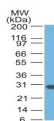


Bcl-x Monoclonal Mouse Antibody (BX006)



Product Description

Bcl-x belongs to the Bcl-2 family of proteins and plays a crucial role in apoptosis. Bcl-X has two isoforms: Bcl-XL (long), a 26 kDa protein which suppresses cell death, and Bcl-XS (short), a 178 amino acid protein lacking a 63 amino acid domain which functions as a dominant inhibitor of Bcl-2 and promotes apoptosis. This MAb reacts with both Bcl-XS and Bcl-XL proteins, and shows no cross-reaction with Bcl-2 or Bax proteins.

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 0006, Anti-Bcl-x (BX006)

Product attributes

Antibody number	#0006
Antibody reactivity (target)	Bcl-x
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BX006
Isotype	IgG2a
Molecular weight of antigen	27 kDa
Synonyms	Bcl-x; Bcl-2-like protein 1; BCL2L1; Apoptosis regulator Bcl-X
Human gene symbol	BCL2L1
Entrez gene ID	598
SwissProt	Q07817
Unigene	516966
Immunogen	A synthetic peptide, aa 3-14 (Cys-QSNRELVDVFLS) of human Bcl-X protein.
Species reactivity	Human, Mouse, Pig, Rat
Verified antibody applications	IHC (FFPE) (verified), WB (verified)
Antibody target cellular localization	Cytoplasmic, Endoplasmic reticulum, Mitochondria
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. 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/million cells/0.1 mL. Western blotting 0.5-1 ug/mL. Optimal dilution for a specific application should be determined by user
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)
Antibody research areas	Apoptosis, Cancer
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
Positive control	Jurkat, K562, HL-60, or HeLa Cells. Reed-Sternberg cells in Hodgkin's lymphoma.
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

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 June 3, 2026 at 08:52:28 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-bcl-x-bx006/>