

Jagged1, Rabbit Monoclonal Antibody

Publication Part no. MAN0004945

Rev. Date: 21 July 2011

Catalog Number: A13984

Store at -20°C

Clonality: Monoclonal
Volume: 100 µL

Host/Class: Rabbit IgG
Reactivity: Human Jagged1
Predicted Reactivity: Human

Product Description

Notch signaling is activated upon engagement of the Notch receptor with its ligands, the DSL (Delta, Serrate, Lag2) proteins of the single-pass type I membrane protein family. Five DSL proteins have been identified in mammals: Jagged1, Jagged2, Delta-like (DLL) 1, 3 and 4. The proteins contain multiple EGF-like repeats and a DSL domain that is required for binding to Notch.

Product Specifications

Immunogen: Synthetic peptide corresponding to residues surrounding Glu¹¹⁴⁰ (intracellular region) of human Jagged1
Alternate Names: Jagged-1
Apparent MW: 180 kDa
Gene ID: 182
Protein Accession No.: P78504
Sequence Identity: Human
Clone/PAD: 28H8
Lot: See product label

Product Applications

Application	Species	Test Material	Concentration
Western blotting	Human	HepG2 cells	1:1000

Additional applications for which this antibody has been used include: Immunoprecipitation (1:50 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, 50% glycerol and less than 0.02% sodium azide.

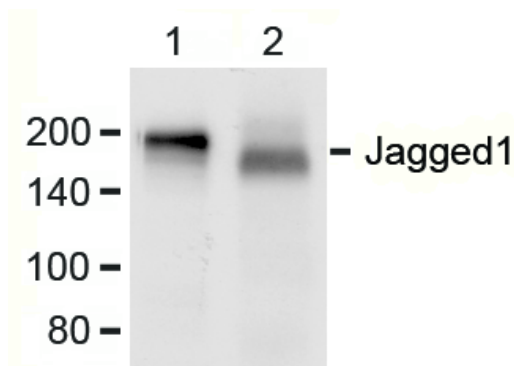


Figure 1 Western blot analysis of total cell lysates from HepG2 (lane1), and LNCaP (lane2) cells, using Jagged1 Rabbit Monoclonal Antibody.

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.









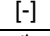
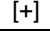


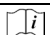
Certificate of Analysis

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.

For support visit www.invitrogen.com/support or email techsupport@invitrogen.com

www.lifetechnologies.com

 life
technologies™