

XIAP Monoclonal Antibody (clone 3B6), Rabbit

Store at -20°C

Catalog Number: A16522

Pub. No. MAN0009154 Rev. 1.00

Clonality: Monoclonal
Reactivity: Human XIAP

Amount: 100 µL
Species Reactivity: Human, Monkey

Host/Class: Rabbit IgG

Product Description

XIAP is a member of the inhibitor of apoptosis protein (IAP) family, consisting of an evolutionarily conserved group of apoptosis inhibitors containing a conserved 70 amino acid baculovirus inhibitor repeat (BIR) domain (1,2). Human members of this family include c-IAP1, c-IAP2, XIAP, survivin, livin, and NAIP. In general, IAP proteins function through direct interactions to inhibit the activity of caspases, including caspase-3, caspase-7, and caspase-9 (3,4). In addition, binding of IAP family members to the mitochondrial protein Smac blocks their interaction with caspase-9, thereby allowing the processing and activation of the caspase (2).

Product Specifications

Immunogen:	A synthetic peptide corresponding to a region surrounding Ser245 of human XIAP
Apparent MW:	53 kDa
Gene ID:	331
Protein Accession No.:	P98170
Clone/PAD:	3B6
Isotype:	IgG
Lot:	See product label

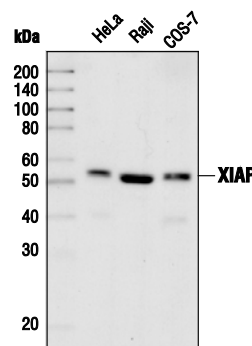


Figure 1 Western blot analysis of extracts from HeLa, Raji and COS cell lines, using XIAP Monoclonal Antibody.

Product Applications

Applications reported for this antibody include western blot. Because conditions may vary, it is recommended that each investigator determine the optimal amount of antibody to be used for each application.

Application	Recommended Dilution
Western blotting	1:1000

Storage and Handling

Store reagents at -20°C. Avoid repeated freezing and thawing. Do not aliquot the antibody.

Stability

When stored as instructed, expires one year from date of receipt unless otherwise indicated on product label.

Storage Buffer

10 mM sodium HEPES (pH 7.5), 150 mM NaCl, 100 µg/mL BSA, 50% glycerol, <0.02% sodium azide.

References

1. Deveraux, Q.L. and Reed, J.C. (1999) Genes Dev 13, 239-52.
2. Deveraux, Q.L. et al. (1998) EMBO J 17, 2215-23.
3. Kasof, G.M. and Gomes, B.C. (2001) J Biol Chem 276, 3238-46.
4. Deveraux, Q.L. et al. (1997) Nature 388, 300-4.

Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit www.lifetechnologies.com/support.

Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

Explanation of symbols

Symbol	Description	Symbol	Description	Symbol	Description
	Manufacturer		Catalog number		Batch code
	Use by		Temperature limitation		
	Consult instructions for use		Caution, consult accompanying documents		

Limited Use Label License No. 508: Cell Signaling Technology and Epitomics Technology for Antibody Production

This antibody is developed, validated, and produced by Cell Signaling Technology (CST) using in part technology under license (granting certain rights including those under U.S. Patents No. 5,675,063 and 7,429,487) from Epitomics, Inc.

DISCLAIMER: LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) DISCLAIM ALL WARRANTIES WITH RESPECT TO THIS DOCUMENT, EXPRESSED OR IMPLIED, INCLUDING BUT NOT LIMITED TO THOSE OF MERCHANTABILITY, FITNESS FOR A PARTICULAR PURPOSE, OR NON-INFRINGEMENT. TO THE EXTENT ALLOWED BY LAW, IN NO EVENT SHALL LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) BE LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY, OR UNDER ANY STATUTE OR ON ANY OTHER BASIS FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING BUT NOT LIMITED TO THE USE THEREOF.

©2013 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation and/or its affiliates or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com

www.lifetechnologies.com

28 August 2013

