Bcl-2 Monoclonal Mouse Antibody (8C8)



Product Description

This antibody recognizes a protein of 25-26 kDa, identified as Bcl-2. This antibody shows no cross-reaction with Bcl-x or Bax proteins. Bcl-2 suppresses apoptosis in a variety of cell systems including factor-dependent lymphohematopoietic and neural cells. It regulates cell death by controlling the mitochondrial membrane permeability. Bcl-2 appears to function in a feedback loop system with caspases. It inhibits caspase activity either by preventing the release of cytochrome c from the mitochondria and/or by binding to the apoptosis-activating factor (APAF-1). Expression of Bcl-2 inhibits programmed cell death (apoptosis). In most follicular lymphomas, neoplastic germinal centers express high levels of Bcl-2 protein, whereas the normal or hyperplastic germinal centers are negative. Consequently, this antibody is valuable when distinguishing between reactive and neoplastic follicular proliferation in lymph node biopsies. It may also be used in distinguishing between those follicular lymphomas that express Bcl-2 protein and the small number in which the neoplastic cells are Bcl-2 negative.

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 to inquire about stock status and lead times before placing your order.

Catalog number key for antibody number 0005, Anti-Bcl-2 (8C8)

Product attributes

Product attributes						
Antibody number	#0005					
Antibody reactivity (target)	Bcl-2					
Antibody type	Primary					
Host species	Mouse					
Clonality	Monoclonal					
Clone	8C8					
Isotype	lgG1, kappa					
Molecular weight	Mouse Monoclonal 8C8 IgG1, kappa 25-26 kDa Apoptosis regulator Bcl-2; B-cell CLL/lymphoma-2 BCL2 596 P10415 150749 A synthetic peptide, aa41-54 (GAAPAPGIFSSQPG-Cys) of human Bcl-2 protein. IHC (FFPE) (verified), WB (verified) Endoplasmic reticulum, Mitochondria, Nuclear membrane Human, Monkey, Pig Higher concentration may be required for direct detection with secondary antibody conjugates than for indirect detection with secondary antibody conjugates than for indirect detection with react with mouse or rat, others not tested, immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes, Flow Cytometry 0.5-1 ug/mlilion cells/c1 mL, Western blotting 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas. Room temperature					
Synonyms	Apoptosis regulator Bcl-2; B-cell CLL/lymphoma-2					
Human gene symbol	Primary Mouse Monoclonal 3C8 gG1, kappa 25-26 kDa Apoptosis regulator Bcl-2; B-cell CLL/lymphoma-2 3CL2 3CL2 596 P10415 150749 A synthetic peptide, aa41-54 (GAAPAPGIFSSQPG-Cys) of numan Bcl-2 protein. HC (FFPE) (verified), WB (verified) Endoplasmic reticulum, Mitochondria, Nuclear membrane Human, Monkey, Pig Higher concentration may be required for direct detection with secondary antibody conjugates than for indirect detection with secondary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not eact with mouse or rat, others not tested, Immunohistology formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, 3H 9.0, for 10-20 min followed by cooling at RT for 20 minutes, 15-10 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular ymphomas. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Vote: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP			BCL2		
Entrez gene ID	596					
SwissProt	P10415					
Unigene	150749					
Immunogen	A synthetic peptide, aa41-54 (GAAPAPGIFSSQPG-Cys) of human Bcl-2 protein.					
Verified antibody applications	IHC (FFPE) (verified), WB (verified)					
Antibody target cellular localization	Endoplasmic reticulum, Mitochondria, Nuclear membrane					
Species reactivity	Human, Monkey, Pig					
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Does not react with mouse or rat, others not tested, Immunohistology formalin-fixed 0.5-1 ug/mL, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes.					
	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user					
Positive control	0.5-1 ug/mL, Optimal dilution for a specific application should be					
Positive control Shipping condition	0.5-1 ug/mL, Óptimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular					
	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas.					
Shipping condition	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,					
Shipping condition Storage Conditions	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when					
Shipping condition Storage Conditions Shelf life	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended					
Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended For research use only (RUO) 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, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 0.2 mg/mL in PBS/0.05% azide, Purified: 0.2					

Call us: 800-304-5357 Email: btinfo@biotium.com

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 2, 2025 at 02:49:28 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-bcl-2-8c8/