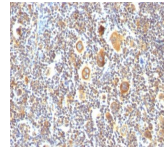


## Bcl-6 Monoclonal Mouse Antibody (BCL6/850)



### Product Description

Bcl-6 is a transcriptional repressor mainly required for germinal center (GC) formation and antibody affinity maturation which has different mechanisms of action specific to the lineage and biological functions. It forms complexes with different corepressors and histone deacetylases to repress the transcriptional expression of different subsets of target genes. Bcl-6 allows GC B-cells to proliferate very rapidly in response to T-cell dependent antigens and tolerate the physiological DNA breaks required for immunoglobulin class switch recombination and somatic hypermutation without inducing a p53/TP53-dependent apoptotic response. Antibodies to Bcl-6 are helpful in a number of diagnostic settings, for example in the differential diagnosis of small B-cell lymphoma, and in distinguishing classical Hodgkin lymphoma from nodular lymphocyte predominant Hodgkin lymphoma (NLPHL). Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors.

**Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email [order@biotium.com](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order. **Catalog number key for antibody number 0850,**

**Anti-Bcl-6 (BCL6/850)**

### References

Garcia JF, et al. 2006. J. Histochem Cytochem. 54:31. | Pasqualucci, L., et al. 2003. Leuk. Lymphoma 44 Suppl 3: S5-12. | Ree, H.J., et al. 2003. Hum. Pathol. 34: 610-616

### Product attributes

Antibody number	#0850
Antibody reactivity (target)	Bcl-6
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BCL6/850
Isotype	IgG1, kappa
Molecular weight	95 kDa
Human gene symbol	BCL6
Synonyms	Bcl-6; B-Cell Lymphoma 6 Protein; B-cell lymphoma 5 protein; Zinc Finger Protein 51; zinc finger transcription factor BCL6S; BCL6 Transcription Repressor
Entrez gene ID	604
SwissProt	P41182
Unigene	478588
Immunogen	Recombinant human bcl-6 protein
Antibody target cellular localization	Nucleus
Species reactivity	Human, Mouse
Antibody application notes	For coating for ELISA, order Ab without BSA. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Optimal dilution and staining procedure for a specific application should be determined by user. Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Positive control	Raji or Ramos cells. Tonsil or Hodgkin's lymphoma.
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Apoptosis, Cancer, Immunology, Transcription factor
Tumor expression	Leukemia/lymphoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF®405S Features</a>
BNC88	CF®488A	490/515	488	GFP, FITC	<a href="#">CF®488A Features</a>
BNC68	CF®568	562/583	532, 561	RFP, TRITC	<a href="#">CF®568 Features</a>
BNC94	CF®594	593/614	561	Texas Red®	<a href="#">CF®594 Features</a>
BNC40	CF®640R	642/662	633-640	Cy®5	<a href="#">CF®640R Features</a>
BNC47	CF®647	650/665	633-640	Cy®5	<a href="#">CF®647 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, LI-COR Bioscience.