
Product Manual

OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence), Trial Size

Catalog Number

STA-342-T

20 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 Trial Size Intracellular ROS Assay Kit provides sufficient reagents to perform up to 20 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

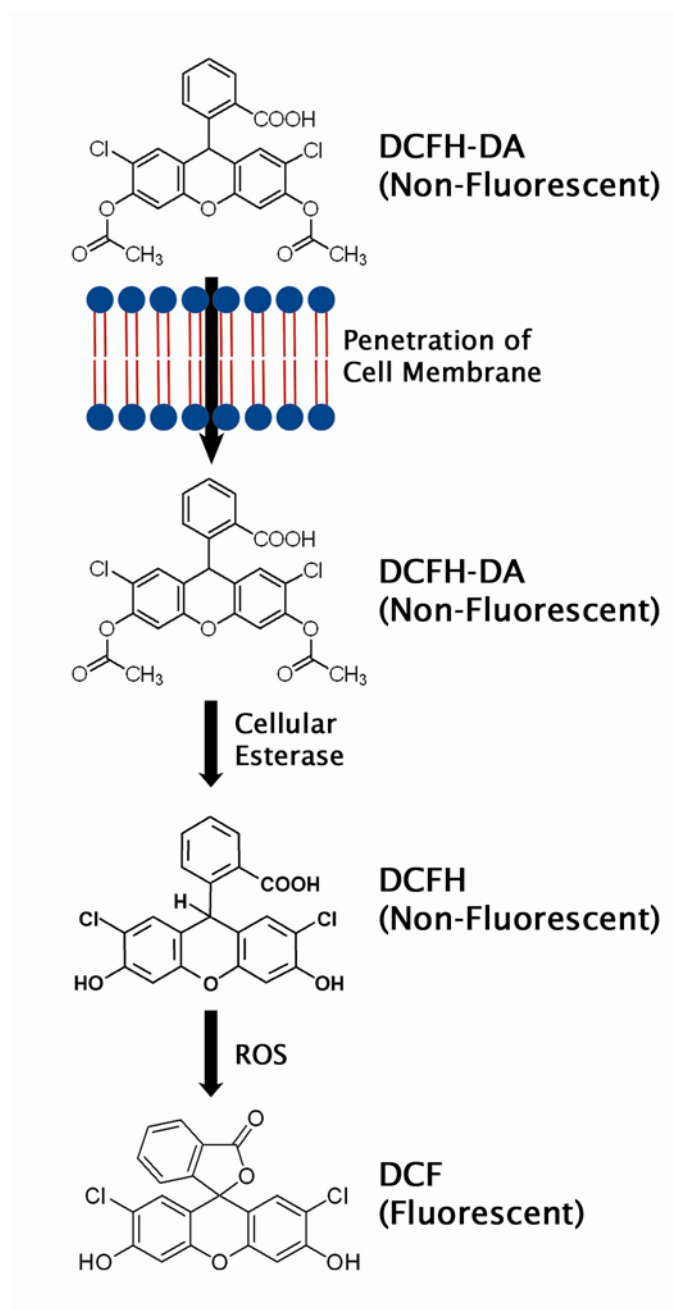


Figure 1. Mechanism of DCF Assay

Kit Components

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

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 μ L of each DCF standard to a 96-well plate suitable for fluorescence measurement. Add 75 μ L 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 μ L 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 μ L 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 μ L of medium to each well. Add 100 μ L of the 2X Cell Lysis Buffer, mix thoroughly and incubate 5 minutes. Transfer 150 μ L 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.

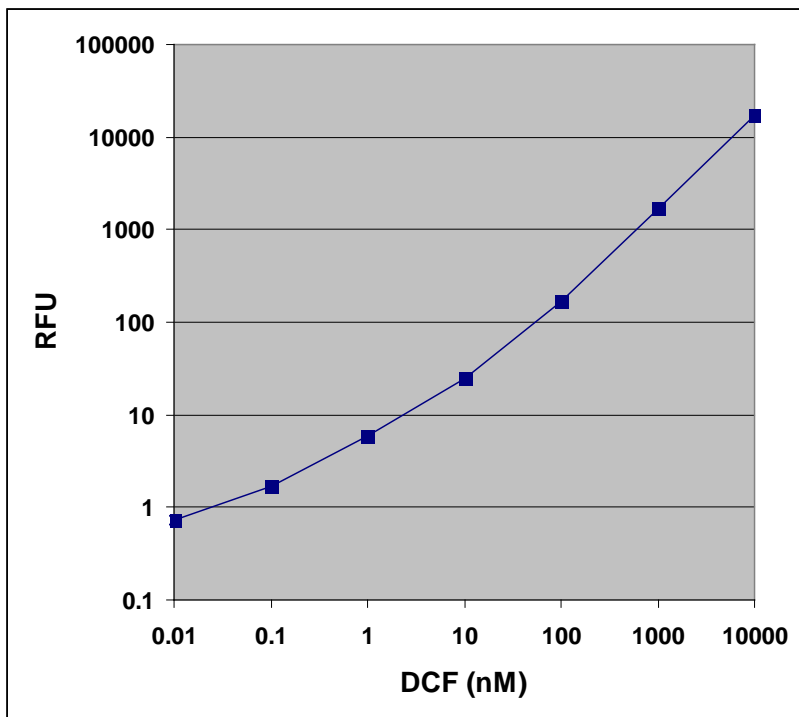


Figure 2. DCF Standard Curve.

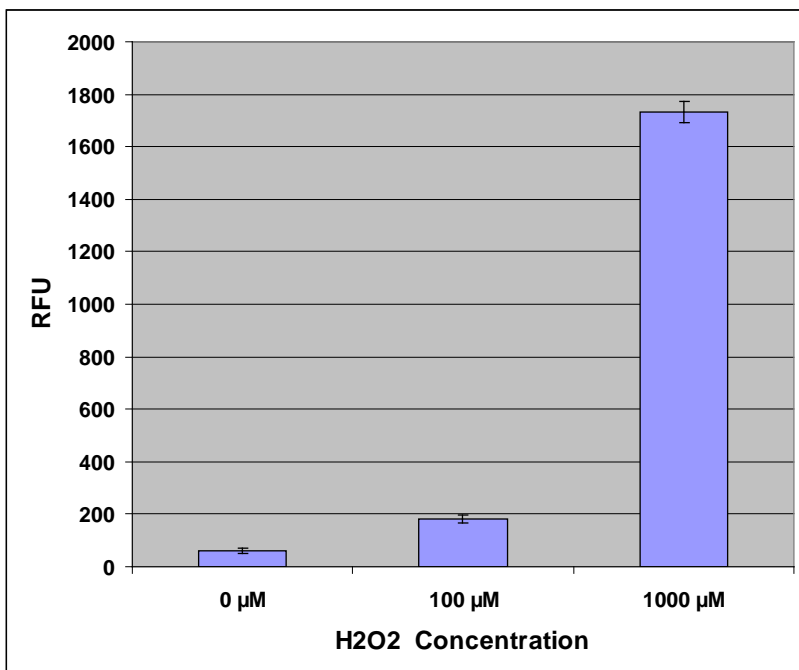


Figure 3. ROS in HeLa cells treated with H₂O₂. 50,000 HeLa cells in a 96-well plate were first pretreated with 1 mM DCFH-DA for 60 minutes at 37°C. Cells were then treated with various concentrations of H₂O₂ for 20 minutes.

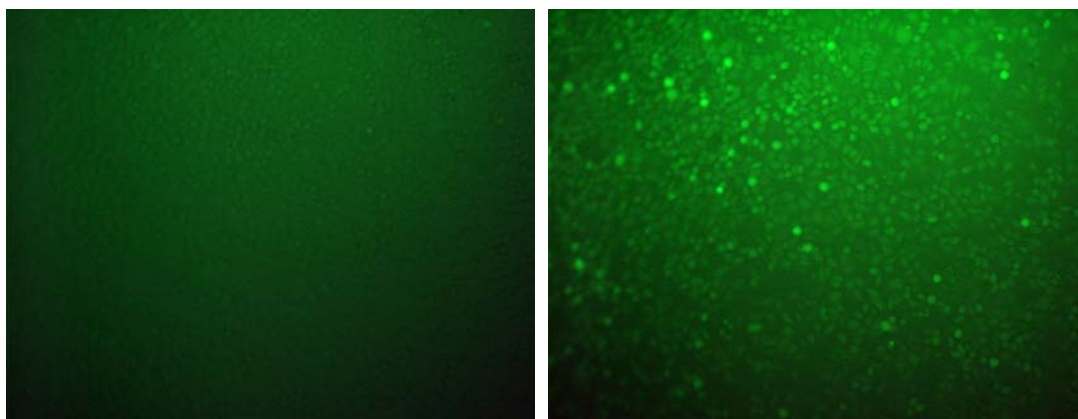


Figure 4. DCF Fluorescence in H₂O₂ treated HeLa cells after 1 hour. Left: 0 μM; Right: 1000 μM.

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. Hirata, T. et al. (2018). 4-Methylthio-3-butenyl isothiocyanate mediates nuclear factor (erythroid-derived 2)-like 2 activation by regulating reactive oxygen species production in human esophageal epithelial cells. *Food Chem Toxicol.* **111**:295-301. doi: 10.1016/j.fct.2017.11.022.
2. Li, L. et al. (2018). Association of anti-HER2 antibody with graphene oxide for curative treatment of osteosarcoma. *Nanomedicine.* **14**(2):581-593. doi: 10.1016/j.nano.2017.11.001.
3. Kim, K.Y. et al. (2017). Mitochondria-targeting self-assembled nanoparticles derived from triphenylphosphonium-conjugated cyanostilbene enable site-specific imaging and anticancer drug delivery. *Nano Res.* <https://doi.org/10.1007/s12274-017-1728-7>
4. Xu, M.J. et al. (2017). Epigallocatechin-3-gallate inhibits TLR4 signaling through the 67-kDa laminin receptor and effectively alleviates acute lung injury induced by H9N2 swine influenza virus. *Int Immunopharmacol.* **52**:24-33. doi: 10.1016/j.intimp.2017.08.023.
5. Hara, K. et al. (2017). Scavenging of reactive oxygen species by astaxanthin inhibits epithelial–mesenchymal transition in high glucose-stimulated mesothelial cells. *PLoS One.* **12**(9): e0184332. doi: 10.1371/journal.pone.0184332
6. Choi, H.N. and Yim, J.E. (2017). Effects of Erythrocyte Membrane Polyunsaturated Fatty Acids in Overweight, Obese, and Morbidly Obese Korean Women. *J Cancer Prev.* **22**(3):182-188. doi: 10.15430/JCP.2017.22.3.182.
7. Sonntag, K.C. et al. (2017). Late-onset Alzheimer's disease is associated with inherent changes in bioenergetics profiles. *Sci Rep.* **7**(1):14038. doi: 10.1038/s41598-017-14420-x.

8. Azimi, I. et al. (2017). Hypoxia-induced reactive oxygen species mediate N-cadherin and SERPINE1 expression, EGFR signalling and motility in MDA-MB-468 breast cancer cells. *Sci Rep.* **7**(1):15140. doi: 10.1038/s41598-017-15474-7.
9. Abd Jalil, A. et al. (2017). Vitamin E-Mediated Modulation of Glutamate Receptor Expression in an Oxidative Stress Model of Neural Cells Derived from Embryonic Stem Cell Cultures. *Evid Based Complement Alternat Med.* 2017:6048936. doi: 10.1155/2017/6048936.
10. 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
11. 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.
12. 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.
13. 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.
14. Benedetti, F. et al. (2017). Anti-inflammatory effects of H₂S during acute bacterial infection: a review. *J. Transl. Med.* **15**(1):100.
15. 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.
16. 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.
17. 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.
18. 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.
19. 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.
20. 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.
21. 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.
22. 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.
23. 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.
24. 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>

25. 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.
26. 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
27. 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.
28. 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.
29. 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.
30. Yao, J. et al. (2016). Pinoresinol diglucoside alleviates oxLDL-induced dysfunction in human umbilical vein endothelial cells. *Evid. Based Complement Alternat. Med.* 2016:3124519.

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

©2014-2018: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.