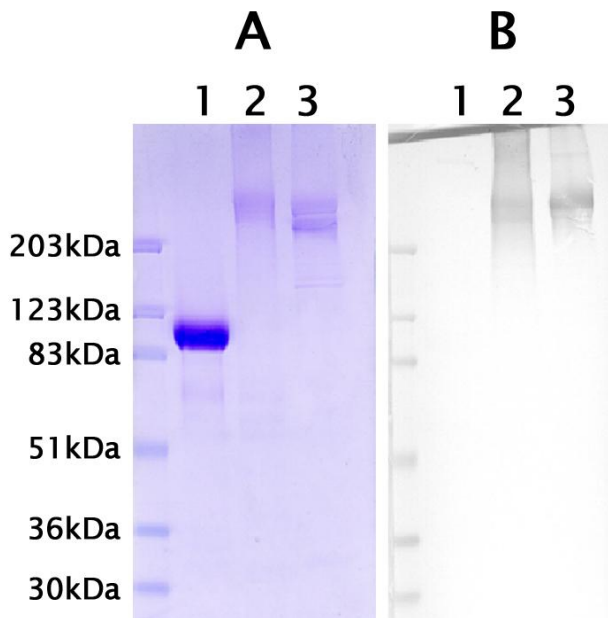


# Sheep Anti-Human Apolipoprotein (a) Polyclonal Antibody

<b>CATALOG NUMBER:</b>	STA-131	<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>	Sheep		
<b>IMMUNOGEN:</b>	Human apolipoprotein (a) purified from lipoprotein (a)		
<b>SPECIFICITY:</b>	Human lipoprotein (a) and apolipoprotein (a); minor reactivity with plasminogen; no reactivity with human apo-B-100/48		
<b>APPLICATION:</b>	Immunoblot (1:200 to 1:80,000) ELISA (1:200 to 1:80,000)		



**Figure 1. Immunoblot using Sheep Anti-Human Apolipoprotein(a) Polyclonal Antibody. A:** Human Plasminogen (1), Apolipoprotein(a) (2), and Lipoprotein(a) (3) were run on a 4-20% SDS Gel at 30 mA for 60 minutes. **B:** Western blot using primary antibody at 1:1000 dilution and HRP-labeled secondary antibody at 1:5000 dilution.

## **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)

©2011-2012: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.