrogen peroxide standards simultaneously with the samples so they may be assayed together. Each sample, including unknown and standard, should be assayed in duplicate or triplicate.
2. Add 90 μL of standard or sample to a microcentrifuge tube, followed by 10 μL of the 10 mM TCEP solution.
3. Vortex tubes thoroughly and incubate them for 30 minutes at room temperature.
4. Transfer 25 μL of the TCEP treated standard or sample to fresh microcentrifuge tubes.
5. Add 250 μL of Lipid Working Reagent to each tube. Vortex the contents thoroughly and incubate on a shaker for 30 minutes at room temperature.
6. Centrifuge the sample and standard tubes at 12,000 x g for 5 minutes to remove any precipitate. Transfer the sample supernatant or standard (~275 μL) to a microtiter plate.
7. Read the plate at 540-600 nm (595 nm optimal). The absorbance should be read the same day.
8. Calculate the concentration of hydrogen peroxide within samples by comparing the sample absorbance to the standard curve.

II. For Cells, Cell Culture Supernatants, Urine or other Biological Fluids

1. Prepare and mix all reagents thoroughly before use. Each sample, including unknown and standard, should be assayed in duplicate or triplicate.
2. Add 25 μL of standard or sample to the microtiter plate wells.
3. Add 250 μL of Aqueous or Lipid Working Reagent to each well. Mix the well contents thoroughly and incubate on a shaker for 30 minutes at room temperature.

4. Read the plate at 540-600 nm (595 nm optimal). The absorbance should be read the same day.
5. Calculate the concentration of hydrogen peroxide within samples by comparing the sample absorbance to the standard curve.

Example of Results

The following figures demonstrate typical Peroxide Assay results. One should use the data below for reference only. This data should not be used to interpret actual results.

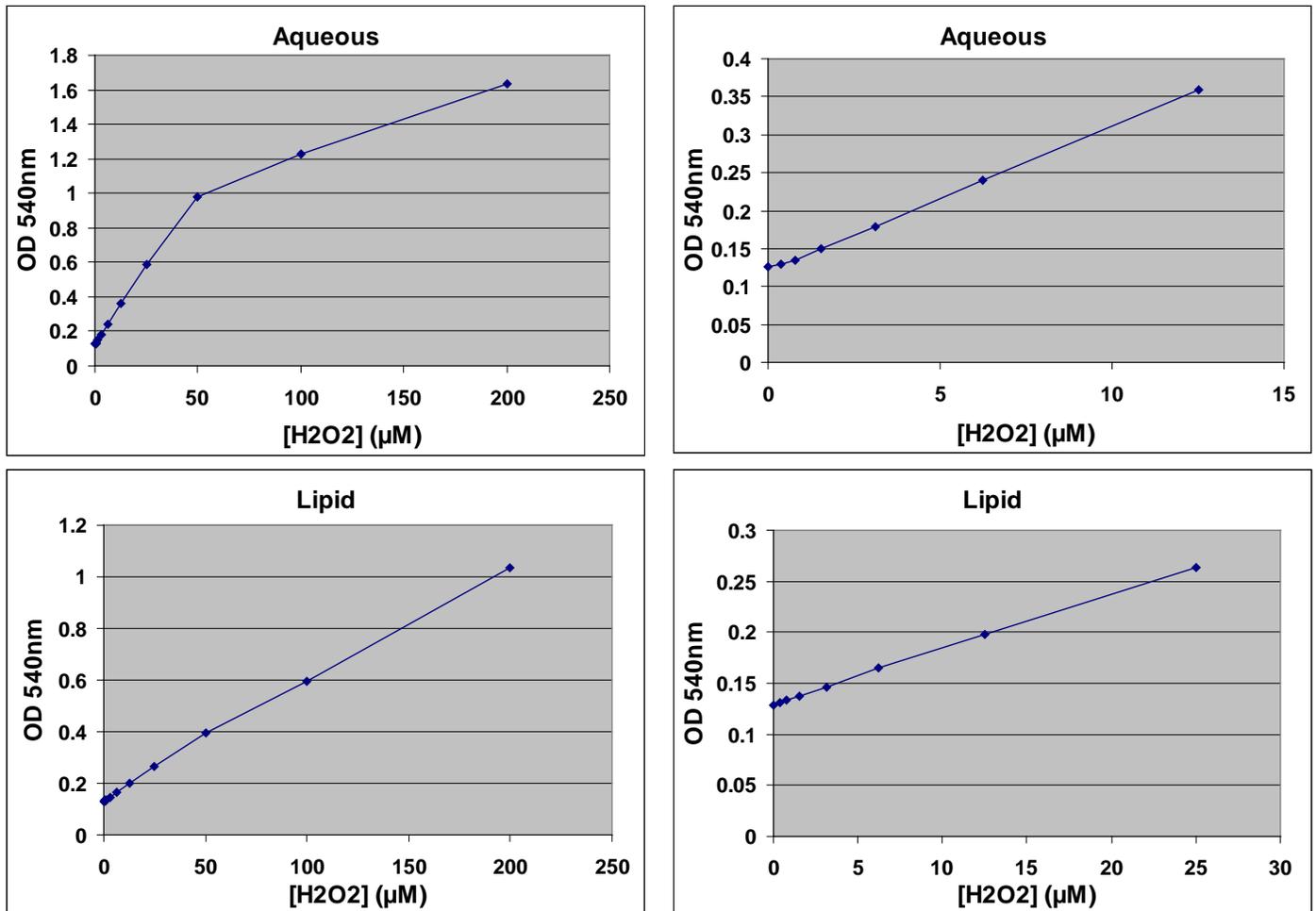


Figure 1. Aqueous and Lipid based H₂O₂ Standard Curves.

References

1. Jiang, Z.Y., Hunt, J.V. and Wolff, S.P. *Anal. Biochem.* (1992) 202: 384-389.

2. Jiang, Z.Y., Woolard, A.C.S. and Wolff, S.P. *FEBS*. (1990) 268(1): 69-71.
3. Jiang, Z.Y., Woolard, A.C.S. and Wolff, S.P. *Lipids*. (1991) 26: 853-856.
4. Nourooz-Zadeh, J. Tajaddini-Sarmadi, J. and Wolff, S.P. *Anal. Biochem.* (1994) 220: 403-409.

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' 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.
7758 Arjons Drive
San Diego, CA 92126
Worldwide: +1 858-271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2010: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.