

## NFκB-p65 (Lys310) acetylate Polyclonal Antibody

Store at -20°C

Catalog Number: A16567

Pub. No. MAN0009272 Rev. 1.00

Clonality: Polyclonal

Amount: 100 μL

Host/Class: Rabbit

Reactivity: NFκB-p65

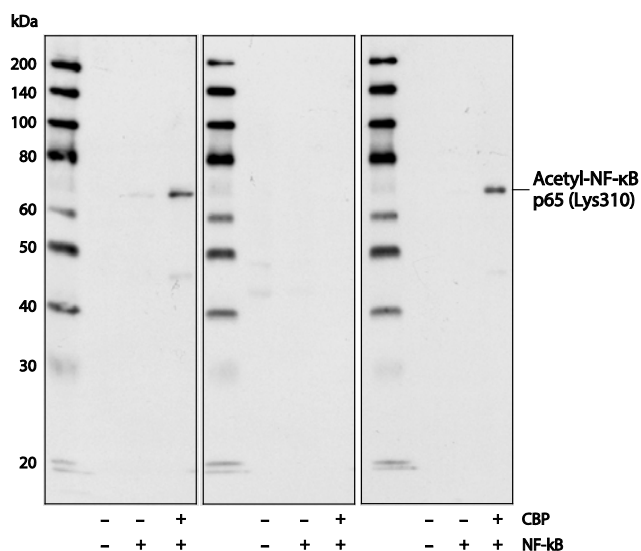
Species Reactivity: Human, Mouse

### Product Description

Transcription factors of the nuclear factor κ B (NF-κB)/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-κB1 (p105/p50), and NF-κB2 (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-κB is sequestered in the cytoplasm by IκB inhibitory proteins (3-5). NF-κB-activating agents can induce the phosphorylation of IκB proteins, targeting them for rapid degradation through the ubiquitin-proteasome pathway and releasing NF-κB to enter the nucleus where it regulates gene expression (6-8). NIK and IKKα (IKK1) regulate the phosphorylation and processing of NF-κB2 (p100) to produce p52, which translocates to the nucleus (9-11).

### Product Specifications

|                               |  |
|-------------------------------|--|
| <b>Immunogen:</b>             | A synthetic acetylated peptide corresponding to residues surrounding Lys310 of NF-κB |
| <b>Predicted Reactivity:</b>  | Rat, Monkey, Bovine, Dog   |
| <b>Apparent MW:</b>           | 65 kDa   |
| <b>Gene ID:</b>               | 5970   |
| <b>Protein Accession No.:</b> | Q04206   |
| <b>Lot:</b>                   | See product label  |



**Figure 1** Western blot analysis of extracts from HeLa cells, mock transfected or transfected with NF-κB or NF-κB plus CBP, using NFκB-p65 (Lys310) acetylate Polyclonal Antibody (left). Western blots were competed with acetylated peptide (center) or non-acetylated peptide (right).

### 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 | 1:50                 |

### 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.

For research use only. Not for use in diagnostic procedures

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

## Storage Buffer

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

## 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.




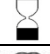
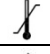
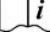

## Product Documentation

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 |   |             |

**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

*life*  
technologies™