
Product Manual

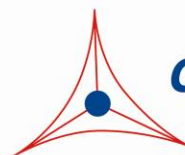
OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit (Colorimetric)

Catalog Number

STA-844

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 (H_2O_2), 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. The colorimetric probe reacts with H_2O_2 and horseradish peroxidase enzyme (HRP) to produce a pink colored product. The probe has less background and greater stability than the commonly used Xylenol Orange (FOX) colorimetric assay for H_2O_2 . The probe can be also used as an ultrasensitive assay for peroxidase activity when H_2O_2 is in excess. The kit has a detection sensitivity limit of $0.8 \mu\text{M}$ (H_2O_2) or 0.16 mU/mL (peroxidase). Each kit provides sufficient reagents to perform up to 500 assays, including standard curve and unknown samples.

Assay Principle

The OxiSelect™ Hydrogen Peroxide/Peroxidase Assay Kit is a sensitive quantitative colorimetric assay for hydrogen peroxide or peroxidase. In the presence of peroxidase, the probe reacts with H_2O_2 in a 1:1 stoichiometry to produce a bright pink colored product. This product can be easily read by a standard colorimetric microplate reader with a filter in the 540-570 nm range. Absorbance values are proportional to the H_2O_2 or peroxidase levels within the samples, depending on the assay employed. The H_2O_2 or peroxidase content in unknown samples is determined by comparison with its respective standard curve.

Related Products

1. STA-330: OxiSelect™ TBARS Assay Kit (MDA Quantitation)
2. STA-342: OxiSelect™ Intracellular ROS Assay Kit (Green Fluorescence)
3. STA-347: OxiSelect™ In Vitro ROS/RNS Assay Kit (Green Fluorescence)
4. STA-832: OxiSelect™ MDA Adduct Competitive ELISA Kit
5. STA-838: OxiSelect™ HNE Adduct Competitive ELISA Kit

Kit Components (shipped at room temperature)

1. Colorimetric Probe (100X) (Part No. 284401): One 250 µL amber tube of solution.
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. 10 µL to 1000 µL adjustable single channel micropipettes with disposable tips
4. 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
5. Standard 96-well microtiter plate
6. Multichannel micropipette reservoir
7. Spectrophotometric microplate reader capable of reading in the 540-570 nm absorbance range.
8. Superoxide dismutase (optional)

Storage

Upon receipt, aliquot and store the Colorimetric Probe and HRP at -20°C. Avoid multiple freeze/thaw cycles. Store the remaining kit components at 4°C. The Colorimetric Probe is light sensitive and 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.
- Hydrogen Peroxide Working Solution (Hydrogen Peroxide Assay): If measuring hydrogen peroxide, prepare a working solution by diluting the Colorimetric Probe 1:100 and HRP to a final

concentration of 0.2 U/mL in 1X Assay Buffer (eg. Add 50 μ L Colorimetric Probe stock solution and 10 μ L HRP stock solution to 4.940 mL 1X Assay Buffer). This volume is enough for ~100 assays. The Hydrogen Peroxide Working Solution should be protected from light and used within 4 hours. Prepare only enough for immediate use.

- Peroxidase Working Solution (Peroxidase Assay): If measuring peroxidases, prepare a working solution by diluting the Colorimetric Probe 1:100 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 μ L 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 Probe/1X Assay Buffer (eg. Add 50 μ L Colorimetric Probe 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 Peroxidase Working Solution should be protected from light and used within 4 hours. 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 probe 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 ($\geq 500 \mu$ M final concentration) or peroxidase (≥ 100 mU/mL) can lower the absorbance because excess H₂O₂ or peroxidase can further oxidize the reaction product.*
- *Samples with NADH concentrations above 10 μ M and glutathione concentrations above 50 μ M will oxidize the 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 (Tatyana et al, Ref. 2).*
- *Avoid samples containing DTT or β -mercaptoethanol since the probe is not stable in the presence of thiols (above 10 μ M).*

Preparation of Standard Curves

- **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. 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 (e.g., Add 11.5 µL of H₂O₂ to 988.5 µL 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	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	0.63
6	500 of Tube #5	500	0.32
7	500 of Tube #6	500	0.16
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 (H_2O_2 standard, control or unknown) into an individual microtiter plate well.
3. Add 50 μL of Hydrogen Peroxide 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 absorbance with a microplate reader in the 540-570 nm range.
5. Calculate the concentration of peroxide within samples by comparing the sample absorbance 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 Peroxidase 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 absorbance with a microplate reader in the 540-570 nm range.
5. Calculate the concentration of peroxidase within samples by comparing the sample absorbance 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.

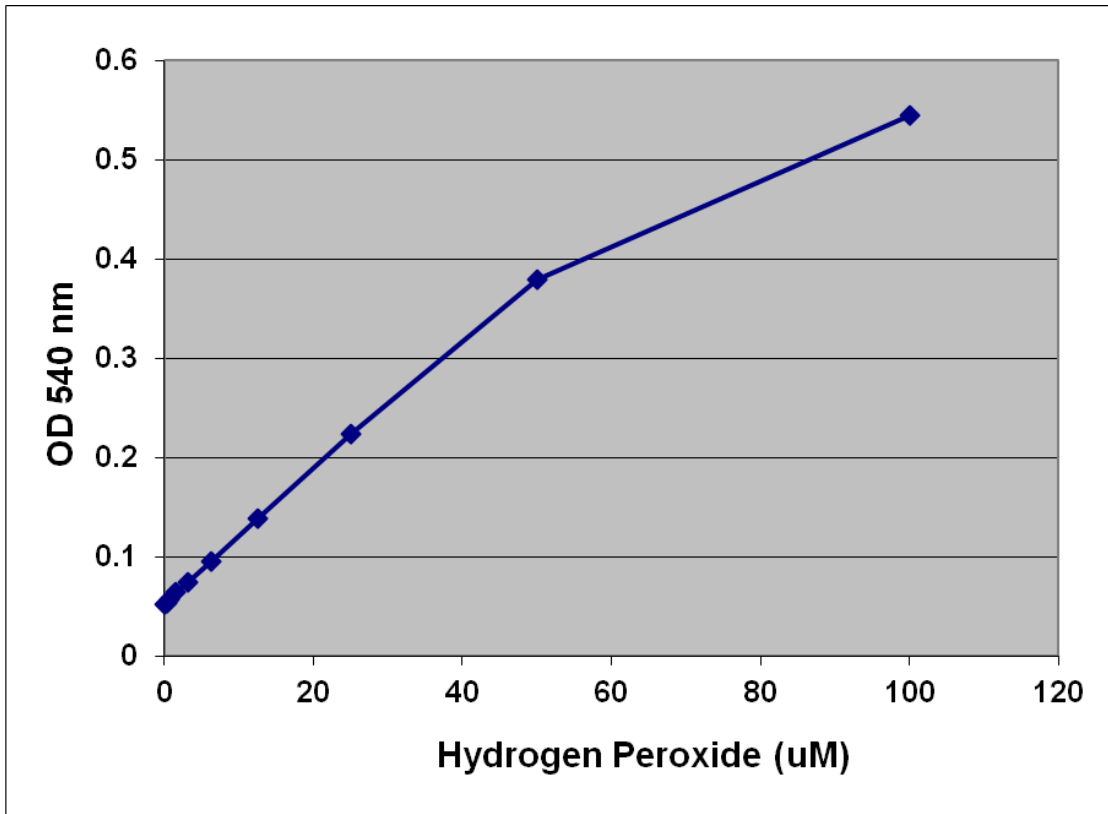


Figure 1. H₂O₂ Standard Curve.

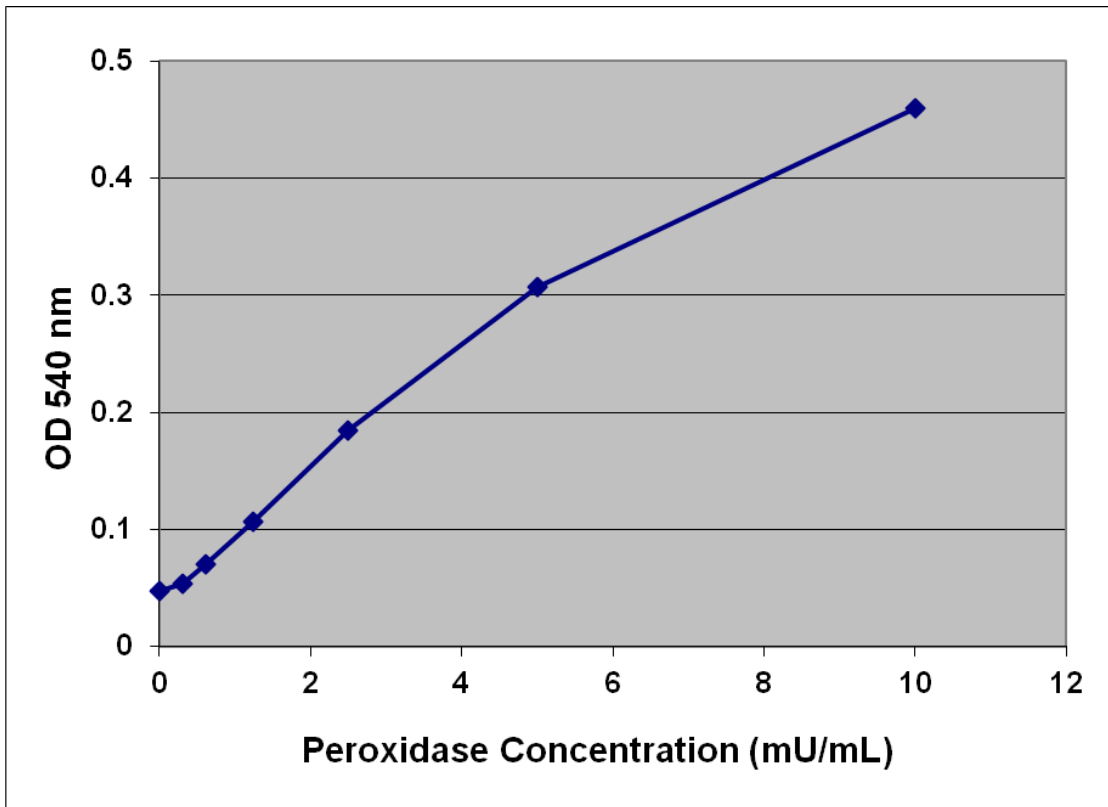


Figure 2. HRP Standard Curve.

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.

Recent Product Citations

1. Miliński, M. et al. (2023). The effect of sulindac on redox homeostasis and apoptosis-related proteins in melanotic and amelanotic cells. *Pharmacol Rep*. doi: 10.1007/s43440-023-00493-1.
2. Coudert, E. et al. (2023). Slow and Fast-Growing Chickens Use Different Antioxidant Pathways to Maintain Their Redox Balance during Postnatal Growth. *Animals (Basel)*. **13**(7):1160. doi: 10.3390/ani13071160.
3. de Oliveira, M.G. et al. (2022). Selective Pharmacological Inhibition of NOX2 by GSK2795039 Improves Bladder Dysfunction in Cyclophosphamide-Induced Cystitis in Mice. *Antioxidants (Basel)*. **12**(1):92. doi: 10.3390/antiox12010092.
4. Ohmori, I. et al. (2022). Thioredoxin deficiency increases oxidative stress and causes bilateral symmetrical degeneration in rat midbrain. *Neurobiol Dis*. doi: 10.1016/j.nbd.2022.105921.
5. Abdrabouh, A.E. (2022). Inflammatory and proapoptotic effects of inhaling gasoline fumes on the lung and ameliorative effects of fenugreek seeds. *Sci Rep*. **12**(1):14446. doi: 10.1038/s41598-022-18607-9.
6. Park, H.J. et al. (2022). Doxorubicin Induces Bone Loss by Increasing Autophagy through a Mitochondrial ROS/TRPML1/TFEB Axis in Osteoclasts. *Antioxidants (Basel)*. **11**(8):1476. doi: 10.3390/antiox11081476.
7. Sul, O.J. et al. (2022). GSPE Protects against Bleomycin-Induced Pulmonary Fibrosis in Mice via Ameliorating Epithelial Apoptosis through Inhibition of Oxidative Stress. *Oxid Med Cell Longev*. doi: 10.1155/2022/8200189.
8. Landerer, S. et al. (2021). UDP-glucuronosyltransferases mediate coffee-associated reduction of liver fibrosis in bile duct ligated humanized transgenic UGT1A mice. *Hepatobiliary Surg Nutr*. **10**(6):766-781. doi: 10.21037/hbsn-20-9.
9. Odatsu, T. et al. (2021). Lactoferrin with Zn-ion protects and recovers fibroblast from H2O2-induced oxidative damage. *Int J Biol Macromol*. **190**:368-374. doi: 10.1016/j.ijbiomac.2021.08.214.
10. Graham, R.J. et al. (2021). Zinc supplementation modulates intracellular metal uptake and oxidative stress defense mechanisms in CHO cell cultures. *Biochem Eng J*. doi: 10.1016/j.bej.2021.107928.
11. Gromotowicz-Poplawska, A. et al. (2021). Hyperglycemia Potentiates Prothrombotic Effect of Aldosterone in a Rat Arterial Thrombosis Model. *Cells*. **10**(2):471. doi: 10.3390/cells10020471.
12. Chiang, C.C. et al. (2020). Aqueous Extract of Kan-Lu-Hsiao-Tu-Tan Ameliorates Collagen-Induced Arthritis in Mice by Inhibiting Oxidative Stress and Inflammatory Responses. *Life (Basel)*. **10**(12):E313. doi: 10.3390/life10120313.
13. Sengupta, S. et al. (2020). Ligand-induced gene activation is associated with oxidative genome damage whose repair is required for transcription. *Proc Natl Acad Sci U S A*. doi: 10.1073/pnas.1919445117.
14. Böhm, A. et al. (2020). Increased mitochondrial respiration of adipocytes from metabolically unhealthy obese compared to healthy obese individuals. *Sci Rep*. **10**(1):12407. doi: 10.1038/s41598-020-69016-9.

15. Sabry, M.M. et al. (2020). Carnitine, apelin and resveratrol regulate mitochondrial quality control (QC) related proteins and ameliorate acute kidney injury: role of hydrogen peroxide. *Arch Physiol Biochem*. doi: 10.1080/13813455.2020.1773504.
16. Ścibior, A. et al. (2020). In vitro effect of vanadyl sulfate on cultured primary astrocytes: cell viability and oxidative stress markers. *J Appl Toxicol*. doi: 10.1002/jat.3939.
17. El-Boshy, M. et al. (2020). Vitamin D3 and calcium cosupplementation alleviates cadmium hepatotoxicity in the rat: Enhanced antioxidative and anti-inflammatory actions by remodeling cellular calcium pathways. *J Biochem Mol Toxicol*. doi: 10.1002/jbt.22440.
18. Oncul, S. et al. (2019). A Kojic Acid Derivative Promotes Intrinsic Apoptotic Pathway of Hepatocellular Carcinoma Cells without Incurring Drug Resistance. *Chem Biol Drug Des*. doi: 10.1111/cbdd.13615.
19. Nishu, S.D. et al. (2019). Nutritional status regulates algicidal activity of *Aeromonas* sp. L23 against cyanobacteria and green algae. *PLoS One*. **14**(3):e0213370. doi: 10.1371/journal.pone.0213370.
20. Sakai, T. et al. (2019). Cytoplasmic OH scavenger TA293 attenuates cellular senescence and fibrosis by activating macrophages through oxidized phospholipids/TLR4. *Life Sci*. **221**:284-292. doi: 10.1016/j.lfs.2019.02.038.
21. Abdrabouh, A.E. (2019). Liver disorders related to exposure to gasoline fumes in male rats and role of fenugreek seed supplementation. *Environ Sci Pollut Res Int*. **26**(9):8949-8957. doi: 10.1007/s11356-019-04307-x.
22. Sakai, T. et al. (2019). Hydrogen Indirectly Suppresses Increases in Hydrogen Peroxide in Cytoplasmic Hydroxyl Radical-Induced Cells and Suppresses Cellular Senescence. *Int J Mol Sci*. **20**(2). pii: E456. doi: 10.3390/ijms20020456.
23. Gamal, M. et al. (2018). Possible involvement of tetrahydrobiopterin in the disturbance of redox homeostasis in sepsis - Induced brain dysfunction. *Brain Res*. **1685**:19-28. doi: 10.1016/j.brainres.2018.02.008.
24. Otreba, M. et al (2017). Prochlorperazine interaction with melanin and melanocytes. *Die Pharmazie - An International Journal of Pharmaceutical Sciences*. **72** (3): 171-176(6).
25. Lee, S.H. et al. (2016). Proteomic analysis indicates activation of reactive oxygen species signaling during seed germination and seedlings growth in *Hordeum vulgare* (barley). *J. Prot. Proteom*. **7**:269-277.

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

©2015-2023: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.