

## RelB Monoclonal Antibody (clone C1E4), Rabbit

Store at -20°C

Catalog Number: A16672

Pub. No. MAN0009141 Rev. 1.00

**Clonality:** Monoclonal  
**Reactivity:** Human RelB

**Amount:** 100 µL  
**Species Reactivity:** Human, Mouse, Rat, Monkey

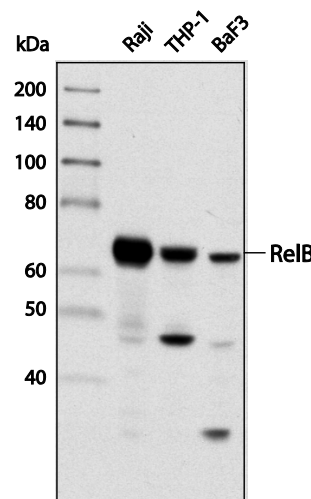
**Host/Class:** Rabbit IgG

### Product Description

Transcription factors of the nuclear factor  $\kappa$  B (NF- $\kappa$ B)/Rel family play a pivotal role in inflammatory and immune responses (1,2). There are five family members in mammals: RelA, c-Rel, RelB, NF- $\kappa$ B1 (p105/p50), and NF- $\kappa$ B2 (p100/p52). Both p105 and p100 are proteolytically processed by the proteasome to produce p50 and p52, respectively. Rel proteins bind p50 and p52 to form dimeric complexes that bind DNA and regulate transcription. In unstimulated cells, NF- $\kappa$ B is sequestered in the cytoplasm by I $\kappa$ B inhibitory proteins (3-5). NF- $\kappa$ B-activating agents can induce the phosphorylation of I $\kappa$ B proteins, targeting them for rapid degradation through the ubiquitin-proteasome pathway and releasing NF- $\kappa$ B to enter the nucleus where it regulates gene expression (6-8). NIK and IKK $\alpha$  (IKK1) regulate the phosphorylation and processing of NF- $\kappa$ B2 (p100) to produce p52, which translocates to the nucleus (9-11).

### Product Specifications

<b>Immunogen:</b>	A synthetic peptide corresponding to residues surrounding Ser424 of human RelB protein
<b>Apparent MW:</b>	70 kDa
<b>Gene ID:</b>	5971
<b>Protein Accession No.:</b>	Q01201
<b>Clone/PAD:</b>	C1E4
<b>Isotype:</b>	IgG
<b>Lot:</b>	See product label



**Figure 1** Western blot analysis of extracts from Raji, THP-1 and BaF3 cells using RelB Monoclonal Antibody.

### Product Applications

Applications reported for this antibody include western blot, and immunoprecipitation.

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
Immunoprecipitation	1:100

### 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.

For research use only. Not for use in diagnostic procedures

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

## References

1. Baeuerle, P.A. and Henkel, T. (1994) Annu Rev Immunol 12, 141-79.
2. Baeuerle, P.A. and Baltimore, D. (1996) Cell 87, 13-20.
3. Haskill, S. et al. (1991) Cell 65, 1281-9.
4. Thompson, J.E. et al. (1995) Cell 80, 573-82.
5. Whiteside, S.T. et al. (1997) EMBO J 16, 1413-26.
6. Traenckner, E.B. et al. (1995) EMBO J 14, 2876-83.
7. Scherer, D.C. et al. (1995) Proc Natl Acad Sci USA 92, 11259-63.
8. Chen, Z.J. et al. (1996) Cell 84, 853-62.
9. Senftleben, U. et al. (2001) Science 293, 1495-9.
10. Coope, H.J. et al. (2002) EMBO J 21, 5375-85.
11. Xiao, G. et al. (2001) Mol Cell 7, 401-9.






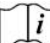

## Product Documentation

To obtain a Certificate of Analysis or Safety Data Sheets (SDSs), visit [www.lifetechnologies.com/support](http://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](http://www.lifetechnologies.com/termsandconditions). If you have any questions, please contact Life Technologies at [www.lifetechnologies.com/support](http://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](http://www.lifetechnologies.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[www.lifetechnologies.com](http://www.lifetechnologies.com)

28 August 2013

