
Product Manual

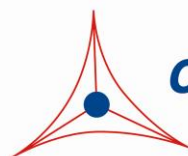
OxiSelect™ Superoxide Dismutase Activity Assay

Catalog Number

STA-340

100 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

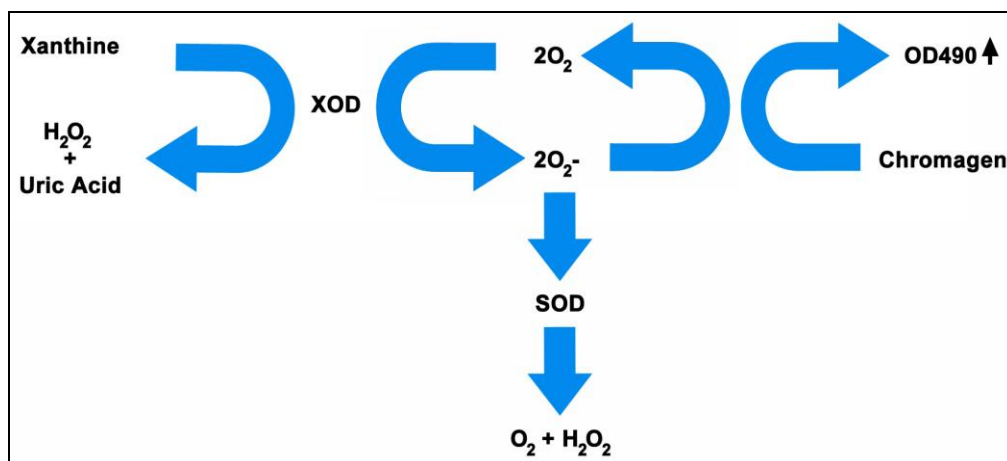
Reactive oxygen species (ROS), such as superoxide (O_2^-) and hydrogen peroxide (H_2O_2), are constantly produced during metabolic processes in all living species. Under normal physiological conditions, cellular ROS generation is counterbalanced by the action of antioxidant enzymes and other redox molecules. However, excessive ROS accumulation will lead to cellular injury, such as damage to DNA, protein, and lipid membrane. Because of their potential harmful effects, excessive ROS must be promptly eliminated from the cells by a variety of antioxidant defense mechanisms. Superoxide dismutase (SOD), which catalyzes the dismutation of the superoxide anion into hydrogen peroxide and molecular oxygen, is one of the most important antioxidative enzymes. SOD enzymes are classified into three groups: cytosolic Cu/Zn-SOD, mitochondrial Mn-SOD, and extracellular Ec-SOD.

Our OxiSelect™ Superoxide Dismutase Activity Assay uses a xanthine/xanthine oxidase (XOD) system to generate superoxide anions. The included chromagen produces a water-soluble formazan dye upon reduction by superoxide anions. The activity of SOD is determined as the inhibition of chromagen reduction (See Figure 1).

The OxiSelect™ Superoxide Dismutase Activity Assay is a fast and reliable kit for the measurement of SOD activity from cell lysate, plasma, serum, tissue homogenates. Each kit provides sufficient reagents to perform up to 100 assays, including blanks, SOD standards and unknown protein samples.

Assay Principle

Superoxide anions (O_2^-) are generated by a Xanthine/Xanthine Oxidase (XOD) system, and then detected with a Chromagen Solution. However, in the presence of SOD, these superoxide anion concentrations are reduced, yielding less colorimetric signal.



Related Products

1. STA-305: OxiSelect™ Nitrotyrosine ELISA Kit
2. STA-310: OxiSelect™ Protein Carbonyl ELISA Kit
3. STA-312: OxiSelect™ Total Glutathione (GSSG/GSH) Assay Kit
4. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
5. STA-341: OxiSelect™ Catalase Activity Assay Kit

Kit Components

1. **SOD Standard** (Part No. 234001): One 100 μ L vial provided at 5 Units/ μ L. Unit Definition: One unit will inhibit the rate of reduction of cytochrome c by 50% in a coupled system, using xanthine and xanthine oxidase, at pH 7.8 at 25°C in a 3.0 ml reaction volume.
2. **Xanthine Solution** (Part No. 234002): One 550 μ L vial.
3. **Xanthine Oxidase Solution, 150X** (Part No. 234003): One 50 μ L vial.
4. **Chromagen Solution** (Part No. 234004): One 550 μ L amber vial.
5. **SOD Assay Buffer, 10X** (Part No. 234005): Two 1.5 mL vials.

Materials Not Supplied

1. 96-well microtiter plate
2. Microplate reader capable of reading at 490 nm

Storage

Store kit components at -20°C. Avoid multiple freeze/thaws by aliquoting. The Chromagen Solution is light sensitive and should be maintained in amber tubes.

Preparation of Reagents

- 1X SOD Assay Buffer: Dilute one vial of 10X SOD Assay Buffer to 1X with deionized water. Mix to homogeneity. Keep the second vial of 10X SOD Assay Buffer undiluted.
- 1X Xanthine Oxidase Solution: Just prior to use, dilute the 150X Xanthine Oxidase Solution to 1X with 1X SOD Assay Buffer. Mix to homogeneity.

Special Precautions

Avoid the use of reducing agents, such as DTT, in the assay due to interference with the Chromagen Solution.

Preparation of Samples

- Suspension Cells: Centrifuge 3-6 x 10⁶ cells at 700 x g for 2 minutes and discard supernatant. Wash cell pellet once with ice-cold PBS, centrifuge, and discard the supernatant. Resuspend cell pellet in 0.5 mL of cold 1X Lysis Buffer (10 mM Tris, pH 7.5, 150 mM NaCl, 0.1 mM EDTA). Lyse cells with sonication or homogenization. Centrifuge at 12000 x g for 10 minutes and collect the cell lysate supernatant.
- Adherent Cells: Wash 1-5 x 10⁶ cells once with 10 mL ice-cold PBS per 100 mm dish. Harvest cells with a cell scraper in 1 mL of cold 1X Lysis Buffer (10 mM Tris, pH 7.5, 150 mM NaCl, 0.1 mM EDTA). Lyse cells with sonication or homogenization. Centrifuge at 12000 x g for 10 minutes and collect the cell lysate supernatant.
- Tissue Lysates: Homogenize tissue sample in 5-10 mL of cold 1X Lysis Buffer (10 mM Tris, pH 7.5, 150 mM NaCl, 0.1 mM EDTA) per gram tissue. Lyse cells with sonication or homogenization. Centrifuge at 12000 x g for 10 minutes and collect the tissue lysate supernatant.
- Plasma: Collect blood with an anticoagulant such as heparin, citrate or EDTA and mix by inversion. Centrifuge the blood at 1000 x g at 4°C for 10 minutes. Collect plasma supernatant

without disturbing the white buffy layer. Sample should be tested immediately or frozen at -80°C for storage.

- Serum: Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage.

Assay Protocol

1. Prepare samples including a blank in a 96-well microtiter plate according to the below table. Allow pre-incubation time if inhibitor is used.

Component	Blank	Sample
SOD Sample	0 μ L	X μ L
Inhibitor (optional)	0 μ L	Y μ L
Xanthine Solution	5 μ L	5 μ L
Chromagen Solution	5 μ L	5 μ L
10X SOD Assay Buffer	10 μ L	10 μ L
DI Water	70 μ L	70-(X+Y) μ L
Total	90 μL	90 μL

2. Finally, add 10 μ L of pre-diluted 1X Xanthine Oxidase Solution (see Preparation of Reagents) to each well. Mix well and incubate for 1 hour at 37°C.
3. Read absorbance at 490 nm on a microplate reader.

Preparation of SOD Standards (Optional)

1. Thaw SOD Standard at 4°C.
2. Freshly prepare a dilution series (1:4 is suggested) of SOD Standard in the concentration range of 5 Units/ μ L – 1.2 mU/ μ L by diluting the SOD Standard in 1X Assay Buffer (see Preparation of Reagents).

Standard Tubes	SOD Standard (μL)	1X Assay Buffer (μL)	SOD (U/ μL)
1	35	0	5
2	10 of Tube #1	30	1.25
3	10 of Tube #2	30	0.312
4	10 of Tube #3	30	0.078
5	10 of Tube #4	30	0.0195
6	10 of Tube #5	30	0.0048
7	10 of Tube #6	30	0.0012
8	0	30	0

Table 1. Suggested preparation of SOD standards

3. Transfer 10 μ L of each dilution to a 96-well microtiter plate, including a 1X Assay Buffer blank.
4. Prepare the following master mixture, adjusting for the required number of wells.

Component	Volume per well
Xanthine Solution	5 μ L
Chromagen Solution	5 μ L
10X SOD Assay Buffer	10 μ L
DI Water	60 μ L
Total	80 μL

- Transfer 80 μ L of the above master mixture to each well.
- Finally, add 10 μ L of pre-diluted 1X Xanthine Oxidase Solution (see Preparation of Reagents) to each well. Mix well and incubate for 1 hour at 37°C.
- Read absorbance at 490 nm on a microplate reader.

Example of Results

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

$$\text{SOD Activity (inhibition \%)} = (\text{OD}_{\text{blank}} - \text{OD}_{\text{sample}}) / (\text{OD}_{\text{blank}}) \times 100$$

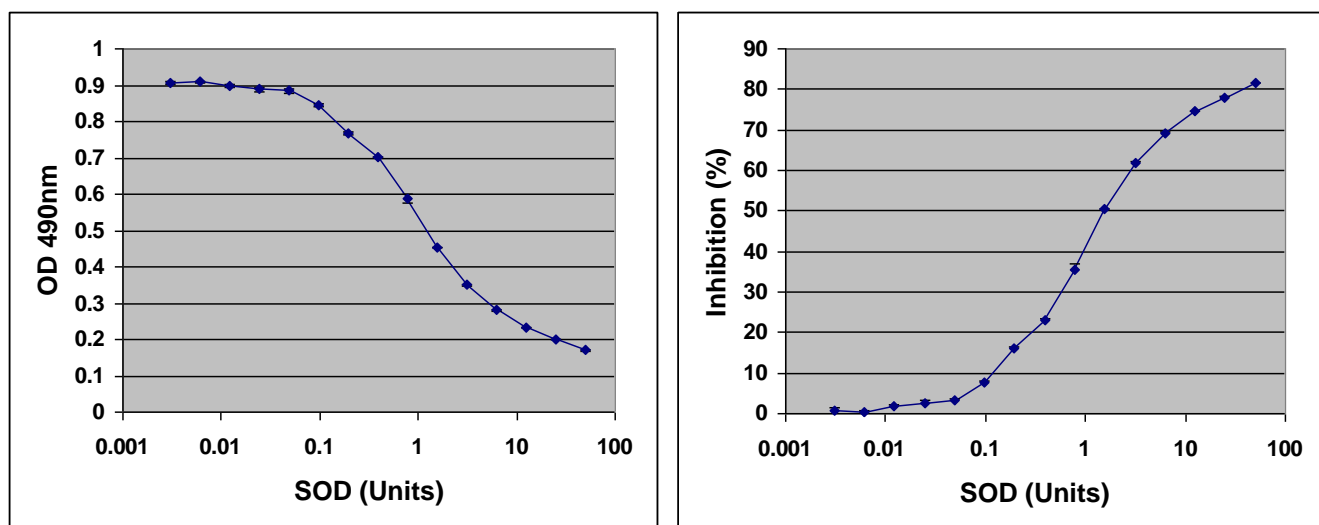


Figure 1: SOD Activity Assay Standard Curve. Left: SOD activity as a function of optical density (OD). **Right:** SOD activity as a function of inhibition percentage.

References

- Lepock, J. R., Arnold, L. D., Torrie, B. H., Andrews, B., and Kruuv, J. (1985) *Arch. Biochem. Biophys.* **241**, 243–251.
- Lepock, J. R., Frey, H. E., and Hallewell, R. A. (1990) *J. Biol. Chem.* **265**, 21612–21618.
- Valentine, J. S., and Hart, P. J. (2003) *Proc. Natl. Acad. Sci. U. S. A.* **100**, 3617-3622.
- Zelko, I. N., Mariani, T. J., and Folz, R. J. (2002) *Free Radic. Biol. Med.* **33**, 337-349.

Recent Product Citations

1. Saha, T.K. et al. (2023). Immuno-physiological effects of dietary reishi mushroom powder as a source of beta-glucan on Rohu, *Labeo rohita* challenged with *Aeromonas veronii*. *Sci Rep.* **13**(1):14652. doi: 10.1038/s41598-023-41557-9.
2. Elrayess, A.R. et al. (2023). Biochemical, Histological and Ultrastructural Studies on the Effect of Citric acid Supplementation on Aflatoxins-intoxicated Japanese Quail. *J. Adv. Vet.* **13**(6):958-964.
3. Hadžić, Z. et al. (2023). Oxidative Stress And C-Reactive Protein As Salivary Biomarkers In Smokers With Periodontitis Stage III And IV After Non-Surgical Periodontal Therapy (A Pilot Study). *Acta Medica Mediterranea.* **39**:947-953. doi: 10.19193/0393-6384_2023_4_131.
4. Sarkar, S. et al. (2023). Microplastic Pollution: Chemical Characterization and Impact on Wildlife. *Int J Environ Res Public Health.* **20**(3):1745. doi: 10.3390/ijerph20031745.
5. Rajab, B.S. et al. (2022). Antioxidative and Anti-Inflammatory Protective Effects of β -Caryophyllene against Amikacin-Induced Nephrotoxicity in Rat by Regulating the Nrf2/AMPK/AKT and NF- κ B/TGF- β /KIM-1 Molecular Pathways. *Oxid Med Cell Longev.* doi: 10.1155/2022/4212331.
6. Emam, K.K. et al. (2022). Assessment of Wheat Germ Oil Role in the Prevention of Induced Breast Cancer in Rats. *ACS Omega.* **7**(16):13942-13952. doi: 10.1021/acsomega.2c00434.
7. Carmona-Rodríguez, L. et al. (2022). Superoxide Dismutase-3 Downregulates Laminin α 5 Expression in Tumor Endothelial Cells via the Inhibition of Nuclear Factor Kappa B Signaling. *Cancers (Basel).* **14**(5):1226. doi: 10.3390/cancers14051226.
8. Refaat, B. & El-Boshy, M. (2022). Protective antioxidative and anti-inflammatory actions of β -caryophyllene against sulfasalazine-induced nephrotoxicity in rat. *Exp Biol Med (Maywood).* doi: 10.1177/15353702211073804.
9. Al-Saleh, I. et al. (2022). Essential metals, vitamins and antioxidant enzyme activities in COVID-19 patients and their potential associations with the disease severity. *Biometals.* doi: 10.1007/s10534-021-00355-4.
10. Zhang, Q. et al. (2021). Micelle silymarin supplementation to sows' diet from day 109 of gestation to entire lactation period enhances reproductive performance, and affects serum hormones and metabolites. *J Anim Sci.* doi: 10.1093/jas/skab354.
11. Nagata, M. et al. (2021). Protective Effects of the Alga Fucoïdan Against Amyloid- β -Induced Neurotoxicity in SH-SY5Y Cells. *BPB Reports.* **4**: 206-213.
12. Hossain, Z. et al. (2021). Heavy metal toxicity in Buriganga river alters the immunology of Nile tilapia (*Oreochromis niloticus* L). *Heliyon.* **7**(11): e08285. doi: 10.1016/j.heliyon.2021.e08285.
13. Widiapradja, A. et al. (2021). Replacement of Lost Substance P Reduces Fibrosis in the Diabetic Heart by Preventing Adverse Fibroblast and Macrophage Phenotype Changes. *Cells.* **10**(10):2659. doi: 10.3390/cells10102659.
14. Ammar, I.M.M. & Salem, M.A.A. et al. (2021). Amelioration of polycystic ovary syndrome-related disorders by supplementation of thymoquinone and metformin. *Middle East Fertil Soc J.* doi: 10.1186/s43043-021-00076-1.
15. Kimura, A.M. et al. (2021). Myricetin prevents high molecular weight A β 1-42 oligomer-induced neurotoxicity through antioxidant effects in cell membranes and mitochondria. *Free Radic Biol Med.* **171**:232-244. doi: 10.1016/j.freeradbiomed.2021.05.019.
16. Sah, D.K. et al. (2021). Sphingosine kinase inhibitor, SKI-II confers protection against the ionizing radiation by maintaining redox homeostasis most likely through Nrf2 signaling. *Life Sci.* doi: 10.1016/j.lfs.2021.119543.

17. Almarhoun, M. et al. (2021). Overexpression of STARD3 attenuates oxidized LDL-induced oxidative stress and inflammation in retinal pigment epithelial cells. *Biochim Biophys Acta Mol Cell Biol Lipids*. **1866**(7):158927. doi: 10.1016/j.bbalip.2021.158927.
18. Zhang, Y. et al. (2021). Neuroprotective effect of the somatostatin receptor 5 agonist L-817,818 on retinal ganglion cells in experimental glaucoma. *Exp Eye Res*. **204**:108449. doi: 10.1016/j.exer.2021.108449.
19. Nour, M.A. et al. (2021). Productive performance, fertility and hatchability, blood indices and gut microbial load in laying quails as affected by two types of probiotic bacteria. *Saudi J Biol Sci*. doi: 10.1016/j.sjbs.2021.07.030.
20. Mukem, S. et al. (2021). Mito-Tempo suppresses autophagic flux via the PI3K/Akt/mTOR signaling pathway in neuroblastoma SH-SY5Y cells. *Heliyon*. doi: 10.1016/j.heliyon. 2021.e07310.
21. de Los Santos-Jiménez, J. et al. (2021). Glutaminase isoforms expression switches microRNA levels and oxidative status in glioblastoma cells. *J Biomed Sci*. **28**(1):14. doi: 10.1186/s12929-021-00712-y.
22. Dong, S. et al. (2021). Leukemia inhibitory factor protects photoreceptor cone cells against oxidative damage through activating JAK/STAT3 signaling. *Ann Transl Med*. **9**(2):152. doi: 10.21037/atm-20-8040.
23. Ra, K. et al. (2021). Comparison of Anti-Oxidative Effect of Human Adipose- and Amniotic Membrane-Derived Mesenchymal Stem Cell Conditioned Medium on Mouse Preimplantation Embryo Development. *Antioxidants (Basel)*. **10**(2):268. doi: 10.3390/antiox10020268.
24. Tsvetkova, A.S. et al. (2020). Melatonin Prevents Early but Not Delayed Ventricular Fibrillation in the Experimental Porcine Model of Acute Ischemia. *Int J Mol Sci*. **22**(1): E328. doi: 10.3390/ijms22010328.
25. Ibnu Rasid, E.N. et al. (2020). Effect of Dioscorea hispida var. Daemona (Roxb) Prain & Burkill on Oxidative Stress and DNA Damage in the Liver of Pregnant Rats. *Int J Biomed Sci*. **16**(3).
26. Alfarisi, H.A.H. et al. (2020). Hepatoprotective Effects of a Novel Trihoney against Nonalcoholic Fatty Liver Disease: A Comparative Study with Atorvastatin. *The Scientific World Journal*. doi: 10.1155/2020/4503253.
27. Hwang, D.K. et al. (2020). Changes in the Systemic Expression of Sirtuin-1 and Oxidative Stress after Intravitreal Anti-Vascular Endothelial Growth Factor in Patients with Retinal Vein Occlusion. *Biomolecules*. **10**(10):1414. doi: 10.3390/biom10101414.
28. Li, T. et al. (2020). Nesfatin-1 Promotes Proliferation, Migration and Invasion of HTR-8/SVneo Trophoblast Cells and Inhibits Oxidative Stress via Activation of PI3K/AKT/mTOR and AKT/GSK3 β Pathway. *Reprod Sci*. doi: 10.1007/s43032-020-00324-1.
29. Kim, J.H. et al. (2020). Comparison of toxic effects of dietary organic or inorganic selenium and prediction of selenium intake and tissue selenium concentrations in broiler chickens using feather selenium concentrations. *Poult Sci*. doi: 10.1016/j.psj.2020.08.061.
30. Chida, J. et al. (2020). Prion protein signaling induces M2 macrophage polarization and protects from lethal influenza infection in mice. *PLoS Pathog*. **16**(8): e1008823. doi: 10.1371/journal.ppat.1008823.

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.
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

©2008-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.