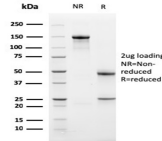


Bcl-x Monoclonal Mouse Antibody (BCL2L1/1762)



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 1762, Bcl-x Monoclonal Mouse Antibody (BCL2L1/1762)**

Product attributes

Antibody number	#1762
Antibody reactivity (target)	Bcl-x
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BCL2L1/1762
Isotype	IgG2b, kappa
Molecular weight	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	Recombinant human BCL2L1 protein fragment (around aa 1-135) (exact sequence is proprietary)
Antibody target cellular localization	Cytoplasmic, Endoplasmic reticulum, Mitochondria
Species reactivity	Human
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
Positive control	Jurkat, K562, HL-60, or HeLa Cells. Reed-Sternberg cells in Hodgkin's lymphoma.
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
Regulatory status	For research use only (RUO)
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody research areas	Apoptosis, Cancer

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel
BNC04	CF®405S	404/431	405	DAPI (microscopy), AF405
BNC88	CF®488A	490/515	488	GFP, FITC
BNC68	CF®568	562/583	532, 561	RFP, TRITC
BNC94	CF®594	593/614	561	Texas Red®
BNC40	CF®640R	642/662	633-640	Cy®5
BNC47	CF®647	650/665	633-640	Cy®5
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

Dye Features

[CF®405S Features](#)

[CF®488A Features](#)

[CF®568 Features](#)

[CF®594 Features](#)

[CF®640R Features](#)

[CF®647 Features](#)

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.

This datasheet was generated on June 30, 2024 at 06:16:21 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/bcl-x-mono-clonal-mouse-antibody-bcl2l1-1762/>