
Product Manual

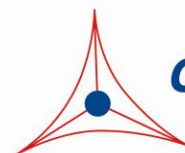
OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit (Fluorometric)

Catalog Number

STA-344

500 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures

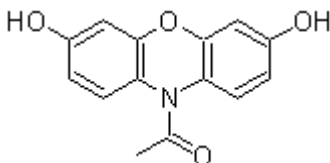


CELL BIOLABS, INC.
Creating Solutions for Life Science Research

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), are some 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/Peroxidase Assay Kit is a simple HTS-compatible assay for measuring hydrogen peroxide concentrations or peroxidase activities in biological samples without any need for pretreatment. In the presence of H₂O₂ and horseradish peroxidase (HRP), non-fluorescent ADHP (10-Acetyl-3, 7-dihydroxyphenoxazine) is oxidized to the highly fluorescent Resorufin. The probe has less background and greater stability. The ADHP-based H₂O₂ detection is at least one order of magnitude more sensitive than the commonly used Xylenol Orange (FOX) colorimetric assay for H₂O₂. The probe can be also used as an ultrasensitive assay for peroxidase activity when H₂O₂ is in excess. The kit has a detection sensitivity limit of 50 nM (H₂O₂) or 0.1 mU/mL (Peroxidase). Each kit provides sufficient reagents to perform up to 500 assays, including standard curve and unknown samples.



ADHP (10-Acetyl-3, 7-dihydroxyphenoxazine)

Assay Principle

The OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit is a sensitive quantitative fluorometric assay for hydrogen peroxide or peroxidase. In the presence of HRP, ADHP reacts with H₂O₂ in a 1:1 stoichiometry to produce highly fluorescent Resorufin. The Resorufin product can be easily read by a fluorescence microplate reader with an excitation of 530-560 nm and an emission of 590 nm. Fluorescence values are proportional to the H₂O₂ or peroxidase levels within the samples. The H₂O₂ or peroxidase content in unknown samples is determined by comparison with its respective 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-342: OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)
5. STA-345: OxiSelect™ ORAC Activity Assay Kit

Kit Components (shipped at room temperature)

1. ADHP Probe (Part No. 234401): One 250 µL amber tube of a 10 mM solution in DMSO.
2. HRP (Part No. 234402): One 100 µL tube of a 100 U/mL solution in glycerol*.
3. Hydrogen Peroxide (Part No. 234102): One 100 µL amber tube of an 8.8 M solution.
4. 10X Assay Buffer (Part No. 234403): One 25 mL bottle.

**Note: One unit is defined as the amount of enzyme that will form 1.0 mg purpurogallin from pyrogallol in 20 seconds at pH 6.0 and 20°C.*

Materials Not Supplied

1. Distilled or deionized water
2. 1X PBS for sample dilutions
3. Standard 96-well fluorescence black microtiter plate and/or black cell culture microplate

Storage

Upon receipt, aliquot and store the ADHP probe and HRP at -20°C. Avoid multiple freeze/thaw cycles. Store the remaining kit components at 4°C. ADHP is light sensitive, must be stored accordingly.

Preparation of Reagents

Note: All reagents must be brought to room temperature prior to use.

- 1X Assay Buffer: Dilute the stock 10X Assay Buffer 1:10 with deionized water for a 1X solution. Stir or vortex to homogeneity.
- ADHP/HRP Working Solution (Hydrogen Peroxide Assay): If measuring hydrogen peroxide, prepare an ADHP/HRP Working Solution by adding ADHP to a final concentration of 100 µM and HRP to a final concentration of 0.2 U/mL in 1X Assay Buffer (eg. Add 50 µL ADHP stock solution and 10 µL HRP stock solution to 4.940 mL 1X Assay Buffer). This volume is enough for ~100 assays. The ADHP/HRP Working Solution is stable for 1 day. Prepare only enough for immediate use.
- ADHP/H₂O₂ Working Solution (Peroxidase Assay): If measuring peroxidases, prepare the ADHP/H₂O₂ Working Solution by adding ADHP to a final concentration of 100 µM and H₂O₂ to a final concentration of 2 mM in 1X Assay Buffer. First perform a 1:1000 dilution of the stock H₂O₂ in 1X Assay Buffer. Use only enough for immediate applications (eg. Add 5 µL of H₂O₂ to 4.995 mL 1X Assay Buffer). This solution has a concentration of 8.8 mM. Use this 8.8 mM H₂O₂ solution to prepare a 2 mM H₂O₂ solution in ADHP/1X Assay Buffer (eg. Add 50 µL ADHP stock

solution and 1.14 mL of the prepared 8.8 mM H₂O₂ solution to 3.81 mL 1X Assay Buffer). This volume is enough for ~100 assays. The Working Solution is stable for 1 day. Prepare only enough for immediate use.

Preparation of Samples

- Cell culture supernatant: To remove insoluble particles, centrifuge at 10,000 rpm for 5 min. The supernatant can be assayed directly or diluted as necessary. Prepare the H₂O₂ standard curve in the same non-conditioned media. Serum should be avoided, as it interferes with the assay.

Note: Maintain pH between 7 and 8 for optimal working conditions as the ADHP is unstable at high pH (>8.5).

- Cell lysate: Resuspend cells at 1-2 x 10⁶ cells/mL in PBS or 1X Assay Buffer. Homogenize or sonicate the cells on ice. Centrifuge to remove debris. Cell lysates can be assayed undiluted or titrated as necessary.
- Plasma, Serum or Urine: To remove insoluble particles, centrifuge at 10,000 rpm for 5 min. The supernatant can be assayed directly or diluted as necessary. Undiluted serum or plasma may interfere with the assay.

Notes:

- *All samples should be assayed immediately or stored at -80°C for up to 1-2 months. Run proper controls as necessary. Optimal experimental conditions for samples must be determined by the investigator. Always run a standard curve with samples.*
- *A serial dilution will be necessary depending on the total H₂O₂ or peroxidase present. Extremely high levels of H₂O₂ (≥ 500 μM final concentration) or peroxidase (≥ 100 mU/mL) can lower the fluorescence because excess H₂O₂ or peroxidase can further oxidize the reaction product, Resorufin, to nonfluorescent product Resazurin.*
- *Samples with NADH concentrations above 10 μM and glutathione concentrations above 50 μM will oxidize the ADHP probe and could result in erroneous readings. To minimize this interference, it is recommended that superoxide dismutase (SOD) be added to the reaction at a final concentration of 40 U/mL (Votyakova and Reynolds, Ref. 2).*
- *Avoid samples containing DTT or β-mercaptoethanol since Resorufin is not stable in the presence of thiols (above 10 μM).*

Preparation of Standard Curves

1. **H₂O₂ Standard:** To prepare the H₂O₂ standards, first perform a 1:1000 dilution of the stock H₂O₂ in 1X Assay Buffer. Prepare only enough for immediate use (e.g., Add 5 μL of H₂O₂ to 4.995 mL 1X Assay Buffer). This solution has a concentration of 8.8 mM.
2. Use this 8.8 mM H₂O₂ solution to prepare standards in the concentration range of 0 μM – 100 μM by further diluting in 1X Assay Buffer; see Table 1 below. H₂O₂ diluted solutions and standards should be prepared fresh.

Standard Tubes	8.8 mM H ₂ O ₂ Standard (μL)	1X Assay Buffer (μL)	H ₂ O ₂ (μM)
1	11.5	988.5	100
2	500 of Tube #1	500	50
3	500 of Tube #2	500	25
4	500 of Tube #3	500	12.5
5	500 of Tube #4	500	6.25
6	500 of Tube #5	500	3.125
7	500 of Tube #6	500	1.56
8	500 of Tube #7	500	0.78
9	500 of Tube #8	500	0.39
10	500 of Tube #9	500	0.195
11	500 of Tube #10	500	0.098
12	0	500	0

Table 1. Preparation of H₂O₂ Standards

- Peroxidase Standard:** To prepare the peroxidase standards, first perform a 1:1000 dilution of the stock HRP in 1X Assay Buffer (e.g., Add 5 μL of HRP stock to 4.995 mL 1X Assay Buffer). Prepare only enough for immediate use. This solution has a concentration of 100 mU/mL. Use this 100 mU/mL solution to prepare standards in the concentration range of 0 mU/mL – 10 mU/mL by further diluting in 1X Assay Buffer (see Table 2 below). HRP diluted solutions and standards should be prepared fresh.

Standard Tubes	100 mU/mL HRP Standard (μL)	1X Assay Buffer (μL)	HRP (mU/mL)
1	100	900	10
2	500 of Tube #1	500	5
3	500 of Tube #2	500	2.5
4	500 of Tube #3	500	1.25
5	500 of Tube #4	500	0.625
6	500 of Tube #5	500	0.3125
7	500 of Tube #6	500	0.1563
8	0	500	0

Table 2. Preparation of HRP Standards

Assay Protocol

I. Hydrogen Peroxide

1. Prepare and mix all reagents thoroughly before use. Each sample, including unknowns and standards, should be assayed in duplicate or triplicate.
2. Add 50 μL of each sample (H₂O₂ standard, control or unknown) into an individual microtiter plate well.
3. Add 50 μL of ADHP/HRP Working Solution to each well. Mix the well contents thoroughly and incubate for 30 minutes at room temperature protected from light.

Note: This assay is continuous (not terminated) and therefore may be measured at multiple time points to follow the kinetics of the reactions.

4. Read the plate with a fluorescence microplate reader equipped for excitation in the 530-570 nm range and for emission in the 590-600 nm range.

5. Calculate the concentration of peroxide within samples by comparing the sample RFU to the standard curve. Subtract the value from the zero H_2O_2 control.

II. Peroxidase

1. Prepare and mix all reagents thoroughly before use. Each sample, including unknowns and standards, should be assayed in duplicate or triplicate.
2. Add 50 μL of each sample (HRP standard, control or unknown) into an individual microtiter plate well.
3. Add 50 μL of ADHP/ H_2O_2 Working Solution to each well. Mix the well contents thoroughly and incubate for 30 minutes at room temperature protected from light.
Note: This assay is continuous (not terminated) and therefore may be measured at multiple time points to follow the kinetics of the reactions.
4. Read the plate with a fluorescence microplate reader equipped for excitation in the 530-570 nm range and for emission in the 590-600 nm range.
5. Calculate the concentration of peroxidase within samples by comparing the sample RFU to the standard curve. Subtract the value from the zero HRP control.

Example of Results

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

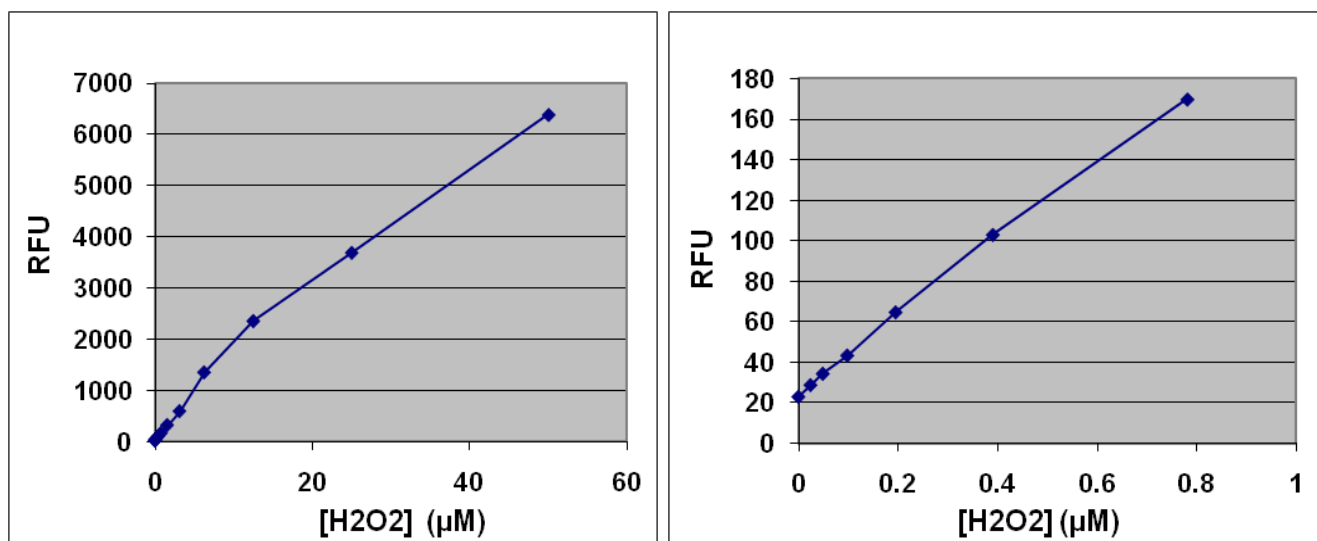


Figure 1. H_2O_2 Standard Curve.

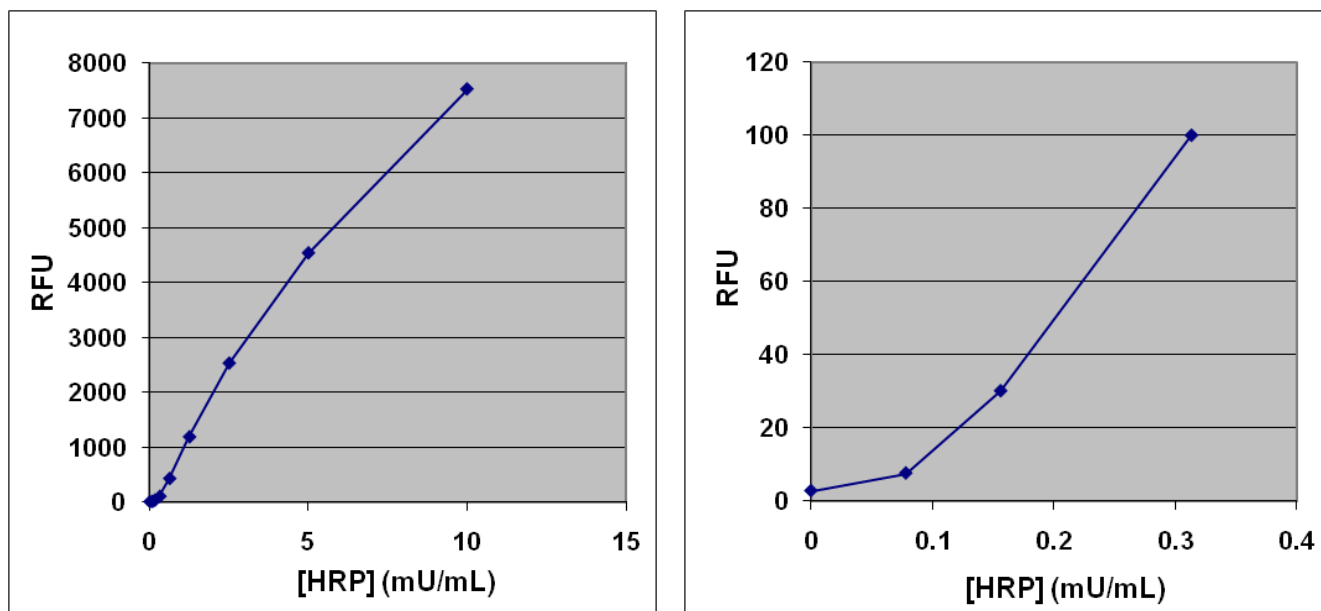


Figure 2. HRP Standard Curve.

References

1. Mohanty, J.G., et al. (1997) *J. Immunol. Methods* 202: 133.
2. Votyakova TV, and Reynolds IJ (2001) *Neurochem.* **79**:266.
3. Votyakova TV, and Reynolds IJ (2004) *Archives of Biochemistry and Biophysics.* **431**: 138-144.
4. Zhang, J., et al. (2001) *Antioxid. Redox Signal.* **3**: 493-504.

Recent Product Citations

1. Saenen, N.D. et al. (2023). Polystyrene Microplastics of Varying Sizes and Shapes Induce Distinct Redox and Mitochondrial Stress Responses in a Caco-2 Monolayer. *Antioxidants (Basel)*. **12**(3):739. doi: 10.3390/antiox12030739.
2. Alghabari, F. et al. (2023). Physio-Chemical and Agronomic-Based Characterization of Synthetic Hexaploid Wheat Germplasm under Field Imposed Conditions of Drought and Heat Stress. *Agronomy*. **13**(2):458. doi: 10.3390/agronomy13020458.
3. Damal Villivalam, S. et al. (2021). A necessary role of DNMT3A in endurance exercise by suppressing ALDH1L1-mediated oxidative stress. *EMBO J.* doi: 10.15252/embj.2020106491.
4. Kassem, S. et al. (2022). In vivo study of dose-dependent antioxidant efficacy of functionalized core-shell yttrium oxide nanoparticles. *Naunyn Schmiedebergs Arch Pharmacol.* doi: 10.1007/s00210-022-02219-1.
5. Han, G. et al. (2022). Nrf2 expands the intracellular pool of the chaperone AHSP in a cellular model of β -thalassemia. *Redox Biol.* **50**:102239. doi: 10.1016/j.redox.2022.102239.
6. Bononi, I. et al. (2022). Antioxidant Activity of Resveratrol Diastereomeric Forms Assayed in Fluorescent-Engineered Human Keratinocytes. *Antioxidants*. **11**(2):196. doi: 10.3390/antiox11020196.
7. Alghabari, F. et al. (2021). Biochemical and Physiological Responses of Thermostable Wheat Genotypes for Agronomic Yield under Heat Stress during Reproductive Stages. *Agronomy*. **11**(10):2080. doi: 10.3390/agronomy11102080.

8. Zunica, E.R.M. et al. (2021). Breast cancer growth and proliferation is suppressed by the mitochondrial targeted furazano[3,4-b]pyrazine BAM15. *Cancer Metab.* **9**(1):36. doi: 10.1186/s40170-021-00274-5.
9. Balmant, K.M. et al. (2021). Guard cell redox proteomics reveals a role of lipid transfer protein in plant defense. *J Proteomics.* doi: 10.1016/j.jprot.2021.104247.
10. 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.
11. Nitta, Y. et al. (2020). Catalase is required for peroxisome maintenance during adipogenesis. *Biochim Biophys Acta Mol Cell Biol Lipids.* doi: 10.1016/j.bbalip.2020.158726.
12. Ansar, M. et al. (2020). Increased Lung Catalase Activity Confers Protection Against Experimental RSV Infection. *Sci Rep.* **10**(1):3653. doi: 10.1038/s41598-020-60443-2.
13. Yoshimoto, S. et al. (2020). Riboflavin Plays a Pivotal Role in the UVA-Induced Cytotoxicity of Fibroblasts as a Key Molecule in the Production of H₂O₂ by UVA Radiation in Collaboration with Amino Acids and Vitamins. *Int J Mol Sci.* **21**(2). pii: E554. doi: 10.3390/ijms21020554.
14. Anand, S. et al. (2019). Agastache honey has superior antifungal activity in comparison with important commercial honeys. *Sci Rep.* **9**(1):18197. doi: 10.1038/s41598-019-54679-w.
15. Pei, J.F. et al. (2019). Diurnal oscillations of endogenous H₂O₂ sustained by p66Shc regulate circadian clocks. *Nat Cell Biol.* **21**(12):1553-1564. doi: 10.1038/s41556-019-0420-4.
16. Mehta, N. et al. (2019). Follistatin Protects against Glomerular Mesangial Cell Apoptosis and Oxidative Stress to Ameliorate Chronic Kidney Disease. *Antioxid Redox Signal.* doi: 10.1089/ars.2018.7684.
17. Lee, D.H. et al. (2019). Peroxiredoxin 6 confers protection against non-alcoholic fatty liver disease through maintaining mitochondrial function. *Antioxid Redox Signal.* doi: 10.1089/ars.2018.7544.
18. Nguyen, K. H. et al. (2018). Overexpression of GmNAC085 enhances drought tolerance in Arabidopsis by regulating glutathione biosynthesis, redox balance and glutathione-dependent detoxification of reactive oxygen species and methylglyoxal. *Environmental and Experimental Botany.* doi: 10.1016/j.envexpbot.2018.12.021.
19. Mu, H.N. et al. (2018). Caffeic acid attenuates rat liver injury after transplantation involving PDIA3-dependent regulation of NADPH oxidase. *Free Radic Biol Med.* **129**:202-214. doi: 10.1016/j.freeradbiomed.2018.09.009.
20. Bucekova, M. et al (2018). Microwave processing of honey negatively affects honey antibacterial activity by inactivation of bee-derived glucose oxidase and defensin-1. *Food Chemistry.* **240**: 1131-1136.
21. Badosa, E. (2017). Control of fire blight infections with synthetic peptides that elicit plant defense responses. *Journal of Plant Pathology.* **99** (Special issue), 65-73.
22. Ohta, Y., et al. (2017). Compound 48/80, a mast cell degranulator, causes oxidative damage by enhancing vitamin C synthesis via reduced glutathione depletion and lipid peroxidation through neutrophil infiltration in rat livers. *J. Clin. Biochem. Nutr.*
23. Endesfelder, S. (2017). Neuroprotection by Caffeine in Hyperoxia-Induced Neonatal Brain Injury. *Int J Mol Sci.* **18**(1). pii: E187. doi: 10.3390/ijms18010187.
24. Son, D.J. et al. (2016). Novel synthetic (E)-2-methoxy-4-(3-(4-methoxyphenyl) prop-1-en-1-yl) phenol inhibits arthritis by targeting signal transducer and activator of transcription 3. *Sci. Rep.* **6**:36852.

25. Begieneman, M.P.V. et al. (2016). Dopamine induces lipid accumulation, NADPH oxidase-related oxidative stress, and a proinflammatory status of the plasma membrane in H9c2 cells. *Am. J. Physiol. Heart Circ. Physiol.* **311**:H1097-H1107.
26. Ye, F. et al. (2016). High glucose induces reactivation of latent Kaposi's sarcoma-associated herpesvirus. *J Virol.* doi:10.1128/JVI.01049-16.
27. Horniackova, M. et al. (2016). Effect of gamma radiation on the antibacterial and antibiofilm activity of honeydew honey. *Eur Food Res Technol.* doi:10.1007/s00217-016-2725-x.
28. Michaels, D. L. et al. (2016). Cellular microbiology of *Mycoplasma canis*. *Infect Immun.* doi: 10.1128/IAI.01440-15.
29. Ruiz-Ojeda, F. J. et al. (2016). Impact of 3-amino-1, 2, 4-triazole (3-AT)-derived increase in hydrogen peroxide levels on inflammation and metabolism in human differentiated adipocytes. *PLoS One.* **11**: e0152550.
30. Douglas, D. N. et al. (2016). Oxidative stress attenuates lipid synthesis and increases mitochondrial fatty acid oxidation in hepatoma cells infected with hepatitis C virus. *J Biol Chem.* **291**:1974-1990.

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.
5628 Copley Drive
San Diego, CA 92111
Worldwide: +1 858 271-6500
USA Toll-Free: 1-888-CBL-0505
E-mail: tech@cellbiolabs.com
www.cellbiolabs.com

©2010-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.