

## BCL-6 Polyclonal Antibody

Store at -20°C

Catalog Number: A16633

Pub. No. MAN0009243 Rev. 1.00

Clonality: Polyclonal

Amount: 100 µL

Host/Class: Rabbit

Reactivity: Human Bcl-6

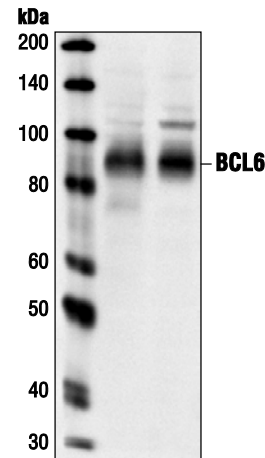
Species Reactivity: Human, Mouse

### Product Description

The BCL-6 is selectively expressed in mature B cells and encodes a nuclear phosphoprotein that belongs to the BTB/POZ zinc finger family of transcription factors (1,2). BCL-6 protein can bind to Stat6 target DNA sequences, and modulate the expression of interleukin-4-induced genes (3). Furthermore, BCL-6 restrains p53-dependent senescence (4). The mitogen-activated protein kinases, Erk1 and Erk2, but not JNK, phosphorylate BCL6 at multiple sites. Phosphorylation of BCL-6 at Ser333 and Ser343 results in degradation of BCL-6 by the ubiquitin/proteasome pathway in B cells (5,6). In addition, BCL-6 is acetylated and its transcriptional repressor function is inhibited by the transcriptional co-activator p300 (7).

### Product Specifications

<b>Immunogen:</b>	A synthetic peptide corresponding to residues near the carboxy-terminal sequence of human BCL6
<b>Predicted Reactivity:</b>	Rat
<b>Apparent MW:</b>	87-98 kDa
<b>Gene ID:</b>	604
<b>Protein Accession No.:</b>	P41182
<b>Lot:</b>	See product label



**Figure 1** Western blot analysis of Ramos and A20 cell extracts using BCL-6 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.

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. Onizuka, T. et al. (1995) Blood 86, 28-37.
2. Zollman, S. et al. (1994) Proc. Natl. Acad. Sci. USA 91, 10717-10721.
3. Harris, M. B. et al. (1999) Mol. Cell. Biol. 19, 7264-7275.
4. Shvarts, A. et al. (2002) Genes Dev. 16, 681-686.
5. Moriyama, M. et al. (1997) Oncogene 14, 2465-2474.
6. Niu, H. et al. (1998) Genes Dev. 12, 1953-1961.
7. Bereshchenko, O. R. et al. (2002) Nat. Genet. 32, 606-613.








## 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

