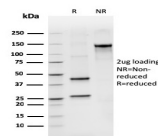


# Bcl-2-like Protein 2 Monoclonal Mouse Antibody (CPTC-BCL2L2-2)



## Product Description

Bcl-2-like protein 2 (BCL2L2) is a member of the Bcl-2 protein family. Bcl-2-like protein 2 promotes cell survival and blocks dexamethasone-induced apoptosis. It also mediates survival of postmitotic Sertoli cells by suppressing death-promoting activity of BAX.

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 2265, BCL2-like 2 Monoclonal Mouse Antibody (CPTC-BCL2L2-2)**

## Product attributes

Antibody number	#2265
Antibody reactivity (target)	Bcl-2-like protein 2
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	CPTC-BCL2L2-2
Isotype	IgG1, kappa
Molecular weight	22-30 kDa
Synonyms	Bcl-2-like Protein 2; BCL2L2; BCL-W; BCL2-L-2; BCL-W
Human gene symbol	BCL2L2
Entrez gene ID	599
SwissProt	Q92843
Unigene	410026
Immunogen	Recombinant human full-length BCL2L2 protein
Verified antibody applications	WB (verified)
Antibody target cellular localization	Mitochondria
Species reactivity	Human
Antibody application notes	Western blot 1-2 ug/mL; Optimal dilution for a specific application should be determined.
Positive control	PC-3, HEK293, HePG2, MCF-7, U87, Human colon extract; human spleen.
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
Validated in protein array	Monospecific
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Antibody research areas	Apoptosis, Cancer
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	<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>
BNC74	CF®740	742/767	633-685	775/50	<a href="#">CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 4, 2025 at 12:31:34 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/bcl2-like-2-monoclonal-mouse-antibody-cptc-bcl2l2-2/>