

Human High Density Lipoprotein (HDL)

CATALOG NUMBER:	STA-243
STORAGE:	-20°C; aliquot to avoid freeze/thaw cycles
QUANTITY AND CONCENTRATION:	100 µg at 1.0 mg/mL in Tris-HCl, pH 8.0, containing 1 mM EDTA, 0.02% NaN ₃ and 50% glycerol
SHELF LIFE:	1 year from date of receipt under proper storage conditions
SOURCE:	Following ultracentrifugation, HDL is isolated from human plasma tested negative for Hepatitis surface antigens and HCV, HIV-I and HIV-II antibodies

Recent Product Citation

Zhu, L. et al. (2018). CETP Inhibition Improves HDL Function but Leads to Fatty Liver and Insulin Resistance in CETP-Expressing Transgenic Mice on a High-Fat Diet. *Diabetes*. **67**(12):2494-2506. doi: 10.2337/db18-0474.

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
www.cellbiolabs.com

©2011-2019: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.