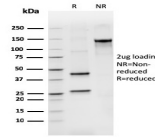


# Bcl-2-like Protein 2 Monoclonal Mouse Antibody (CPTC-BCL2L2-2)



## 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-L2; 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

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.