

A20/TNFAIP3 Antibody

MAN0004657

Rev. Date: 21 June 2011

Catalog Number: A14028

Store at -20°C

Clonality: Polyclonal
Volume: 100 µL

Host/Class: Rabbit IgG
Reactivity: Human A20/TNFAIP3
Predicted Reactivity: Human, Monkey

Product Description

A20, also referred to as TNF- α -induced protein 3 (TNFAIP3), is cytokine inducible protein which functions to inhibit apoptosis and activate NF- κ B. It is a 790-amino acid protein containing seven Cys2/Cys2-zinc finger motifs. Constitutive expression of A20 is observed in lymphoid tissues, but it is transiently expressed in a variety of cell types in response to inflammatory signals.

Product Specifications

Immunogen: Synthetic peptide corresponding to residues surrounding Gly267 of human A20/TNFAIP3
Alternate Names: TNFAIP3
Apparent MW: 82 kDa
Gene ID: 7128
Protein Accession No.: P21580
Sequence Identity: Human
Lot: See product label

Product Applications

Application	Species	Test Material	Concentration
Western Blotting	Human	THP-1 cells	1:1000

Additional applications for which this antibody has been used include: Immunoprecipitation (1:25 dilution).

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.

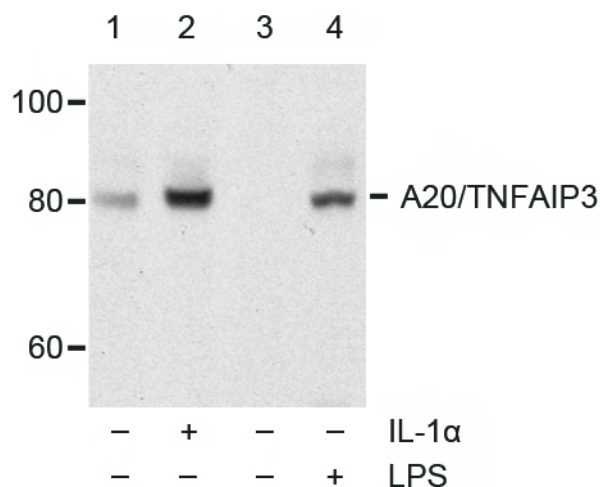


Figure 1 Western blot analysis of extracts from HeLa cells, untreated (lane 1), or treated (lane 2) with IL-1 α (20 ng/mL for 2 hours), and from differentiated THP-1 cells (treated with 80 nM TPA overnight), untreated (lane 3), or treated (lane 4) with LPS (1 µg/mL overnight), using A20/TNFAIP3 Antibody.

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.



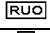
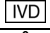



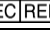
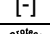
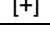



Certificate of Analysis

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		In vitro 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.