

eNOS [pSer113] Antibody

MAN0004672

Rev. Date: 21 June 2011

Catalog Number: A14043

Store at -20°C

Clonality: Polyclonal
Volume: 100 µL

Host/Class: Rabbit IgG
Reactivity: Human eNOS
Predicted Reactivity: Human, Bovine

Product Description

Endothelial nitric-oxide synthase (eNOS) is an important enzyme in the cardiovascular system. It catalyzes the production of nitric oxide (NO), a key regulator of blood pressure, vascular remodeling and angiogenesis. The activity of eNOS is regulated by phosphorylation at multiple sites. eNOS is activated by VEGF, and this activation is associated with dephosphorylation of eNOS at serine 113. Cyclosporin A blocks this dephosphorylation of eNOS upon VEGF stimulation. The eNOS [pSer113] Antibody detects endogenous levels of eNOS only when phosphorylated at serine 113.

Product Specifications

Immunogen: Synthetic phosphopeptide (KLH-coupled) corresponding to residues surrounding Ser¹¹³ of human eNOS
Apparent MW: 140 kDa
Sequence Identity: Human
Lot: See product label

Product Applications

Application	Species	Test Material	Concentration
Western Blotting	Bovine	Aortic endothelial cells	1:1000

Storage and Handling

Store the antibody at -20°C. Do not make aliquots of the antibody.

Stability

Expires one year from date of receipt when stored as instructed.

Storage Buffer

10 mM sodium HEPES (pH 7.5), 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.

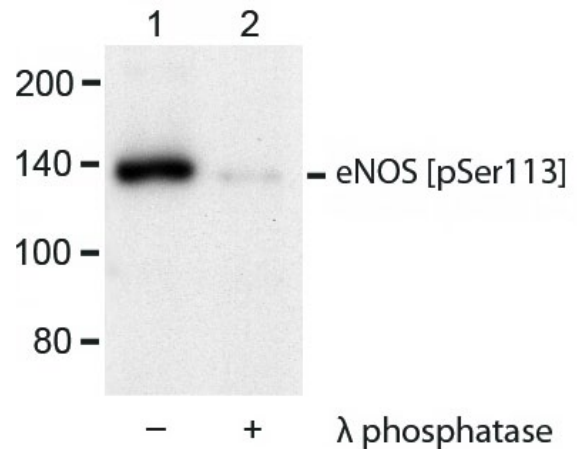


Figure 1 Western blot analysis of extracts from bovine aortic endothelial cells (BAEC), untreated (lane 1), or treated (lane 2) with λ phosphatase, using eNOS [pSer113] Antibody.

For research use only. Not for human or animal therapeutic or diagnostic use.

Manufacturing Site • 7335 Executive Way • Frederick • MD 21704 • E-mail: techsupport@invitrogen.com

Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at www.invitrogen.com/sds.









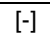
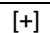


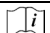
Certificate of Analysis

The Certificate of Analysis provides detailed quality control and product qualification information for each product. Certificates of Analysis are available on our website. Go to www.invitrogen.com/support and search for the Certificate of Analysis by product lot number, which is printed on the box.

Related Products

Product Name	Quantity	Catalog No.
iBlot® Dry Blotting System	1 unit	IB1001
WesternBreeze® Chromogenic Kit Anti-Rabbit	1 kit	WB7105
WesternBreeze® Chemiluminescent Kit, Anti-Rabbit	1 kit	WB7106
Goat anti-mouse (H+L), HRP conj.	1 mg	G21040
Goat anti-rabbit (H+L), HRP conj.	1 mg	G21234
Goat anti-mouse (H+L), AP conj.	1 mg	G21060
Goat anti-rabbit (H+L), AP conj.	1 mg	G21079
Nitrocellulose, 0.2 µm	20/pack	LC2000

Explanation of symbols

Symbol	Description	Symbol	Description
	Catalogue Number		Batch code
	Research Use Only		<i>In vitro</i> diagnostic medical device
	Use by		Temperature limitation
	Manufacturer		European Community authorised representative
	Without, does not contain		With, contains
	Protect from light		Consult accompanying documents
	Directs the user to consult instructions for use (IFU), accompanying the product.		

Limited Use Label License: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.