

---

**Product Manual**

# **Total Cholesterol Assay Kit (Fluorometric)**

**Catalog Number**

**STA-390**

**192 assays**

**FOR RESEARCH USE ONLY  
Not for use in diagnostic procedures**

---

## **Introduction**

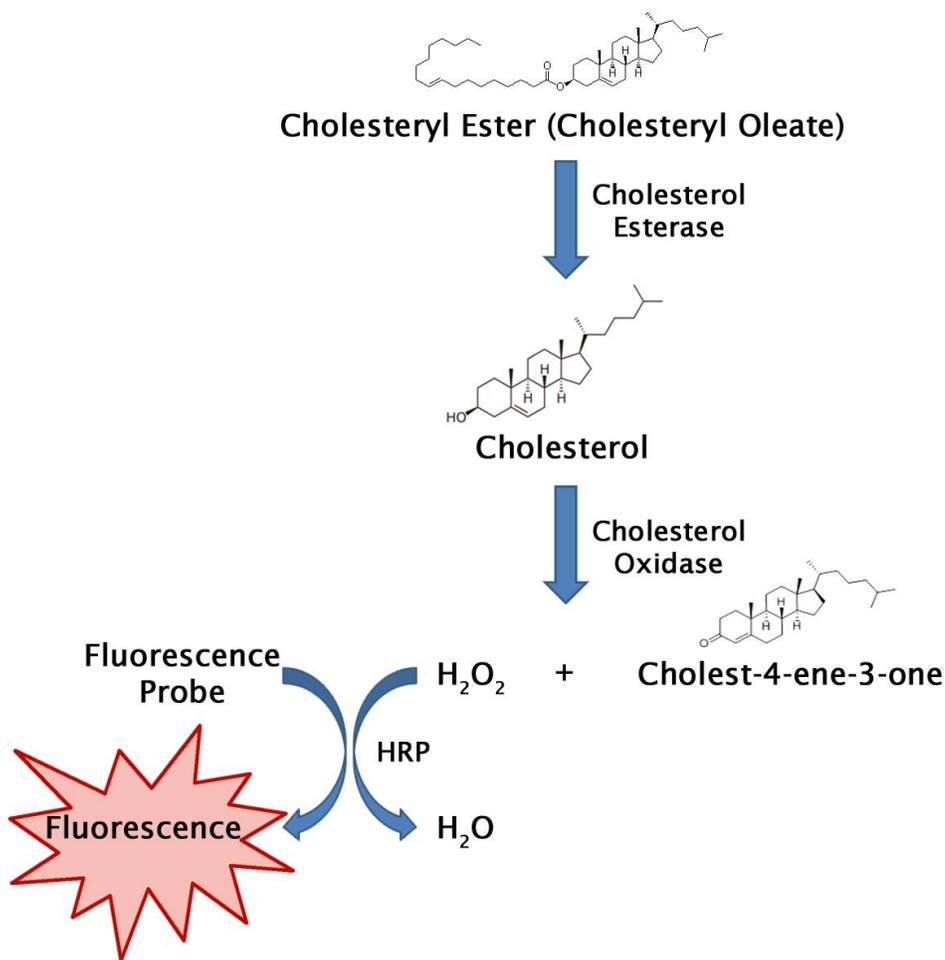
Cholesterol is a lipid sterol that is produced in and transported throughout the bloodstream in eukaryotes. Cholesterol is a critical compound used in the structure of cell membranes, hormones, and cell signaling. It is an essential component of animal cell structure in order to maintain permeability 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 fluorometric 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 fluorescence 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 within the 96-well microtiter plate format. Samples and standards are incubated for 45 minutes and then read with a standard 96-well fluorometric plate reader (Figure 1).



**Figure 1.** Cholesterol Assay Principle

**Related Products**

1. STA-241: Human Low Density Lipoprotein
2. STA-367: Human ApoE ELISA Kit
3. STA-369: OxiSelect™ Human Oxidized LDL ELISA Kit (MDA-LDL Quantitation)
4. STA-391: HDL and LDL/VLDL Cholesterol Assay Kit

## **Kit Components**

### **Box 1 (shipped at room temperature)**

1. 96-well Microtiter Plate (Part No. 234501): Two 96-well clear bottom black plates.
2. Cholesterol Standard (Part No. 239001): One 50  $\mu$ L tube of a 10 mM cholesterol solution in ethanol.
3. Assay Diluent (5X) (Part No. 239002): One 100 mL bottle.
4. Fluorescence Probe (Part No. 239005): One 200  $\mu$ L tube in DMSO.
5. HRP (Part No. 234402): Two 100  $\mu$ L 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  $\mu$ L tube.

## **Materials Not Supplied**

1. Distilled or deionized water
2. 1X PBS
3. Fluorescence microplate reader capable of reading excitation in the 530-570 nm range and emission in the 590-600 nm range.
4. Superoxide dismutase (optional)

## **Storage**

Upon receipt, store the Cholesterol Standard, Fluorescence Probe, HRP, Cholesterol Oxidase, and Cholesterol Esterase at  $-20^{\circ}\text{C}$ . The Fluorescence Probe is light sensitive and must be stored accordingly. Avoid multiple freeze/thaw cycles. Store the remaining kit components at  $4^{\circ}\text{C}$ .

## **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  $4^{\circ}\text{C}$  up to six months.
- Cholesterol Esterase: Reconstitute the powder with 200  $\mu$ L of 1X Assay Diluent. Vortex vigorously until dissolved. Prepare aliquots and store at  $-20^{\circ}\text{C}$  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, Fluorescence Probe 1:50, and Cholesterol Esterase 1:250 in 1X Assay Diluent. (eg. For 100 assays, combine 100  $\mu$ L of Cholesterol Oxidase, 100  $\mu$ L of HRP, 100  $\mu$ L Fluorescence Probe, and 20  $\mu$ 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 Fluorescence 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 µL 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 µL 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 µL of extracted sample per assay. Next, adjust the volume to 50 µL 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<sup>6</sup> cells, extract with 200 µL 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 µL 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 µL of extracted sample per assay. Next, adjust the volume to 50 µL 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 1:200 to 1:400 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 1:200 to 1:400 with Assay Diluent. This will provide values within the range of the standard curve.

Notes:

1. Samples with NADH concentrations above 10  $\mu\text{M}$  and glutathione concentrations above 50  $\mu\text{M}$  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  $\beta$ -mercaptoethanol since the fluorescence probe is not stable in the presence of thiols (above 10  $\mu\text{M}$ ).

### **Preparation of Cholesterol Standard Curve**

1. Prepare fresh cholesterol standards by diluting in 1X Assay Diluent. First, dilute the stock Cholesterol Standard 10 mM solution 1:50 in 1X Assay Diluent for a 200  $\mu\text{M}$  solution. (eg. add 20  $\mu\text{L}$  of the stock 10 mM standard to 980  $\mu\text{L}$  of 1X Assay Diluent). Vortex thoroughly.
2. Use this 200  $\mu\text{M}$  solution to prepare a series of the remaining cholesterol standards according to Table 1 below.

<b>Tubes</b>	<b>200 <math>\mu\text{M}</math> Cholesterol Standard (<math>\mu\text{L}</math>)</b>	<b>1X Assay Diluent (<math>\mu\text{L}</math>)</b>	<b>Resulting Cholesterol Concentration (<math>\mu\text{M}</math>)</b>
1	60	940	12
2	50	950	10
3	40	960	8
4	30	970	6
5	20	980	4
6	10	990	2
7	5	995	1
8	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  $\mu\text{L}$  of the diluted cholesterol standards or samples to the 96-well microtiter plate.
2. Add 50  $\mu\text{L}$  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 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 cholesterol within samples by comparing the sample RFU 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.

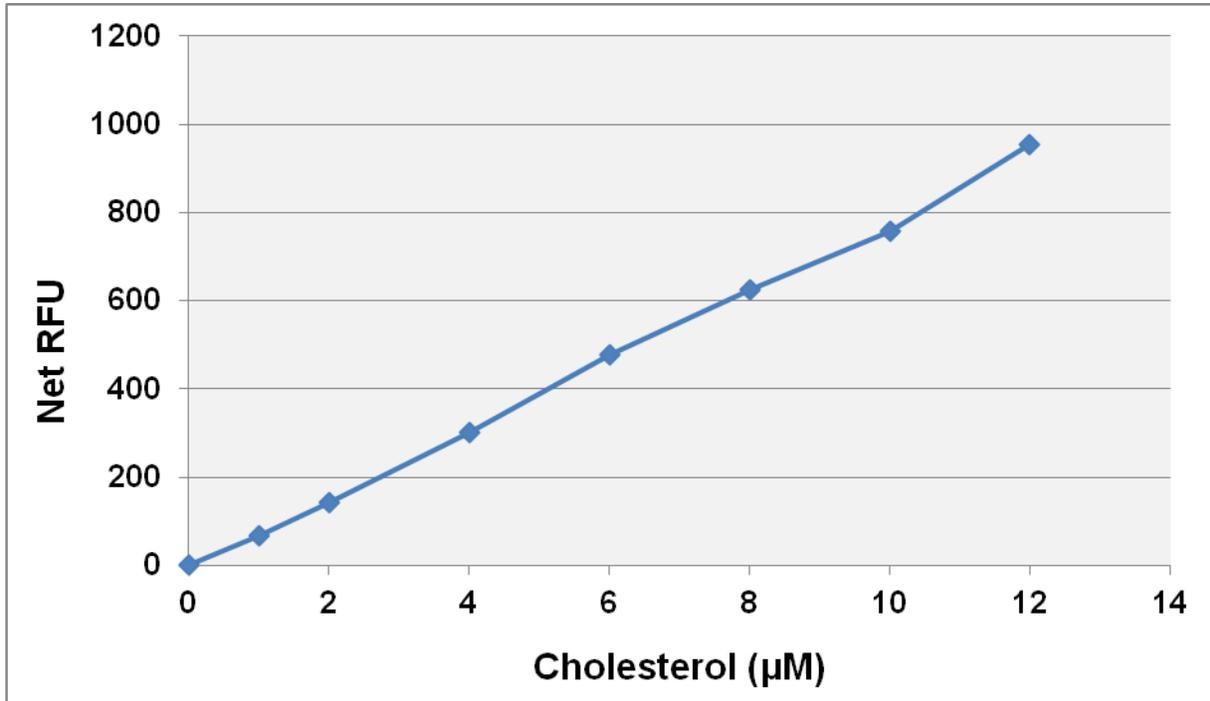


Figure 2: Cholesterol Standard Curve.

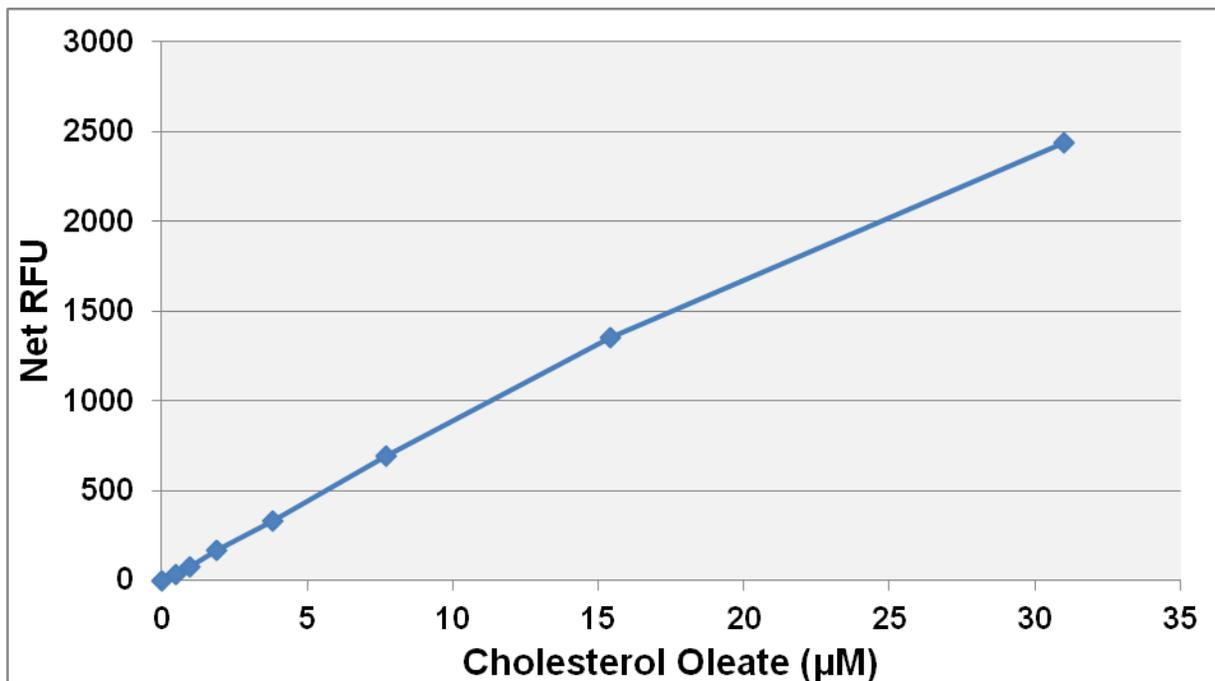


Figure 3: Cholesterol Oleate tested in Total Cholesterol Assay Kit.

## Calculation of Results

1. Calculate the average fluorescence 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 fluorescence.
2. Plot the corrected fluorescence 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 fluorescence values for each sample. Remember to account for dilution factors.

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

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

*Note: For the conversion of results from  $\mu\text{M}$  to  $\text{mg/dl}$ , divide the cholesterol concentration ( $\mu\text{M}$ ) 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. Tran, G.B. et al. (2023). Caffeine supplementation and FOXM1 inhibition enhance the antitumor effect of statins in neuroblastoma. *Cancer Res.* doi: 10.1158/0008-5472.CAN-22-3450.
2. Zahid, S. et al. (2023). The Geroprotective Drug Candidate CMS121 Alleviates Diabetes, Liver Inflammation, and Renal Damage in db/db Leptin Receptor Deficient Mice. *Int J Mol Sci.* **24**(7):6828. doi: 10.3390/ijms24076828.
3. Wang, Y. et al. (2021). N1-methyladenosine methylation in tRNA drives liver tumorigenesis by regulating cholesterol metabolism. *Nat Commun.* **12**(1):6314. doi: 10.1038/s41467-021-26718-6.
4. Deng, Q. et al. (2020). Dietary Lactic Acid Bacteria Modulate Yolk Components and Cholesterol Metabolism by Hmgr Pathway in Laying Hens. *Braz. J. Poult.* **22**(3): eRBCA-2020-1261. doi: 10.1590/1806-9061-2020-1261.
5. Singhal, A. et al. (2020). 2-Hydroxypropyl-gamma-cyclodextrin overcomes NPC1 deficiency by enhancing lysosome-ER association and autophagy. *Sci Rep.* **10**(1):8663. doi: 10.1038/s41598-020-65627-4.
6. Kidnapillai, S. et al. (2019). Drugs Used in the Treatment of Bipolar Disorder and their Effects on Cholesterol Biosynthesis- A Possible Therapeutic Mechanism. *World J Biol Psychiatry.* doi: 10.1080/15622975.2019.1669823.

7. Le Menn, G. et al. (2019). Decrease in  $\alpha\beta/\gamma\delta$  T-cell ratio is accompanied by a reduction in high-fat diet-induced weight gain, insulin resistance, and inflammation. *FASEB J.* **33**(2):2553-2562. doi: 10.1096/fj.201800696RR.
8. Martiskainen, H. et al. (2017). DHCR24 exerts neuroprotection upon inflammation-induced neuronal death. *J Neuroinflammation.* **14**(1):215. doi: 10.1186/s12974-017-0991-6.
9. Hsieh, H.Y. et al. (2017). Quantification of Endogenous Cholesterol in Human Serum on Paper Using Direct Analysis in Real Time Mass Spectrometry. *Anal Chem.* **89**(11):6146-6152.
10. Joseph, B. K. et al. (2015). Inhibition of AMP kinase by the protein phosphatase 2A heterotrimer, PP2A<sup>ppp2r2d</sup>. *J Biol Chem.* doi:10.1074/jbc.M114.626259.
11. Rodriguez-Jorquera, I. A. et al. (2015). Transcriptional and physiological response of fathead minnows (*Pimephales promelas*) exposed to urban waters entering into wildlife protected areas. *Environ Pollut.* **199**:155-165.
12. Liu, Z. H. et al. (2014). Abnormal lipid metabolism down-regulates adenosine triphosphate synthase  $\beta$ -subunit protein expression in corpus cavernosum smooth muscle in vitro and in vivo. *Andrologia.* **46**:487-494.
13. Ananth, S. et al. (2014). Regulation of the cholesterol efflux transporters ABCA1 and ABCG1 in retina in hemochromatosis and by the endogenous siderophore 2, 5-dihydroxybenzoic acid. *Biochim Biophys Acta.* **1842**:603-612.

## **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](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://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.