

IκB-beta Polyclonal Antibody

Store at -20°C

Catalog Number: A16709

Pub. No. MAN0009199 Rev. 1.00

Clonality: Polyclonal

Amount: 100 μL

Host/Class: Rabbit

Reactivity: Human IκB beta

Species Reactivity: Human, Mouse, Rat, Monkey

Product Description

Cytosolic NF-κB/Rel transcription factors are in an inactive state, complexed with inhibitory IκB proteins (1-3). Activation occurs via phosphorylation of IκBα at Ser32 and Ser36 followed by proteasome-mediated degradation that results in the release and nuclear translocation of active NF-κB (3-7). IκBα phosphorylation and resulting Rel-dependent transcription are activated by a highly diverse group of extracellular signals including inflammatory cytokines, growth factors, and chemokines. Kinases that phosphorylate IκB at these activating sites have been identified (8).

Product Specifications

Immunogen:	A synthetic peptide corresponding to residues near the middle of human and mouse IκBβ
Apparent MW:	48 kDa
Gene ID:	4793
Protein Accession No.:	Q15653
Lot:	See product label

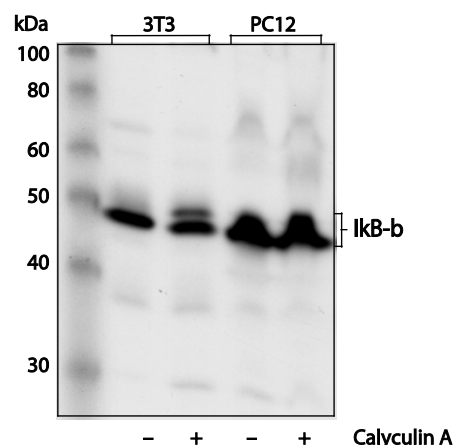


Figure 1 Western blot analysis of extracts from NIH-3T3 and PC12 cells, untreated or treated with calyculin A (50 nM for 15 minutes), using IκB-beta Polyclonal Antibody.

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.

References

1. Baeuerle, P.A. and Baltimore, D. (1988) Science 242, 540-6.
2. Beg, A.A. and Baldwin, A.S. (1993) Genes Dev 7, 2064-70.
3. Finco, T.S. et al. (1994) Proc Natl Acad Sci USA 91, 11884-8.
4. Brown, K. et al. (1995) Science 267, 1485-8.
5. Brockman, J.A. et al. (1995) Mol Cell Biol 15, 2809-18.
6. Traenckner, E.B. et al. (1995) EMBO J 14, 2876-83.
7. Chen, Z.J. et al. (1996) Cell 84, 853-62.
8. Karin, M. and Ben-Neriah, Y. (2000) Annu Rev Immunol 18, 621-63.








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		

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

