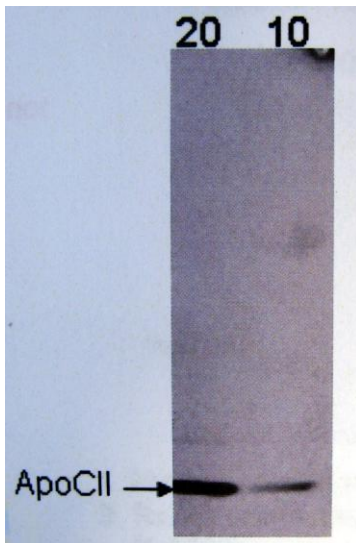


# Rabbit Anti-Human Apolipoprotein CII Polyclonal Antibody

<b>CATALOG NUMBER:</b>	STA-136	<b>STORAGE:</b>	-20°C
<b>QUANTITY AND CONCENTRATION:</b>	100 µg of affinity purified antibody at 1 mg/mL in PBS, pH 7.2, containing 0.5 mM EDTA and 0.02% NaN <sub>3</sub>		
<b>SHELF LIFE:</b>	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
<b>HOST SPECIES:</b>	Rabbit		
<b>IMMUNOGEN:</b>	Purified human apolipoprotein CII		
<b>SPECIFICITY:</b>	Human apolipoprotein CII		
<b>APPLICATION:</b>	Immunoblot (1:200 to 1:8,000) ELISA (1:200 to 1:8,000)		

## 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 CII.** Immunoblot using Rabbit Anti-Human Apolipoprotein CII 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'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: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.