

## Copper (Cu<sup>++</sup>) Oxidized Human Low Density Lipoprotein (LDL)

---

|                                    |  |
|------------------------------------|--|
| <b>CATALOG NUMBER:</b>             | STA-214  |
| <b>STORAGE:</b>                    | 4°C; DO NOT FREEZE   |
| <b>QUANTITY AND CONCENTRATION:</b> | 100 µg at 1.0 mg/mL in PBS, pH 7.5, containing 0.5 mM EDTA and 0.02% NaN <sub>3</sub>  |
| <b>SHELF LIFE:</b>                 | 6 months from date of receipt under proper storage conditions  |
| <b>SOURCE:</b>                     | Isolated from human plasma tested negative for Hepatitis surface antigens and HCV, HIV-I and HIV-II antibodies; purified by ultracentrifugation and oxidized with Cu <sup>++</sup> |
| <b>PURITY:</b>                     | 98% by SDS-PAGE  |

### Recent Product Citations

1. Kaçar, C. et al. (2015). Amperometric magnetoimmunoassay for the determination of lipoprotein (a). *Microchim Acta*. doi: 10.1007/s00604-015-1472-1.
2. Haller, E. et al. (2014). Gold nanoparticle–antibody conjugates for specific extraction and subsequent analysis by liquid chromatography–tandem mass spectrometry of malondialdehyde-modified low density lipoprotein as biomarker for cardiovascular risk. *Anal Chim Acta*. **857**:53-63.

### 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.

*This product is for RESEARCH USE ONLY; not for use in diagnostic procedures.*

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

©2011-2017: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.