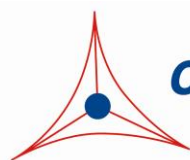

Product Manual

OxiSelect™ *In Vitro* ROS/RNS Assay Kit (Green Fluorescence)

Catalog Number

STA-347	96 assays
STA-347-5	5 x 96 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) and reactive nitrogen species (RNS) are well-established molecules responsible for the deleterious effects of oxidative stress. Accumulation of free radicals coupled with an increase in oxidative stress has been implicated in the pathogenesis of several disease states. The role of oxidative stress in vascular diseases, diabetes, renal ischemia, atherosclerosis, pulmonary pathological states, inflammatory diseases, cancer, as well as ageing has been well established. Free radicals and other reactive species are constantly generated *in vivo* and cause oxidative damage to biomolecules, a process held in check by the existence of multiple antioxidant and repair systems as well as the replacement of damaged nucleic acids, proteins and lipids. Measuring the effect of antioxidant therapies and ROS/RNS activity is crucial to suppressing or treating oxidative stress inducers.

The OxiSelect™ *In Vitro* ROS/RNS Assay Kit is an assay for measuring the total free radical presence of a sample. The assay employs a proprietary quenched fluorogenic probe, dichlorodihydrofluorescein DiOxyQ (DCFH-DiOxyQ), which is a specific ROS/RNS probe that is based on similar chemistry to the popular 2', 7'-dichlorodihydrofluorescein diacetate. The DCFH-DiOxyQ probe is first primed with a quench removal reagent, and subsequently stabilized in the highly reactive DCFH form. In this reactive state, ROS and RNS species can react with DCFH, which is rapidly oxidized to the highly fluorescent 2', 7'-dichlorodihydrofluorescein (DCF) (Figure 1). Fluorescence intensity is proportional to the total ROS/RNS levels within the sample. The DCFH-DiOxyQ probe can react with hydrogen peroxide (H₂O₂), peroxy radical (ROO·), nitric oxide (NO), and peroxynitrite anion (ONOO⁻). These free radical molecules are representative of both ROS and RNS, thus allowing for measurement of the total free radical population within a sample. OxiSelect™ *In Vitro* ROS/RNS Assay Kit can also be used to evaluate antioxidant's effect on free radicals. The kit has a detection sensitivity limit of 10 pM for DCF and 40 nM for H₂O₂ respectively. Each kit provides sufficient reagents to perform up to 96 assays, including standard curve and unknown samples.

Assay Principle

The OxiSelect™ *In Vitro* ROS/RNS Assay Kit is an *in vitro* assay for measuring total ROS/RNS free radical activity. Unknown ROS or RNS samples or standards are added to the wells with a catalyst that helps accelerate the oxidative reaction. After a brief incubation, the prepared DCFH probe is added to all wells and the oxidation reaction is allowed to proceed (Figure 1). Samples are measured fluorometrically against a hydrogen peroxide or DCF standard. The assay is performed in a 96-well fluorescence plate format that can be read on a standard fluorescence plate reader. The free radical content in unknown samples is determined by comparison with the predetermined DCF or hydrogen peroxide standard curve.

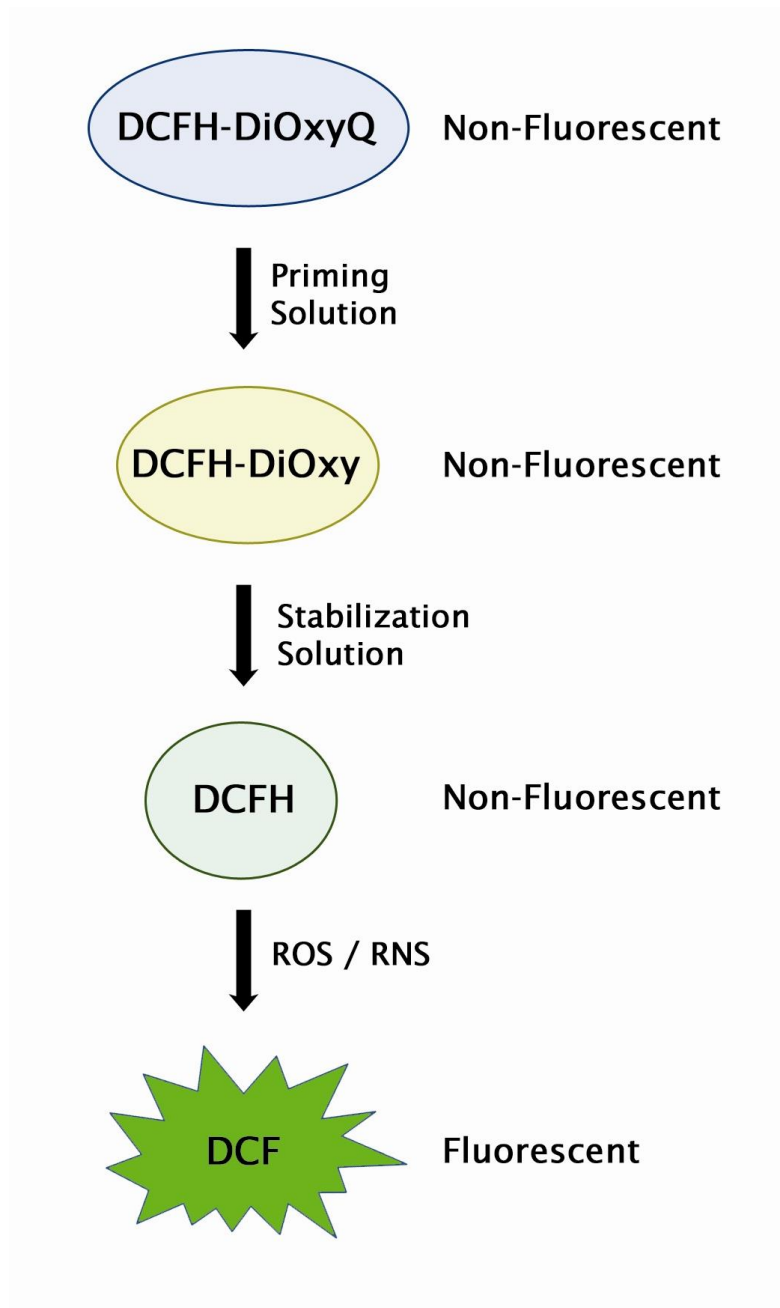


Figure 1. Mechanism of *In Vitro* ROS/RNS Assay.

Related Products

1. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
2. STA-340: OxiSelect™ Superoxide Dismutase Activity Assay
3. STA-341: OxiSelect™ Catalase Activity Assay Kit
4. STA-342: OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)

5. STA-832: OxiSelect™ MDA Adduct Competitive ELISA Kit
6. STA-838: OxiSelect™ HNE Adduct Competitive ELISA Kit

Kit Components

1. Priming Reagent (Part No. 234701): One 250 µL tube of solution.
2. Stabilization Solution (10X) (Part No. 234702): One 1.5 mL tube of solution.
3. Catalyst (250X) (Part No. 234703): One 20 µL tube of solution.
4. DCF-DiOxyQ (Part No. 234704): One 50 µL amber tube of solution in methanol.
5. DCF Standard (Part No. 234202): One 100 µL amber tube of a 1 mM solution in DMSO.
6. Hydrogen Peroxide (Part No. 234102): One 100 µL amber tube of an 8.821 M solution.

Materials Not Supplied

1. 10 µL to 1000 µL adjustable single channel micropipettes with disposable tips
2. 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
3. Multichannel micropipette reservoir
4. Phosphate Buffered Saline for sample preparations and dilutions
5. 96-well black or fluorescence microtiter plate
6. Fluorescent microplate reader capable of reading 480 nm (excitation) and 530 nm (emission)

Storage

Upon receipt, store the DCF-DiOxyQ and DCF Standard at -20°C. Avoid multiple freeze/thaw cycles. Store all other components at 4°C.

Preparation of Reagents

- 1X Stabilization Solution: Dilute the 10X Stabilization Solution 1:10 by adding 1.5 mL of solution to 13.5 mL of deionized water. Stir or vortex to homogeneity. Store the solution at 4°C.
- 1X Catalyst: Prior to use, dilute the 250X Catalyst 1:250 in PBS. Vortex thoroughly. Prepare only enough for immediate applications (eg. add 10 µL of Catalyst to 2.49 mL PBS for 50 wells).
- DCFH Solution: Prepare only enough DCFH Solution for immediate applications in an amber tube or aluminum foil covered tube. Prepare DCFH Solution by diluting the stock solution of DCF-DiOxyQ 1:5 with Priming Reagent (eg. for 50 assays, add 25 µL DCF-DiOxyQ to 100 µL Priming Reagent). Vortex to homogeneity. Incubate the solution for 30 minutes at room temperature. Next, dilute the reaction 1:40 with 1X Stabilization Solution (eg. for 50 assays, add 125 µL DCF-DiOxyQ/ Priming Reagent reaction to 4.875 mL of Stabilization Solution). Vortex to homogeneity. Protect the solution from light. This solution is now stable in the DCFH form and ready to use. The solution may be stored at -20°C for up to one week when protected from light.

Note: Due to light-induced auto-oxidation, the stock DCF-DiOxyQ solution and all subsequent DCF-DiOxy and DCFH solutions must be protected from light.

Preparation of Samples

All samples should be assayed immediately or stored at -80°C for up to 1-2 months. The assay may be used on cell or tissue lysates, cell culture supernatants, serum, plasma, urine, and other biological fluids. Always run a standard curve with samples. Use PBS for dilution and preparation of samples.

Some common detergents and denaturants have been tested for compatibility in the assay (below table). Dilution of samples, and interfering substances, may be necessary for assay compatibility.

Substance	Compatible Concentration
Triton X-100	≤1%
NP-40	≤1%
SDS	≤0.1%
Deoxycholate	≤1%
Tween-20	≤0.1%
EDTA	≤10 mM
EGTA	≤10 mM
Glycerol	≤10%

Table 1. Substance Compatibility Table

- Cells or Tissues: Resuspend cells at $1-2 \times 10^7$ cells/mL or tissues at 10-50 mg/mL in PBS. Homogenize or sonicate on ice. To remove insoluble particles, spin at 10,000 g for 5 min. The homogenate can be assayed directly or stored at -80°C as necessary.
- Serum, Plasma, Urine or 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.

Preparation of the DCF Standard Curve

1. Prepare a 1:10 dilution series of DCF standards in the concentration range of 0 μ M – 10 μ M by diluting the 1mM DCF stock in 1X PBS (see Table 2).

Standard Tubes	DCF Standard (μ L)	PBS (μ L)	DCF (nM)
1	10	990	10,000
2	100 of Tube #1	900	1000
3	100 of Tube #2	900	100
4	100 of Tube #3	900	10
5	100 of Tube #4	900	1
6	0	1000	0

Table 2. Preparation of DCF Standards

- Transfer 200 μL of each DCF standard to a 96-well plate suitable for fluorescence measurement.
- Read the relative fluorescence with a fluorescence plate reader at 480 nm excitation / 530 nm emission.

Preparation of the H_2O_2 Standard Curve

- To prepare the Hydrogen Peroxide standards, first perform a 1:4400 dilution of the stock Hydrogen Peroxide in deionized water. Use only enough for immediate applications (eg. Add 5 μL of Hydrogen Peroxide to 22 mL deionized water). This solution has a concentration of 2 mM.
- Use the 2 mM H_2O_2 solution to prepare standards in the concentration range of 0 μM – 20 μM by further diluting in PBS (see Table 3). H_2O_2 diluted solutions and standards should be prepared fresh. Use the table below as a reference guide only. The volumes and concentrations of the standard may be adjusted by the user.

Standard Tubes	2 mM H_2O_2 Standard (μL)	PBS (μL)	H_2O_2 (μM)
1	10	990	20
2	500 of Tube #1	500	10
3	500 of Tube #2	500	5
4	500 of Tube #3	500	2.5
5	500 of Tube #4	500	1.25
6	500 of Tube #5	500	0.625
7	500 of Tube #6	500	0.313
8	500 of Tube #7	500	0.156
9	500 of Tube #8	500	0.078
10	500 of Tube #9	500	0.039
11	0	1000	0

Table 3. Preparation of H_2O_2 Standards

Assay Protocol

- Prepare and mix all reagents thoroughly before use. Each sample, including unknown(s) and standard(s), should be assayed in duplicate or triplicate.
- Add 50 μL of unknown sample or hydrogen peroxide standard to wells of a 96-well plate suitable for fluorescence measurement.
- Add 50 μL of Catalyst to each well. Mix well and incubate 5 minutes at room temperature.
- Add 100 μL of DCFH solution to each well. Cover the plate reaction wells to protect them from light and incubate at room temperature for 15-45 minutes.
- Read the fluorescence with a fluorescence plate reader at 480 nm excitation / 530 nm emission.

Example of Results

The following figures demonstrate typical Free Radical ROS/RNS Assay results. Fluorescence measurement was performed on SpectraMax Gemini XS Fluorometer (Molecular Devices) with a

485/538 nm filter set and 530 nm cutoff. One should use the data below for reference only. This data should not be used to interpret actual results.

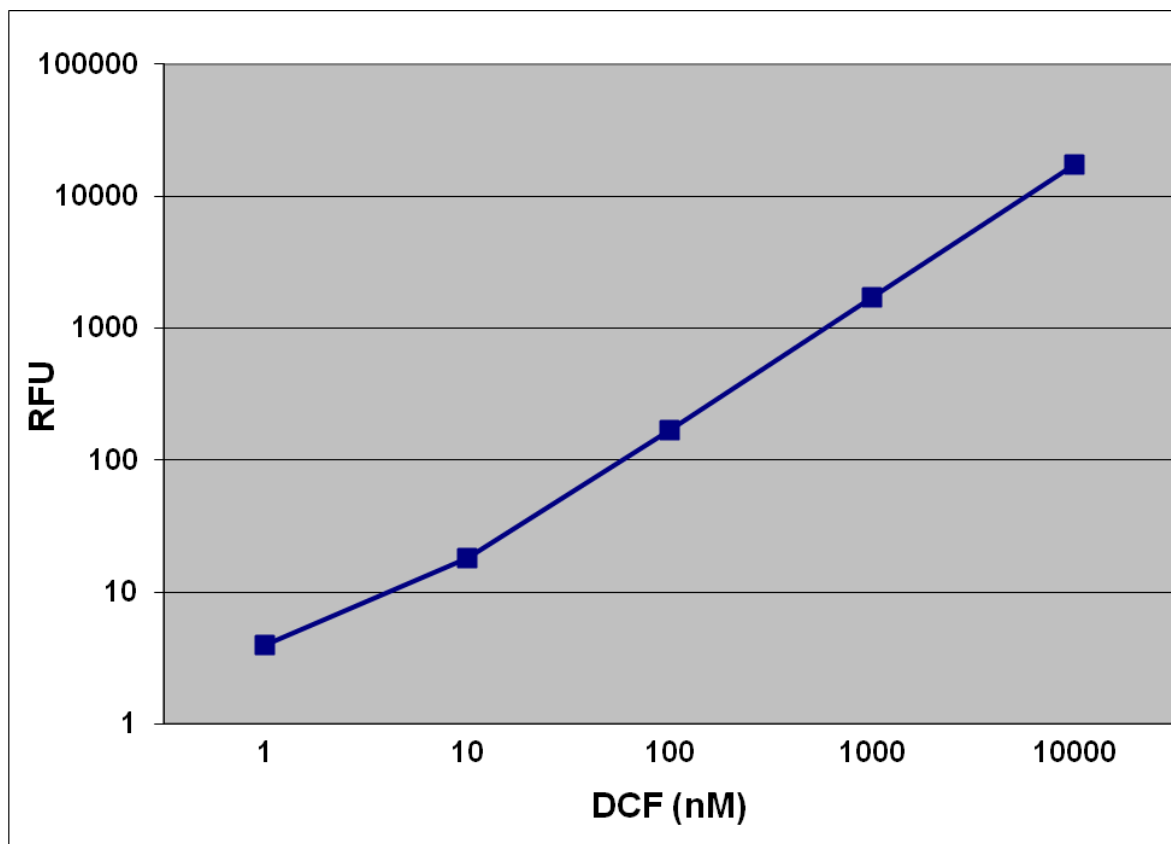


Figure 2. DCF Standard Curve.

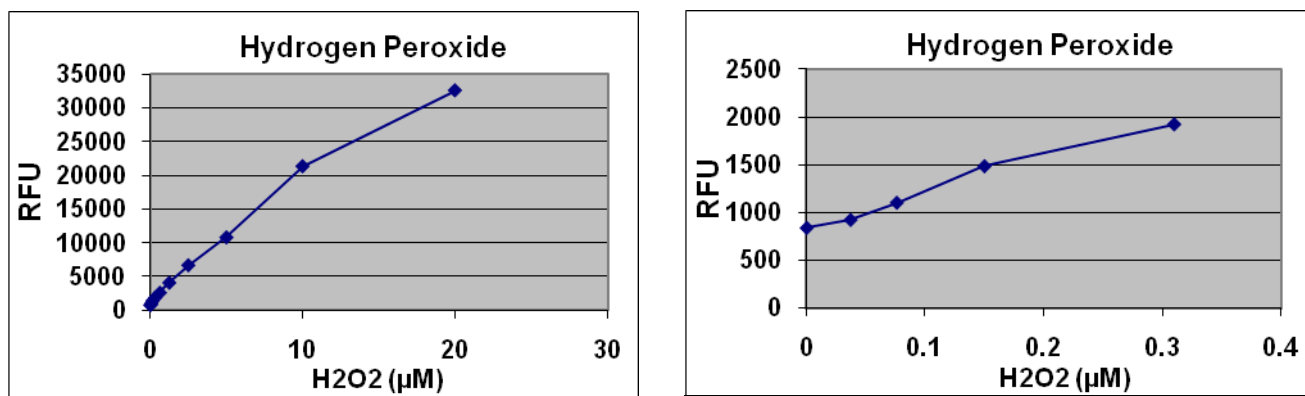


Figure 3. Hydrogen Peroxide Standard Curve.

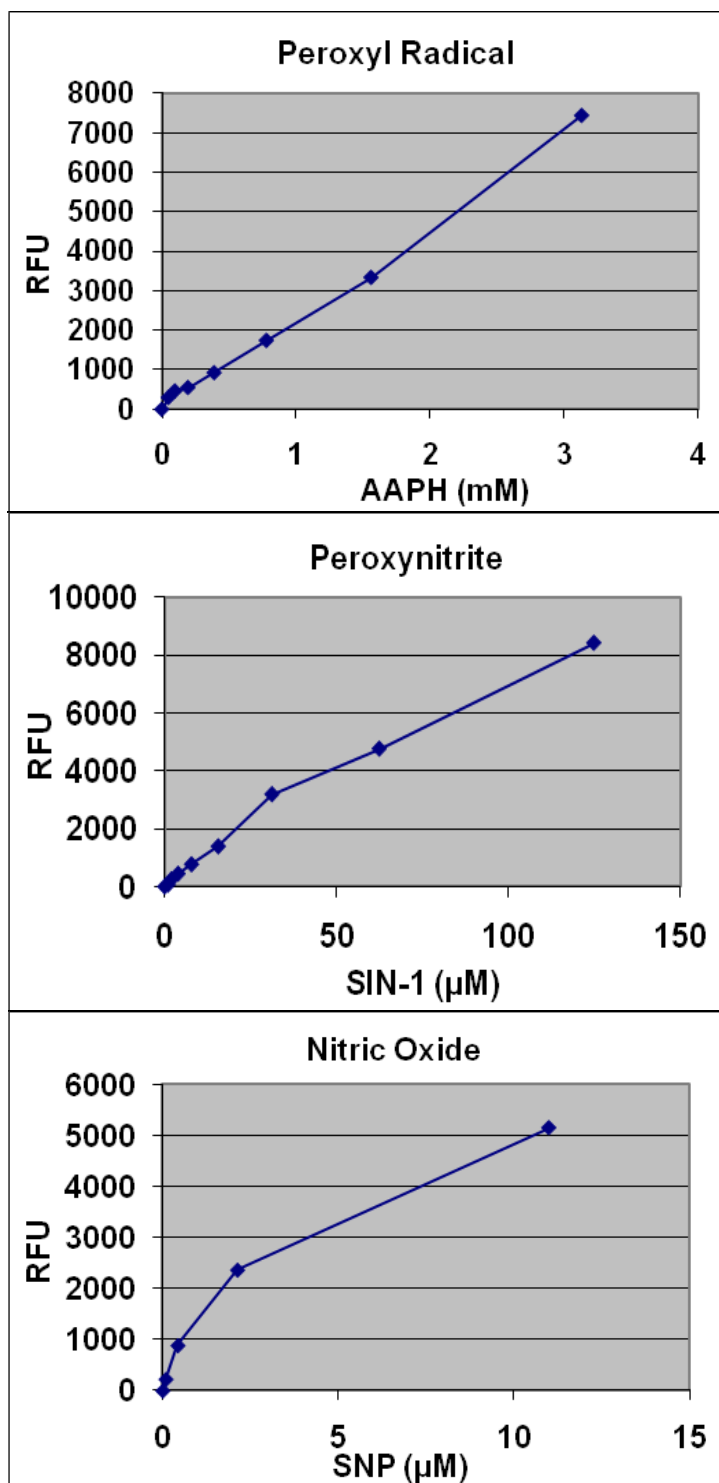


Figure 4. Detection of various free radical species using OxiSelect™ *In Vitro* ROS/RNS Assay Kit. DCF Fluorescence curves for AAPH (peroxyl radical generator, top), SIN-1 (peroxynitrite generator, center), and SNP (nitric oxide generator, bottom).

References

1. Bass DA, Parce JW, Dechatelet LR, Szejda P, Seeds MC, Thomas M. Flow cytometric studies of oxidative product formation by neutrophils: A graded response to membrane stimulation. *J Immunol.* 1983; 130:1910-1917.
2. Brandt R, Keston AS. Synthesis of diacetyldichlorofluorescein: A stable reagent for fluorometric analysis. *Anal Biochem.* 1965; 11:6-9.
3. Keston AS, Brandt R. The fluorometric analysis of ultramicro quantities of hydrogen peroxide. *Anal Biochem.* 1965; 11:1-5.

Recent Product Citations

1. Aljaser, F. et al. (2021). Effect of trace elements on the seminal oxidative status and correlation to sperm motility in infertile Saudi males. *Saudi J Biol Sci.* doi: 10.1016/j.sjbs.2021.04.042.
2. ALTamimi, J.Z. et al. (2021). Ellagic acid improved diabetes mellitus-induced testicular damage and sperm abnormalities by activation of Nrf2. *Saudi J Biol Sci.* doi: 10.1016/j.sjbs.2021.04.005.
3. Kim, D.H. et al. (2021). Comparison of therapeutic effects between topical 8-oxo-2'-deoxyguanosine and corticosteroid in ocular alkali burn model. *Sci Rep.* **11**(1):6909. doi: 10.1038/s41598-021-86440-7.
4. Shao, A. et al. (2021). Melatonin Ameliorates Hemorrhagic Transformation via Suppression of ROS-Induced NLRP3 Activation after Cerebral Ischemia in Hyperglycemic Rats. *Oxid Med Cell Longev.* **2021**:6659282. doi: 10.1155/2021/6659282.
5. ALTamimia, J.Z. et al. (2021). Ellagic acid protects against diabetic nephropathy in rats by regulating the transcription and activity of Nrf2. *J Funct Foods.* doi: 10.1016/j.jff.2021.104397.
6. Cuervo, W. et al. (2021). Oxidative Stress Compromises Lymphocyte Function in Neonatal Dairy Calves. *Antioxidants (Basel).* **10**(2):255. doi: 10.3390/antiox10020255.
7. BinMowyna, M.N. et al. (2021). Kaempferol suppresses acetaminophen-induced liver damage by upregulation/activation of SIRT1. *Pharm Biol.* **59**(1):146-156. doi: 10.1080/13880209.2021.1877734.
8. Lee, S.H. (2021). Human Adipose-Derived Stem Cells' Paracrine Factors in Conditioned Medium Can Enhance Porcine Oocyte Maturation and Subsequent Embryo Development. *Int J Mol Sci.* **22**(2):E579. doi: 10.3390/ijms22020579.
9. Grau, M. et al. (2021). Sub-Fractions of Red Blood Cells Respond Differently to Shear Exposure Following Superoxide Treatment. *Biology (Basel).* **10**(1):E47. doi: 10.3390/biology10010047.
10. Nakanishi, K. et al. (2021). High-Dose Vitamin C Administration Inhibits the Invasion and Proliferation of Melanoma Cells in Mice Ovary. *Biol Pharm Bull.* **44**(1):75-81. doi: 10.1248/bpb.b20-00637.
11. Wang, H. et al. (2020). Comprehensive Subchronic Inhalation Toxicity Assessment of an Indoor School Air Mixture of PCBs. *Environ Sci Technol.* doi: 10.1021/acs.est.0c04470.
12. Katerji, M. et al. (2020). Oxidative stress markers in patient-derived non-cancerous cervical tissues and cells. *Sci Rep.* **10**(1):19044. doi: 10.1038/s41598-020-76159-2.
13. Chan, K.C. et al. (2020). Effects of fermented red bean extract on nephropathy in streptozocin-induced diabetic rats. *Food Nutr Res.* doi: 10.29219/fnr.v64.4272.
14. Yeh, W.J. et al. (2020). *Hylocereus polyrhizus* Peel Extract Retards Alcoholic Liver Disease Progression by Modulating Oxidative Stress and Inflammatory Responses in C57BL/6 Mice. *Nutrients.* **12**(12):3884. doi: 10.3390/nu12123884.

15. Xu, Y. et al. (2020). Hepatocyte-specific Expression of Human Carboxylesterase 2 Attenuates Non-alcoholic Steatohepatitis in Mice. *Am J Physiol Gastrointest Liver Physiol*. doi: 10.1152/ajpgi.00315.2020.
16. Warowicka, A. et al. (2020). Alternations in mitochondrial genome in carcinogenesis of HPV positive cervix. *Exp Mol Pathol*. doi: 10.1016/j.yexmp.2020.104530.
17. Tonogawa, U. et al. (2020). Abnormal increases in reactive oxygen species in dying insects infected with nematodes. *Arch Insect Biochem Physiol*. doi: 10.1002/arch.21758.
18. Katerji, M. et al. (2020). Oxidative stress markers in patient-derived non-cancerous cervical tissues and cells. *Sci Rep*. **10**(1):19044. doi: 10.1038/s41598-020-76159-2.
19. Olejnik, A. et al. (2020). Ameliorating Effect of Klotho Protein on Rat Heart during I/R Injury. *Oxid Med Cell Longev*. doi: 10.1155/2020/6427284.
20. García-Laorden, M.I. et al. (Systemic Effects Induced by Hyperoxia in a Preclinical Model of Intra-abdominal Sepsis. *Mediators Inflamm*. doi: 10.1155/2020/5101834.
21. Itam, M. et al. (2020). Metabolic and physiological responses to progressive drought stress in bread wheat. *Sci Rep*. doi: 10.1038/s41598-020-74303-6.
22. Siregar, A.S. et al. (2020). Dipeptide YA is Responsible for the Positive Effect of Oyster Hydrolysates on Alcohol Metabolism in Single Ethanol Binge Rodent Models. *Mar. Drugs*. **18**(10):512. doi: 10.3390/md18100512.
23. Handa, K. et al. (2020). Bone loss caused by dopaminergic degeneration and levodopa treatment in Parkinson's disease model mice. *Sci Rep*. **9**(1):13768. doi: 10.1038/s41598-019-50336-4.
24. Opgenorth, J. et al. (2020). Colostrum supplementation with n-3 fatty acids alters plasma polyunsaturated fatty acids and inflammatory mediators in newborn calves. *J Dairy Sci*. doi: 10.3168/jds.2019-18045.
25. Barroso, E. et al. (2020). SIRT3 deficiency exacerbates fatty liver by attenuating the HIF1 α -LIPIN 1 pathway and increasing CD36 through Nrf2. *Cell Commun Signal*. **18**(1):147. doi: 10.1186/s12964-020-00640-8.
26. Seon, G. et al. (2020). Effect of post-treatment process of microalgal hydrolysate on bioethanol production. *Sci Rep*. **10**(1):16698. doi: 10.1038/s41598-020-73816-4.
27. ALTamimi, J.Z. et al. (2020). Curcumin reverses diabetic nephropathy in streptozotocin-induced diabetes in rats by inhibition of PKC β /p⁶⁶Shc axis and activation of FOXO-3a. *J Nutr Biochem*. doi: 10.1016/j.jnutbio.2020.108515.
28. Hiramoto, K. et al. (2020). Innate immune activation and antitumor effects of Lactobacillus-fermented Sparassis crispa extract in mice. *J Funct Foods*. doi: 10.1016/j.jff.2020.104215.
29. Walter, L. et al. (2020). Matrix metalloproteinase 9 (MMP9) limits reactive oxygen species (ROS) accumulation and DNA damage in colitis-associated cancer. *Cell Death Dis*. **11**(9):767. doi: 10.1038/s41419-020-02959-z.
30. Langbøl, M. et al. (2020). Increased Antioxidant Capacity and Pro-Homeostatic Lipid Mediators in Ocular Hypertension-A Human Experimental Model. *J Clin Med*. **9**(9):E2979. doi: 10.3390/jcm9092979.

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

©2010-2021: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.