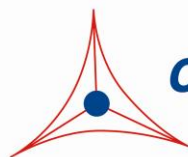

Product Manual

OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)

Catalog Number

STA-342	96 assays
STA-342-5	5 x 96 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.
Creating Solutions for Life Science Research

Introduction

Accumulation of reactive oxygen species (ROS) 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, and cancer 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 activity intracellularly is crucial to suppressing or treating oxidative stress inducers.

Cell Biolabs' OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence) is a cell-based assay for measuring hydroxyl, peroxy, or other reactive oxygen species activity within a cell. The assay employs the cell-permeable fluorogenic probe 2', 7'-Dichlorodihydrofluorescein diacetate (DCFH-DA). In brief, DCFH-DA is diffused into cells and is deacetylated by cellular esterases to non-fluorescent 2', 7'-Dichlorodihydrofluorescein (DCFH), which is rapidly oxidized to highly fluorescent 2', 7'-Dichlorodihydrofluorescein (DCF) by ROS (Figure 1). The fluorescence intensity is proportional to the ROS levels within the cell cytosol. The effect of antioxidant or free radical compounds on DCFH-DA can be measured against the fluorescence of the provided DCF standard. The kit has a DCF detection sensitivity limit of 10 pM. Each kit provides sufficient reagents to perform up to 96 assays, including standard curve and unknown samples.

Assay Principle

The OxiSelect™ Intracellular ROS Assay Kit is a cell-based assay for measuring antioxidant or ROS activity. Cells are cultured in a 96-well cell culture plate and then pre-incubated with DCFH-DA, which is cell-permeable (Figure 1). The unknown antioxidant or ROS samples are then added to the cells. After a brief incubation, the cells can be read on a standard fluorescence plate reader. The ROS or antioxidant content in unknown samples is determined by comparison with the predetermined DCF standard curve.

Related Products

1. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
2. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
3. STA-341: OxiSelect™ Catalase Activity Assay Kit
4. STA-343: OxiSelect™ Hydrogen Peroxide Assay Kit
5. STA-344: OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit
6. STA-345: OxiSelect™ ORAC Activity Assay Kit
7. STA-347: OxiSelect™ In Vitro ROS/RNS Assay Kit (Green Fluorescence)
8. STA-350: OxiSelect™ Comet Assay Kit (3-Well Slides), 15 assays
9. STA-832: OxiSelect™ MDA Adduct Competitive ELISA Kit
10. STA-838: OxiSelect™ HNE Adduct Competitive ELISA Kit

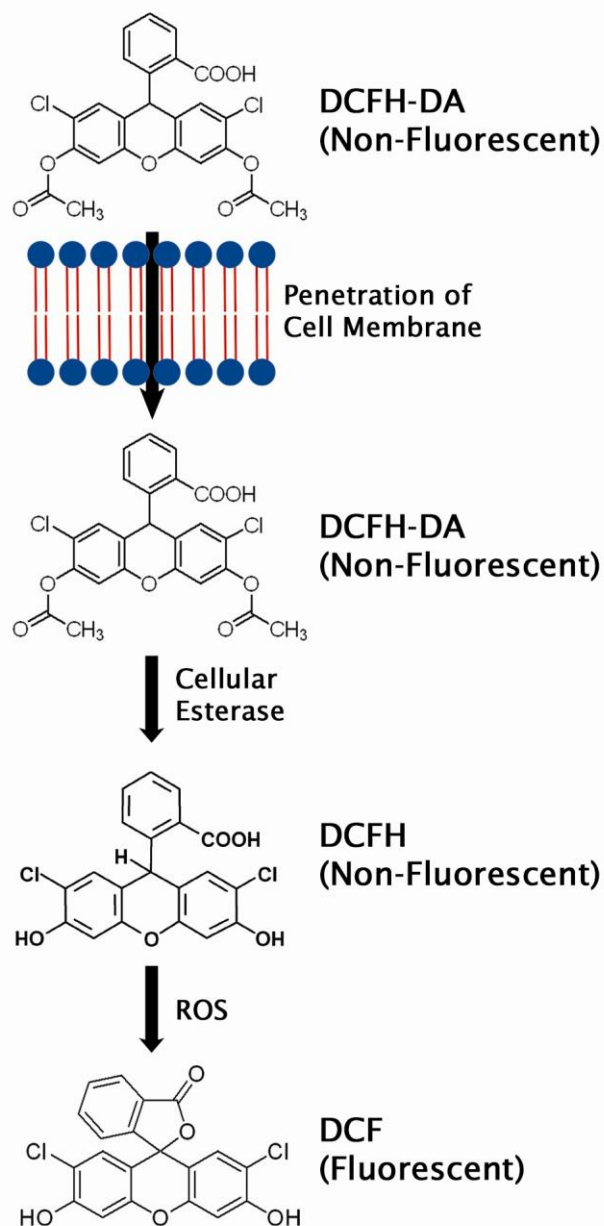


Figure 1. Mechanism of DCF Assay

Kit Components

1. 20X DCFH-DA (Part No. 234201): One 500 μ L amber tube of a 20 mM solution in methanol.
2. DCF Standard (Part No. 234202): One 100 μ L amber tube of a 1 mM solution in DMSO.
3. Hydrogen Peroxide (Part No. 234102): One 100 μ L amber tube of an 8.821 M solution.
4. 2X Cell Lysis Buffer (Part No. 234203): One 20 mL bottle.

Materials Not Supplied

1. Sterile DPBS for washes and buffer dilutions
2. Hank's Balanced Salt Solution (HBSS)
3. Cell culture medium (ie: DMEM +/-10% FBS)
4. 96-well black or fluorescence microtiter plate
5. Fluorescent microplate reader capable of reading 480 nm (excitation) and 530 nm (emission)

Storage

Upon receipt, store the DCFH-DA and DCF Standard at -20°C. Avoid multiple freeze/thaw cycles. Store the Cell Lysis Buffer and Hydrogen Peroxide at 4°C.

Preparation of Reagents

- 1X DCFH-DA: Dilute the 20X DCFH-DA stock solution to 1X in cell culture media, preferably without FBS. Stir or vortex to homogeneity. Prepare only enough for immediate applications.

Notes:

- *1X DCFH-DA/media solution contains 5% methanol. For cells that are sensitive to methanol, we recommend instead preparing a 0.1X (100 µM) solution of DCFH-DA in cell culture media.*
- *Due to light-induced auto-oxidation, DCFH-DA solutions at any concentration must be protected from light.*
- Hydrogen Peroxide (H₂O₂): Prepare H₂O₂ dilutions in DMEM or DPBS as necessary. Do not store diluted solutions. Hydrogen Peroxide may be used as a positive control in the assay, or as a cell treatment.

Preparation of Standard Curve

1. Prepare a 1:10 dilution series of DCF standards in the concentration range of 0 µM – 10 µM by diluting the 1 mM DCF stock in cell culture media (see Table 1).

Standard Tubes	DCF Standard (µL)	Culture Medium (µL)	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	100 of Tube #5	900	0.1
7	100 of Tube #6	900	0.01
8	0	1000	0

Table 1. Preparation of DCF Standards

2. Transfer 75 μL of each DCF standard to a 96-well plate suitable for fluorescence measurement. Add 75 μL of the 2X Cell Lysis Buffer.
3. Read the fluorescence with a fluorescence plate reader at 480 nm excitation /530 nm emission.

Assay Protocol

I. DCF Dye Loading

1. Prepare and mix all reagents thoroughly before use. Each unknown sample should be assayed in duplicate or triplicate.
2. Culture cells in either a clear or black 96-well cell culture plate.
Note: If using a black plate, choose an appropriate plate based on your fluorometer's reader (i.e. choose a clear bottom black plate for bottom readers).
3. Remove media from all wells and discard. Wash cells gently with DPBS or HBSS 2-3 times. Remove the last wash and discard.
4. Add 100 μL of 1X DCFH-DA/media solution to cells. Incubate at 37°C for 30-60 minutes.
5. Remove solution. Repeat step three using multiple washes with DPBS or HBSS. Remove the last wash and discard.
6. Treat DCFH-DA loaded cells with desired oxidant or antioxidant in 100 μL medium.

II. Quantitation of Fluorescence

- Fluorescence microscopy or Flow cytometry: Fluorescence can be analyzed on an inverted fluorescence microscope or by flow cytometry using excitation and emission wavelengths of 480 nm and 530 nm, respectively.
- Fluorescence Plate Reader:
 - Assays performed in black cell culture fluorometric plates: Plate may be read immediately for kinetic analysis or after 1 hour for static analysis. Plates read for kinetic analysis may be read in increments of 1 and 5 minutes up to 1 hour or more as necessary. Read the fluorescence with a fluorometric plate reader at 480 nm/530 nm.
 - Assays performed in clear cell culture plates: After treatment with desired oxidant or antioxidant, carefully remove treatment media from all wells and discard. Wash cells gently with DPBS or HBSS 2-3 times. Remove the last wash and discard. Add 100 μL of medium to each well. Add 100 μL of the 2X Cell Lysis Buffer, mix thoroughly and incubate 5 minutes. Transfer 150 μL of the mixture to a 96-well plate suitable for fluorescence measurement. Read the fluorescence with a fluorometric plate reader at 480 nm/530 nm.

Example of Results

The following figures demonstrate typical ROS 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.

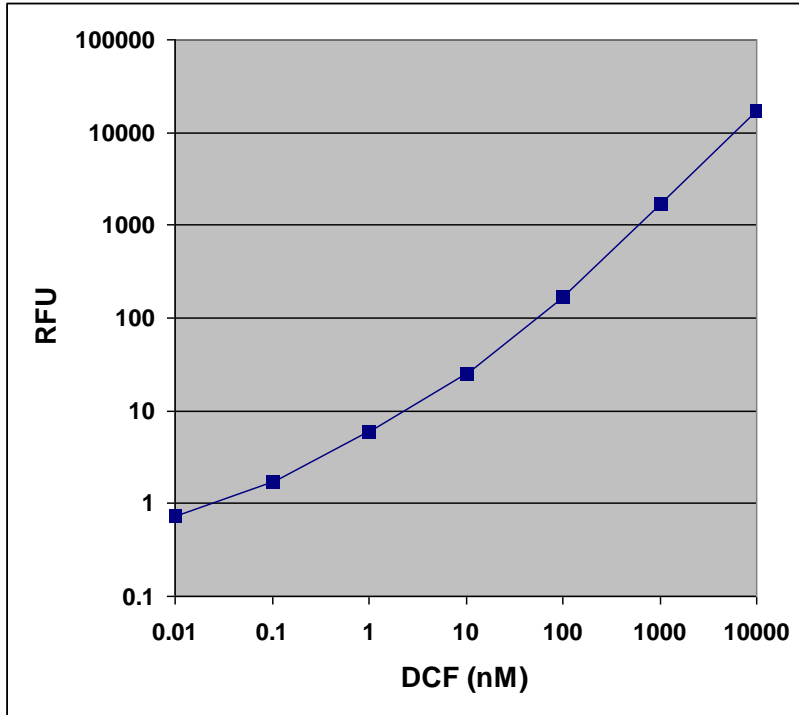


Figure 2. DCF Standard Curve.

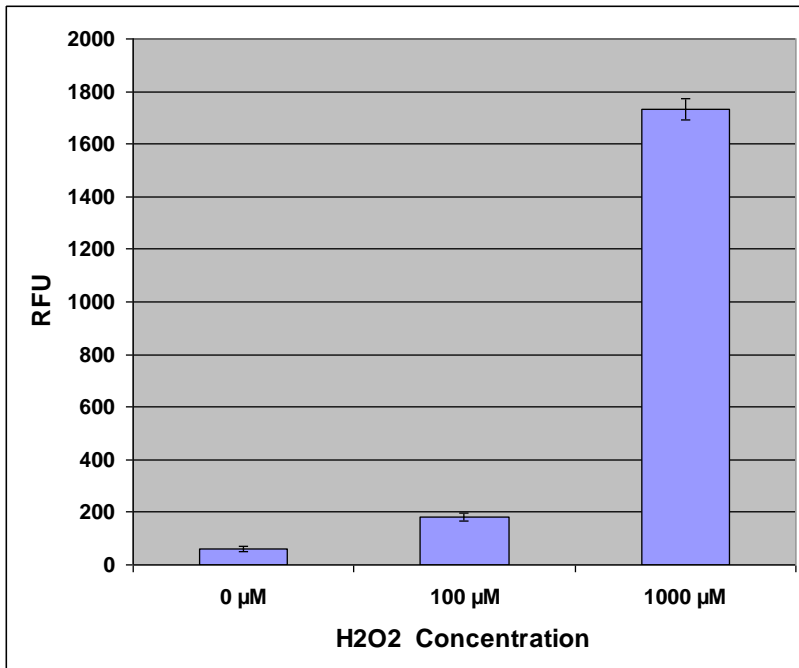


Figure 3. ROS in HeLa cells treated with H₂O₂. 50,000 HeLa cells in a 96-well plate were first pretreated with 1 mM DCFH-DA for 60 minutes at 37°C. Cells were then treated with various concentrations of H₂O₂ for 20 minutes.

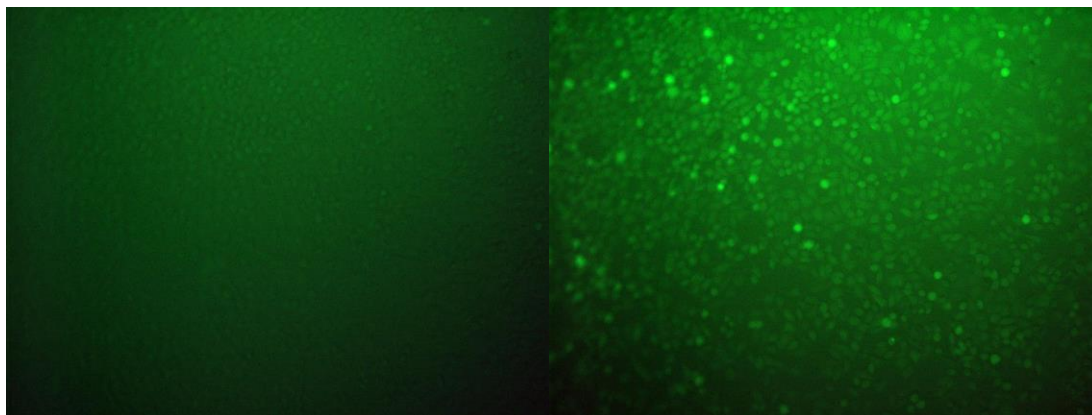


Figure 4. DCF Fluorescence in H₂O₂ treated HeLa cells after 1 hour. Left: 0 μM; Right: 1000 μM.

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. Li, X.M., Zhang, X.J., Dong, M.X. (2017). Isorhynchophylline Attenuates MPP⁺-Induced Apoptosis Through Endoplasmic Reticulum Stress- and Mitochondria-Dependent Pathways in PC12 Cells: Involvement of Antioxidant Activity. *Neuromolecular Med.* doi: 10.1007/s12017-017-8462-x
2. Kim, J. et al. (2017). Aldehyde dehydrogenase 2*2 knock-in mice show increased reactive oxygen species production in response to cisplatin treatment. *J Biomed Sci.* **24(1)**:33.
3. Osamura, T. et al. (2017). Specific expression and function of the A-type cytochrome c oxidase under starvation conditions in *Pseudomonas aeruginosa*. *PLoS One* **12(5)**:e0177957.
4. Palao-Suay, R. et al. (2017). Photothermal and photodynamic activity of polymeric nanoparticles based on α -tocopheryl succinate-RAFT block copolymers conjugated to IR-780. *Acta Biomater.* **57**:70-84.
5. Benedetti, F. et al. (2017). Anti-inflammatory effects of H₂S during acute bacterial infection: a review. *J. Transl. Med.* **15(1)**:100.
6. Yin, S. et al. (2017). Environmental temperature affects physiology and survival of nanosecond pulsed electric field-treated cells. *J. Cell. Physiol.* doi:10.1002/jcp.25984.
7. Stebbins, K.J. et al. (2017). In vitro and in vivo pharmacology of NXT629, a novel and selective PPAR α antagonist. *Eur. J. Pharmacol.* **809**:130-140.
8. Sweet, C. et al. (2017). Two-photon fluorescent molybdenum disulfide dots for targeted prostate cancer imaging in the biological II window. *ACS Omega* **2(5)**:1826-1835.
9. Haga, H., et al. (2017). Extracellular vesicles from bone marrow derived mesenchymal stem cells protect against murine hepatic ischemia-reperfusion injury. *Liver Transpl.* doi: 10.1002/lt.24770.

10. Vijayarathna, S. et al. (2017). Polyalthia longifolia Methanolic Leaf Extracts (PLME) induce apoptosis, cell cycle arrest and mitochondrial potential depolarization by possibly modulating the redox status in hela cells. *Biomed Pharmacother.* **89**:499-514. doi: 10.1016/j.biopha.2017.02.075.
11. Kavitha, N. et al. (2017). Phaleria macrocarpa (Boerl.) fruit induce G0/G1 and G2/M cell cycle arrest and apoptosis through mitochondria-mediated pathway in MDA-MB-231 human breast cancer cell. *J Ethnopharmacol.* **201**:42-55. doi: 10.1016/j.jep.2017.02.041.
12. Xu, Z. and Kong, X.Q. (2017). Bixin ameliorates high fat diet-induced cardiac injury in mice through inflammation and oxidative stress suppression. *Biomed Pharmacother.* **89**:991-1004. doi: 10.1016/j.biopha.2017.02.052.
13. Kuete, V. et al (2017). In vitro cytotoxicity of compounds isolated from Desbordesia glaucescens against human carcinoma cell lines. *South African Journal of Botany.* **111**: 37–43.
14. Yu C, et al. (2017). Astragaloside IV attenuates lead acetate-induced inhibition of neurite outgrowth through activation of Akt-dependent Nrf2 pathway in vitro. *Biochim Biophys Acta.* pii: S0925-4439(17)30082-0. doi: 10.1016/j.bbadis.2017.03.006.
15. Acevedoa, F. et al (2017). Volatile and non-volatile/semi-volatile compounds and in vitro bioactive properties of Chilean Ulmo (Eucryphia cordifolia Cav.) honey. *Food Research International.* **94**:20-28. <http://dx.doi.org/10.1016/j.foodres.2017.01.021>
16. Crookenden MA, (2017). Effects of precalving body condition and prepartum feeding level on gene expression in circulating neutrophils. *J Dairy Sci.* **100**(3):2310-2322. doi: 10.3168/jds.2016-12105.
17. Golestaneh, N., et al. (2017). Dysfunctional autophagy in RPE, a contributing factor in age-related macular degeneration. *Cell Death and Disease.* **8**(1):e2537. doi:10.1038/cddis.2016.453
18. Ludwig R, et al. (2017). Nanoparticle-based hyperthermia distinctly impacts production of ROS, expression of Ki-67, TOP2A, and TPX2, and induction of apoptosis in pancreatic cancer. *Int J Nanomedicine.* **12**:1009-1018. doi: 10.2147/IJN.S108577.
19. Zhang X, et al (2017). Different virulence of candida albicans is attributed to the ability of escape from neutrophil extracellular traps by secretion of DNase. *Am J Transl Res.* **9**(1):50-62.
20. Paul A, et al (2016). Nanoengineered biomimetic hydrogels for guiding human stem cell osteogenesis in three dimensional microenvironments. *J Mater Chem B Mater Biol Med.* May **28**;4(20):3544-3554.
21. Yao, J. et al. (2016). Pinoresinol diglucoside alleviates oxLDL-induced dysfunction in human umbilical vein endothelial cells. *Evid. Based Complement Alternat. Med.* 2016:3124519.
22. Zhao, Q. et al. (2016). Sulodexide protects contrast-induced nephropathy in Sprague-Dawley rats. *Cell Physiol. Biochem.* **40**:621-632.
23. Cao, X. et al. (2016). TUSC2(FUS1)-erlotinib induced vulnerabilities in Epidermal Growth Factor Receptor (EGFR) wildtype non-small cell lung cancer (NSCLC) targeted by the repurposed drug auranofin. *Sci Rep.* **6**:35741.
24. Golestaneh, N. et al. (2016). Repressed SIRT1/PGC-1 α pathway and mitochondrial disintegration in iPSC-derived RPE disease model of age-related macular degeneration. *J. Transl. Med.* **14**:344.
25. Kuete, V. et al. (2016). Cytotoxicity of Plumbagin, Rapanone and 12 other naturally occurring Quinones from Kenyan flora towards human carcinoma cells. *BMC Pharmacol. Toxicol.* **17**:60.
26. Chayapong, J. et al. (2016). Arsenic trioxide induces ROS activity and DNA damage, leading to G0/G1 extension in skin fibroblasts through the ATM-ATR-associated Chk pathway. *Env. Sci. Pollut. Res. Int.* **24**(6):5316-5325. doi:10.1007/s11356-016-8215-7.

27. Feng, J. et al. (2016). Inhibitor of nicotinamide phosphoribosyltransferase sensitizes glioblastoma cells to temozolomide via activating ROS/JNK signaling pathway. *BioMed Res. Int.* doi:10.1155/2016-1450843.
28. Yao, J. et al. (2016). Pinoresinol diglucoside alleviates oxLDL-induced dysfunction in human umbilical vein endothelial cells. *Evid. Based Complement Alternat. Med.* **2016**:3124519.
29. Zhao, Q. et al. (2016). Sulodexide protects contrast-induced nephropathy in Sprague-Dawley rats. *Cell Physiol. Biochem.* **40**:621-632.
30. Cao, X. et al. (2016). TUSC2(FUS1)-erlotinib induced vulnerabilities in Epidermal Growth Factor Receptor (EGFR) wildtype non-small cell lung cancer (NSCLC) targeted by the repurposed drug auranofin. *Sci Rep.* **6**:35741.
31. Xiaobo, C. et al. (2016). TUSC2 (FUS1)-erlotinib induced vulnerabilities in epidermal growth factor receptor (EGFR) wildtype non-small cell lung cancer (NSCLC) targeted by the repurposed drug auranofin. *Sci. Rep.* **6**:35741.
32. Chen, L. et al. (2016). Neuroprotective effects of vitexin against isoflurane-induced neurotoxicity by targeting the TRPV1 and NR2B signaling pathways. *Mol. Med. Rep.* **14**:5607-5613.
33. Brufau, M.T. et al. (2016). Salmosan®, a β -galactomannan-rich product, in combination with *Lactobacillus plantarum* contributes to restore intestinal epithelial barrier function by modulation of cytokine production. *J. Nutr. Biochem.* doi:10.1016/j.jnutbio.2016.11.011.
34. Liu, H. et al. (2016). Synergistic protective effect of paeoniflorin and β -ecdysterone against rotenone-induced neurotoxicity in PC12 cells. *Apoptosis* doi:10.1007/s10495-016-1293-7.
35. Koo, M. A. et al. (2016). Controlled delivery of extracellular ROS based on hematoporphyrin-incorporated polyurethane film for enhanced proliferation of endothelial cells. *ACS Appl Mater Interfaces.* doi:10.1021/acsami.6b07628.
36. Takayama, K. et al. (2016). Nuclear factor (erythroid-derived)-related factor 2-associated retinal pigment epithelial cell protection under blue light-induced oxidative stress. *Oxid Med Cell Longev.* doi:10.1155/2016/8694641.

Please see the complete list of product citations: <http://www.cellbiolabs.com/reactive-oxygen-species-ros-assay>.

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

©2009-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.