
Product Manual

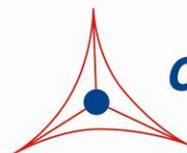
Total Cholesterol Assay Kit (Colorimetric)

Catalog Number

STA-384

192 assays

FOR RESEARCH USE ONLY
Not for use in diagnostic procedures



CELL BIOLABS, INC.

Creating Solutions for Life Science Research

Introduction

Cholesterol is a lipid sterol that is produced in and transported throughout the bloodstream in eukaryotes. Cholesterol is a critical compound used in the structure of cell membranes, hormones, and cell signaling. It is an essential component of animal cell structure in order to maintain permeability and fluidity. Cholesterol is a precursor for steroid hormones including the adrenal gland hormones cortisol and aldosterone, sex hormones progesterone, estrogens, and testosterone, and bile acids and vitamin D. Cholesterol is transported throughout the body within lipoproteins, which have cell-specific signals that direct the lipids they transport to certain tissues. For this reason, lipoproteins exist in different forms within the blood based on their density. These include chylomicrons, very-low density lipoproteins (VLDLs), low-density lipoproteins (LDLs), intermediate-density lipoproteins (IDLs), and high-density lipoproteins (HDLs). The higher the lipid content within a lipoprotein, the lower its density. Cholesterol exists within a lipoprotein as a free alcohol and as a fatty cholesteryl ester, which is the predominant form of cholesterol transport and storage.

Determining circulatory levels of lipoproteins is critical to the diagnosis of lipid transport disorders. High levels of cholesterol and cholesteryl esters (hypercholesterolemia) have been associated with cardiovascular disease such as atherosclerosis and heart disease, although lower levels (hypocholesterolemia) may be associated with cancer, depression, or respiratory diseases.

Cell Biolabs' Total Cholesterol Assay Kit is a simple colorimetric assay that measures the amount of total cholesterol present in plasma, serum, tissue homogenates, or cell lysates in a 96-well microtiter plate format. The assay will detect total cholesterol (cholesteryl esters plus free cholesterol) in the presence of cholesterol esterase or only free cholesterol in the absence of the esterase enzyme. Each kit provides sufficient reagents to perform up to 192 assays, including blanks, cholesterol standards and unknown samples. Sample cholesterol concentrations are determined by comparison with a known cholesterol standard. Cholesteryl esters can be quantified by subtracting the free cholesterol values from the total cholesterol value.

Assay Principle

Cell Biolabs' Total Cholesterol Assay Kit measures the total cholesterol within serum, plasma, lysate, or tissue samples. The assay is based on the enzyme driven reaction that quantifies both cholesterol esters and free cholesterol. Cholesterol esters are hydrolyzed via cholesterol esterase into cholesterol, which is then oxidized by cholesterol oxidase into the ketone cholest-4-en-3-one plus hydrogen peroxide. The hydrogen peroxide is then detected with a highly specific colorimetric probe. Horseradish peroxidase catalyzes the reaction between the probe and hydrogen peroxide, which bind in a 1:1 ratio. Samples are compared to a known concentration of cholesterol standard in a 96-well microtiter plate format. Samples and standards are incubated for 45 minutes and then read with a standard 96-well colorimetric plate reader (Figure 1).

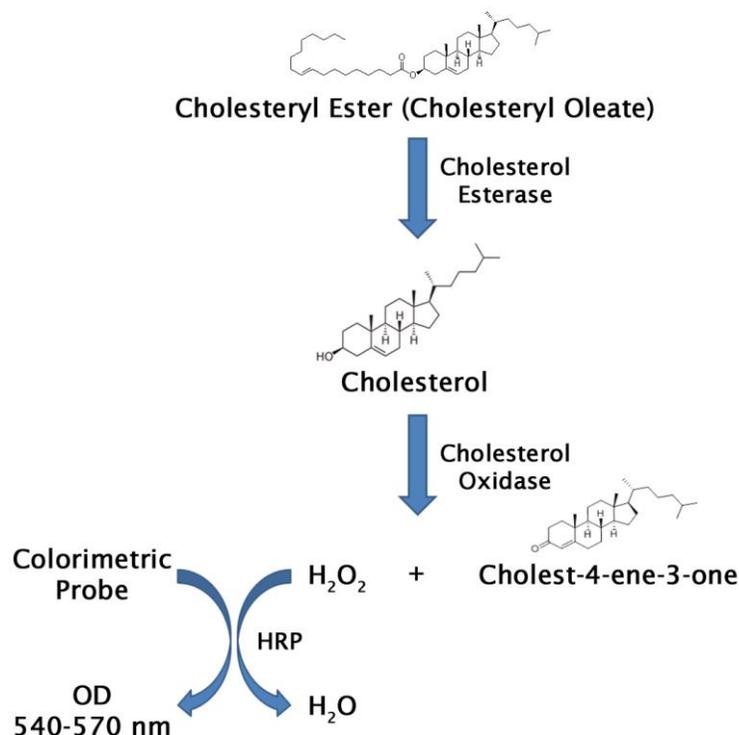


Figure 1. Colorimetric Cholesterol Assay Principle

Related Products

1. STA-367: Human ApoE ELISA Kit
2. STA-369: OxiSelect™ Human Oxidized LDL ELISA Kit (MDA-LDL Quantitation)
3. STA-390: Total Cholesterol Assay Kit (Fluorometric)
4. STA-391: HDL and LDL/VLDL Cholesterol Assay Kit

Kit Components

Box 1 (shipped at room temperature)

1. Cholesterol Standard (Part No. 239001): One 50 µL tube of a 10 mM cholesterol solution in ethanol.
2. Assay Diluent (5X) (Part No. 239002): One 100 mL bottle.
3. 50X Colorimetric Probe (Part No. 238401): One 200 µL tube in DMSO.
4. HRP (Part No. 234402): Two 100 µL tubes of 100 U/mL HRP solution in glycerol.

Box 2 (shipped on blue ice packs)

1. Cholesterol Esterase (Part No. 239003): One tube of 10 Units enzyme in powder.
2. Cholesterol Oxidase (Part No. 239004): One 200 μ L tube.

Materials Not Supplied

1. 96-well microtiter plates
2. Distilled or deionized water
3. 1X PBS
4. 10 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
5. 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
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, store the Assay Diluent at 4°C. Store the remaining kit components at -20°C. The 50X Colorimetric Probe is light sensitive and must be stored accordingly. Avoid multiple freeze/thaw cycles.

Preparation of Reagents

- 1X Assay Diluent: Warm the Assay Diluent (5X) to room temperature prior to using. Dilute the Assay Diluent (5X) with deionized water by diluting the 100 mL Diluent with 400 mL deionized water for 500 mL total. Mix to homogeneity. Store the 1X Assay Diluent at 4°C up to six months.
- Cholesterol Esterase: Reconstitute the powder with 200 μ L of 1X Assay Diluent. Vortex vigorously until dissolved. Prepare aliquots and store at -20°C to avoid multiple freeze thaws of the reconstituted powder.
- Cholesterol Reaction Reagent: Prepare the reagent by diluting the Cholesterol Oxidase 1:50, HRP 1:50, Colorimetric Probe 1:50, and Cholesterol Esterase 1:250 in 1X Assay Diluent. (e.g., For 100 assays, combine 100 μ L of Cholesterol Oxidase, 100 μ L of HRP, 100 μ L Colorimetric Probe, and 20 μ L Cholesterol Esterase with 1X Assay Diluent to 5 mL total solution). Mix thoroughly and protect the solution from light. For best results, place the Cholesterol Reaction Reagent on ice and use within 30 minutes of preparation. Do not store the Cholesterol Reaction Reagent solution.

Notes:

1. *If testing for the concentration of free cholesterol is needed only, omit the addition of Cholesterol Esterase from the Cholesterol Reaction Reagent solution.*
2. *The Colorimetric Probe is light sensitive and must be stored accordingly.*

Preparation of Samples

Samples should be assayed immediately or stored at -80°C prior to performing the assay. Optimal experimental conditions for samples must be determined by the investigator. The following

recommendations are only guidelines and may be altered to optimize or complement the user's experimental design. A set of serial dilutions is recommended for samples to achieve optimal assay results and minimize possible interfering compounds. Run proper controls as necessary. Always run a standard curve with samples.

- **Tissue Lysates:** For 10 mg of tissue, extract with 200 μ L of a mixture of chloroform: isopropanol: NP-40 (7:11:0.1) in a micro-homogenizer. Centrifuge the extract 10 minutes at 15,000 x g. Transfer the liquid (organic phase) to a new tube, taking care to avoid the pellet. Air dry at 50°C to remove the chloroform. Put samples under vacuum for 30 minutes to remove the trace amounts of organic solvent. Dissolve the dried lipids in 200 μ L of 1X Assay Diluent with sonicating and vortexing until the solution is homogenous (the solution may appear cloudy). This extraction procedure may be scaled up if larger sample amounts are desired. Use 1 - 50 μ L of extracted sample per assay. Next, adjust the volume to 50 μ L per well with 1X Assay Diluent. For unknown samples, we suggest testing different amounts of samples to ensure that the readings are within the linear portion of the standard curve.
- **Cell Lysates:** Wash cells 3 times with cold PBS prior to lysis. For 10^6 cells, extract with 200 μ L of a mixture of chloroform: isopropanol: NP-40 (7:11:0.1) in a micro-homogenizer. Centrifuge the extract 10 minutes at 15,000 x g. Transfer the liquid (organic phase) to a new tube, taking care to avoid the pellet. Air dry at 50°C to remove the chloroform. Put samples under vacuum for 30 minutes to remove the trace amounts of organic solvent. Dissolve the dried lipids in 200 μ L of 1X Assay Diluent with sonicating and vortexing until the solution is homogenous (the solution may appear cloudy). This extraction procedure may be scaled up if larger sample amounts are desired. Use 1 - 50 μ L of extracted sample per assay. Next, adjust the volume to 50 μ L per well with 1X Assay Diluent. For unknown samples, we suggest testing different amounts of samples to ensure that the readings are within the linear portion of the standard curve.
- **Serum:** Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the serum layer and store on ice. Avoid disturbing the white buffy layer. Aliquot samples for testing and store at -80°C. Perform dilutions in 1X Assay Diluent. Serum samples must be diluted at least 1:100 to 1:200 with Assay Diluent. This will provide values within the range of the standard curve. Cholesterol levels in serum average about 3% higher in value than in the corresponding plasma pair (Ref. 2).
- **Plasma:** Avoid hemolyzed and lipemic blood samples. Collect blood with heparin or citrate and centrifuge at 2000 x g and 4°C for 10 minutes. Remove the plasma layer and store on ice. Avoid disturbing the white buffy layer. Aliquot samples for testing and store at -80°C. Perform dilutions in 1X Assay Diluent. Plasma samples must be diluted at least 1:100 to 1:200 with Assay Diluent. This will provide values within the range of the standard curve.

Notes:

1. *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.*
2. *Avoid samples containing DTT or β -mercaptoethanol since the colorimetric probe is not stable in the presence of thiols (above 10 μ M).*

Preparation of Cholesterol Standard Curve

1. Prepare fresh cholesterol standards before use by diluting in 1X Assay Diluent. First, dilute the stock Cholesterol Standard 10 mM solution 1:40 in 1X Assay Diluent for a 250 μ M solution. (eg. add 25 μ L of the stock 10 mM standard to 975 μ L of 1X Assay Diluent). Vortex thoroughly. Use the diluted Cholesterol Standards promptly.
2. Use this 250 μ M solution to prepare a series of the remaining cholesterol standards according to Table 1 below.

Tubes	250 μM Cholesterol Standard (μL)	1X Assay Diluent (μL)	Resulting Cholesterol Concentration (μM)
1	1000	0	250
2	500 of Tube #1	500	125
3	500 of Tube #2	500	62.5
4	500 of Tube #3	500	31.3
5	500 of Tube #4	500	15.6
6	500 of Tube #5	500	7.8
7	500 of Tube #6	500	3.9
8	500 of Tube #7	500	1.9
9	500 of Tube #8	500	1.0
10	0	1000	0

Table 1. Preparation of Cholesterol Standards.

Note: Do not store diluted cholesterol standard solutions.

Assay Protocol

Each cholesterol standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

1. Add 50 μ L of the diluted cholesterol standards or samples to a 96-well microtiter plate.
2. Add 50 μ L of the prepared Cholesterol Reaction Reagent to each well and mix the well contents thoroughly.
3. Cover the plate wells to protect the reaction from light. Incubate the plate for 45 minutes at 37°C.
4. IMMEDIATELY read the plate with a spectrophotometric microplate reader in the 540-570 nm range.
5. Calculate the concentration of cholesterol within samples by comparing the sample absorbance values to the cholesterol standard curve.

Example of Results

The following figures demonstrate typical Total Cholesterol Assay results. One should use the data below for reference only. This data should not be used to interpret or calculate actual sample results.

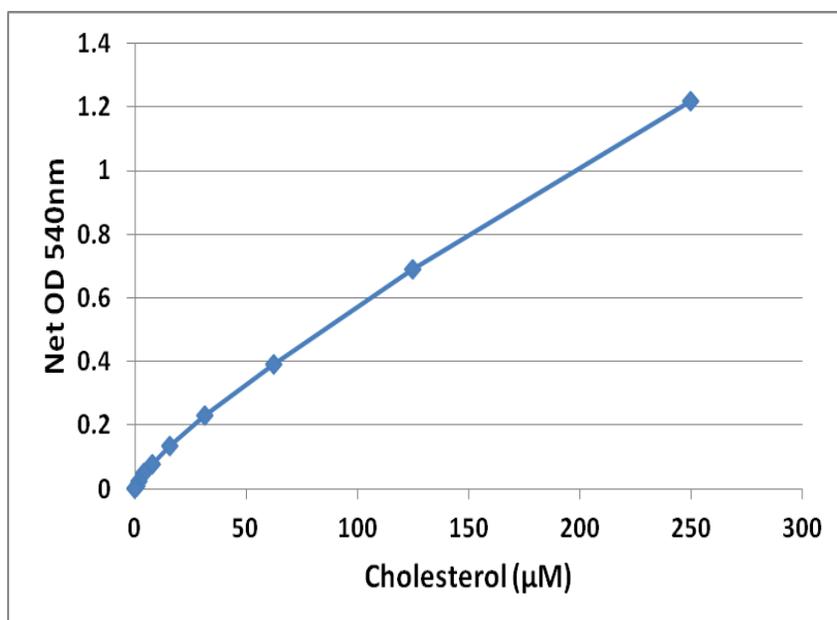


Figure 2: Cholesterol Standard Curve.

Calculation of Results

1. Calculate the average absorbance values for every standard, control, and sample. Subtract the average zero standard value from itself and all standard and sample values. This is the corrected absorbance.
2. Plot the corrected absorbance for the standards against the final concentration of the cholesterol standards from Table 1 to determine the best curve. See Figure 2 for an example standard curve.
3. Determine the cholesterol concentration of the samples with the equation obtained from the linear regression analysis of the standard curve. Substitute the corrected absorbance values for each sample. Remember to account for dilution factors.

$$\text{Total Cholesterol } (\mu\text{M}) = \left[\frac{\text{Sample corrected absorbance}}{\text{Slope}} \right] \times \text{Sample dilution}$$

$$\text{Cholesteryl Ester } (\mu\text{M}) = \text{Total Cholesterol} - \text{Free Cholesterol}$$

Note: For the conversion of results from µM to mg/dl, divide the cholesterol concentration (µM) by 25.9.

References

1. Admundson, D.M., et al. (1999) *J. Biochem. Biophys. Meth.* **38**: 43-52.
2. Cholesterol and Triglyceride concentrations in Serum/Plasma Pairs. (1977) *Clin. Chem.* **23**: 60-63.
3. Fossati, P., et al. (1982) *Clin. Chem.* **28**: 2077-2080.
4. Ledwozyw, A., et al. (1986) *Clin. Chim. Acta.* **155**: 275-284.
5. Lee, S.M. et al. (2008) *Lipids* **43**: 419-429.

Recent Product Citations

1. Tashiro, J. et al. (2023). CYP11A1 silencing suppresses HMGCR expression via cholesterol accumulation and sensitizes CRPC cell line DU-145 to atorvastatin. *J Pharmacol Sci.* **153**(3):104-112. doi: 10.1016/j.jphs.2023.08.002.
2. Alhelf, M. et al. (2023). Long noncoding RNA (taurine upregulated gene 1) and micro RNA-377: emerging players in the development of metabolic syndrome among psoriasis patients. *Beni-Suef Univ J Basic Appl Sci.* **12**(68). doi: 10.1186/s43088-023-00404-6.
3. Nopparat, J. et al. (2023). Probiotics of *Lacticaseibacillus paracasei* SD1 and *Lacticaseibacillus rhamnosus* SD11 attenuate inflammation and β -cell death in streptozotocin-induced type 1 diabetic mice. *PLoS One.* **18**(4):e0284303. doi: 10.1371/journal.pone.0284303.
4. Yang, X. et al. (2023). Sprouty1 has a protective role in atherogenesis and modifies the migratory and inflammatory phenotype of vascular smooth muscle cells. *Atherosclerosis.* **373**:17-28. doi: 10.1016/j.atherosclerosis.2023.04.007.
5. Mou, Y. et al. (2023). Chenodeoxycholic acid rescues axonal degeneration in induced pluripotent stem cell-derived neurons from spastic paraplegia type 5 and cerebrotendinous xanthomatosis patients. *Orphanet J Rare Dis.* **18**(1):72. doi: 10.1186/s13023-023-02666-w.
6. Bashir, K.M.I. et al. (2023). Efficacy Confirmation Test of Black Cumin (*Nigella sativa* L.) Seeds Extract Using a High-Fat Diet Mouse Model. *Metabolites.* **13**(4):501. doi: 10.3390/metabo13040501.
7. Abdelwahed, K.S. et al. (2023). Pseurotin A Validation as a Metastatic Castration-Resistant Prostate Cancer Recurrence-Suppressing Lead via PCSK9-LDLR Axis Modulation. *Mar Drugs.* **21**(4):215. doi: 10.3390/md21040215.
8. Fahrner, A. et al. (2023). microRNA-501 controls myogenin+/CD74+ myogenic progenitor cells during muscle regeneration. *Mol Metab.* **71**:101704. doi: 10.1016/j.molmet.2023.101704.
9. Lin, B. et al. (2023) Vitamin E Supplement Protects Against Gestational Diabetes Mellitus in Mice Through nuclear factor-erythroid factor 2-related factor 2/heme oxygenase-1 Signaling Pathway. *Diabetes Metab Syndr Obes.* **16**:565-574. doi: 10.2147/DMSO.S397255.
10. Begemann, K. et al. (2023). Rest phase snacking increases energy resorption and weight gain in male mice. *Mol Metab.* **69**:101691. doi: 10.1016/j.molmet.2023.101691.
11. Bedsted, A.E. et al. (2023). Detection of Porcine Deltacoronavirus RNA in the Upper and Lower Respiratory Tract and Biliary Fluid and the Effect of Infection on Serum Cholesterol Levels and Blood T Cell Population Frequencies in Gnotobiotic Piglets. *Vet Sci.* **10**(2):117. doi: 10.3390/vetsci10020117.
12. Landowski, M. et al. (2023). Transmembrane protein 135 regulates lipid homeostasis through its role in peroxisomal DHA metabolism. *Commun Biol.* **6**(1):8. doi: 10.1038/s42003-022-04404-7.
13. Moreau, F. et al. (2023). Liver-specific FGFR4 knockdown in mice on an HFD increases bile acid synthesis and improves hepatic steatosis. *J Lipid Res.* **64**(2):100324. doi: 10.1016/j.jlr.2022.100324.
14. Pathak, R. et al. (2022). Prolonged effects of DPP-4 inhibitors on steato-hepatic changes in Sprague-Dawley rats fed a high-cholesterol diet. *Inflamm Res.* doi: 10.1007/s00011-022-01572-4.
15. Bhat, N. et al. (2022). TCF7L2 transcriptionally regulates Fgf15 to maintain bile acid and lipid homeostasis through gut-liver crosstalk. *FASEB J.* **36**(3):e22185. doi: 10.1096/fj.202101607R.
16. Ranasinghe, N. et al. (2022). Cholesterol Accumulation in Livers of Indian Medaka, *Oryzias dancena*, Acclimated to Fresh Water and Seawater. *Front. Mar. Sci.* doi: 10.3389/fmars.2022.891706.

17. Sugiyama, T. et al. (2022). Chemical chaperones ameliorate neurodegenerative disorders in Derlin-1-deficient mice via improvement of cholesterol biosynthesis. *Sci Rep.* **12**(1):21840. doi: 10.1038/s41598-022-26370-0.
18. Pasello, M. et al. (2022). ABCA6 affects the malignancy of Ewing sarcoma cells via cholesterol-guided inhibition of the IGF1R/AKT/MDM2 axis. *Cell Oncol (Dordr).* doi: 10.1007/s13402-022-00713-5.
19. Shang, Y. et al. (2022). A CHCHD6-APP axis connects amyloid and mitochondrial pathology in Alzheimer's disease. *Acta Neuropathol.* doi: 10.1007/s00401-022-02499-0.
20. Hwang, S.M. et al. (2022). Preventive and Therapeutic Effects of Krill Oil on Obesity and Obesity-Induced Metabolic Syndromes in High-Fat Diet-Fed Mice. *Mar Drugs.* **20**(8):483. doi: 10.3390/md20080483.
21. Cho, I.J. et al. (2022). Lemon Balm and Corn Silk Mixture Alleviates Metabolic Disorders Caused by a High-Fat Diet. *Antioxidants (Basel).* **11**(4):730. doi: 10.3390/antiox11040730.
22. Choi, J.Y. et al. (2022). Combination Effects of Metformin and a Mixture of Lemon Balm and Dandelion on High-Fat Diet-Induced Metabolic Alterations in Mice. *Antioxidants (Basel).* **11**(3):580. doi: 10.3390/antiox11030580.
23. Zhao, Y. et al. (2022). ATAD3A oligomerization promotes neuropathology and cognitive deficits in Alzheimer's disease models. *Nat Commun.* **13**(1):1121. doi: 10.1038/s41467-022-28769-9.
24. Alsabaani, N.A. et al. (2022). Maslinic Acid Protects against Streptozotocin-Induced Diabetic Retinopathy by Activating Nrf2 and Suppressing NF- κ B. *J Ophthalmol.* **2022**:3044202. doi: 10.1155/2022/3044202.
25. Chen, T.Y. et al. (2022). Effects of the Water Extract of Fermented Rice Bran on Liver Damage and Intestinal Injury in Aged Rats with High-Fat Diet Feeding. *Plants (Basel).* **11**(5):607. doi: 10.3390/plants11050607.
26. Valdivia, A.O. & Bhattacharya, S.K. (2022). Lyso-Lipid-Induced Oligodendrocyte Maturation Underlies Restoration of Optic Nerve Function. *eNeuro.* **9**(1): ENEURO.0429-21.2022. doi: 10.1523/ENEURO.0429-21.2022.
27. Landowski, M. et al. (2022). A mutation in transmembrane protein 135 impairs lipid metabolism in mouse eyecups. *Sci Rep.* **12**(1):756. doi: 10.1038/s41598-021-04644-3.
28. Fitzpatrick, A.M. et al. (2021). Obesity Is Associated with Sustained Symptomatology and Unique Inflammatory Features in Children with Asthma. *J Allergy Clin Immunol Pract.* doi: 10.1016/j.jaip.2021.10.020.
29. Fujimoto, N. et al. (2021). Interaction of galectin-7 with HMGCS1 in vitro may facilitate cholesterol deposition in cultured keratinocytes. *J Invest Dermatol.* doi: 10.1016/j.jid.2021.04.038.
30. Ruano-Gallego, D. et al. (2021). Type III secretion system effectors form robust and flexible intracellular virulence networks. *Science.* **371**(6534):eabc9531. doi: 10.1126/science.abc9531.

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's 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

©2012-2024: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.