

## Goat Anti-Nitrotyrosine Polyclonal Antibody

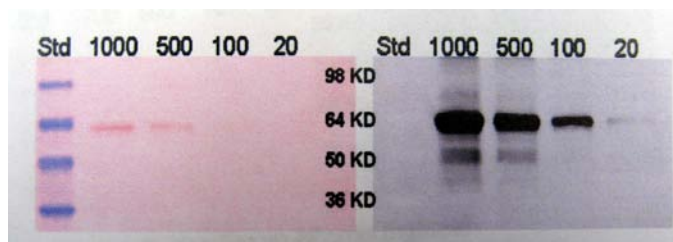
<b>CATALOG NUMBER:</b>	STA-003	<b>STORAGE:</b>	-20°C
<b>QUANTITY AND CONCENTRATION:</b>	100 µg of affinity purified antibody at 1 mg/mL in 75 mM PBS, pH 7.2, 0.5 mM EDTA and 0.02% NaN <sub>3</sub>		
<b>SHELF LIFE:</b>	1 year from date of receipt under proper storage conditions; aliquot to avoid multiple freeze thaw cycles		
<b>HOST SPECIES:</b>	Goat		
<b>IMMUNOGEN:</b>	Nitrated KLH		
<b>SPECIFICITY:</b>	Nitrated proteins		
<b>APPLICATION:</b>	Immunoblot (1:200 to 1:40,000) ELISA (1:200 to 1:40,000)		

### Background

The modification of tyrosine residues in proteins to 3-nitrotyrosine by peroxynitrite (Figure 1) or other potential nitrating agents has been detected in biological systems that are subject to oxidative stress. Detection of nitrotyrosine-containing proteins has been reported in many human and animal diseases or cellular models of disease. While all tyrosine residues in proteins may theoretically be targets for nitration, presumably the efficiency of tyrosine nitration is dependent on various biological conditions such as the local production and concentration of the reactive species, the existence and availability of antioxidants and scavengers, the accumulation of inflammatory cell and the presence of pro-inflammatory cytokines, as well as the proximity and compartmentation of these components.

### 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 Nitrated BSA. Left:** Ponceau S staining. **Right:** Immunoblot using Goat Anti-Nitrotyrosine Polyclonal Antibody at 1:1000 dilution, followed by HRP-conjugated secondary antibody. Numbers indicate ng/lane.

## **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.