

Goat Anti-Human Apolipoprotein E Polyclonal Antibody

CATALOG NUMBER:	STA-138	STORAGE:	-20°C
QUANTITY AND CONCENTRATION:	100 µg of affinity purified antibody at 1 mg/mL in PBS, pH 7.2, containing 0.5 mM EDTA and 0.02% NaN ₃		
SHELF LIFE:	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
HOST SPECIES:	Goat		
IMMUNOGEN:	Purified human apolipoprotein E		
SPECIFICITY:	Human apolipoprotein E		
APPLICATION:	Immunoblot (1:200 to 1:80,000) ELISA (1:200 to 1:80,000)		

Example of Results

The following figures demonstrate typical results. One should use the data below for reference only. This data should not be used to interpret actual results.

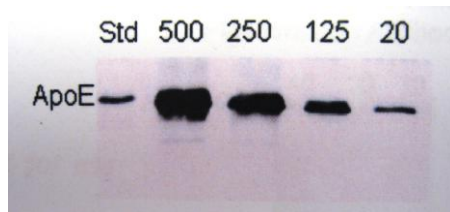


Figure 1. Immunoblot of Human Apolipoprotein E. Immunoblot using Goat Anti-Human Apolipoprotein E Polyclonal Antibody at 1:1000 dilution, followed by HRP-conjugated secondary antibody.

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.

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