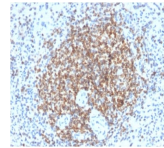


Bcl-2 Recombinant Monoclonal Rabbit Antibody (BCL2/2210R)



This antibody recognizes a protein of 25-26 kDa, identified as the bcl-2 α oncoprotein.

Product Description

This antibody recognizes a protein of 25-26 kDa, identified as the bcl-2 oncoprotein. It shows no cross-reaction with Bcl-x or Bax protein. Expression of bcl-2 oncoprotein inhibits the 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 $\text{\textcircled{R}}$ dyes and other labels. CF $\text{\textcircled{R}}$ dyes offer exceptional brightness and photostability. See the [CF \$\text{\textcircled{R}}\$ Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF $\text{\textcircled{R}}$ 405S and CF $\text{\textcircled{R}}$ 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 $\text{\textcircled{R}}$ 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 2210, Anti-Bcl-2 (BCL2/2210R)**

Antibody #	prefix	Conjugation	Ex/Em	Concentration	Storage Buffer
BNC04	CF $\text{\textcircled{R}}$ 405S		404/431 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC05	CF $\text{\textcircled{R}}$ 405M		408/452 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC06	CF $\text{\textcircled{R}}$ 405L		395/545 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC88	CF $\text{\textcircled{R}}$ 488A		490/515 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC14	CF $\text{\textcircled{R}}$ 514		516/548 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC43	CF $\text{\textcircled{R}}$ 543		541/560 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC55	CF $\text{\textcircled{R}}$ 555		555/565 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC68	CF $\text{\textcircled{R}}$ 568		562/583 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC94	CF $\text{\textcircled{R}}$ 594		593/614 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC40	CF $\text{\textcircled{R}}$ 640R		642/662 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC47	CF $\text{\textcircled{R}}$ 647		650/665 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC60	CF $\text{\textcircled{R}}$ 660C		667/685 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC61	CF $\text{\textcircled{R}}$ 660R		663/682 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC80	CF $\text{\textcircled{R}}$ 680		681/698 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC81	CF $\text{\textcircled{R}}$ 680R		680/701 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC00	CF $\text{\textcircled{R}}$ 700		695/720 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNC70	CF $\text{\textcircled{R}}$ 770		770/797 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCR	R-PE (PE)		496, 546, 565/578 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCA	APC		650/660 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCP	PerCP		482/677 nm	0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCB	Biotin	N/A		0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCAP	Alkaline Phosphatase	N/A		0.1 mg/mL	PBS, 0.1% BSA, 0.05% azide
BNCH	Horse radish Peroxidase	N/A		0.1 mg/mL	PBS, 0.05% BSA, no azide
BNUB	Purified, with BSA	N/A		0.2 mg/mL	PBS, 0.05% BSA, 0.05% azide
BNUM	Purified, BSA-free	N/A		1 mg/mL	PBS, no BSA, no azide

References

Adams, J.M., et al. 1998. The Bcl-2 protein family: arbiters of cell survival. *Science* 281: 1322-1326.

Product attributes

Antibody number	2210
Reactivity (target)	Bcl-2
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	BCL2/2210R
Isotype	IgG
Molecular weight	25-26 kDa
Synonyms	Apoptosis regulator Bcl-2, B-cell CLL/lymphoma-2
Human gene symbol	BCL2
Entrez gene ID	596
SwissProt	P10415
Unigene	150749
Immunogen	Recombinant full-length human bcl-2 protein
Cellular localization	Endoplasmic reticulum, Mitochondria, Nuclear membrane
Species reactivity	Human
Applications	Immunohistology (formalin), Western
Application notes	Immunohistology (formalin): 0.5-1 μ g/mL for 30 minutes at RT. Staining of formalin-fixed tissues is enhanced by boiling tissue sections in 10 mM Tris with 1 mM EDTA Buffer pH 9.0 for 10-20 minutes followed by cooling at RT for 20 minutes. Western blotting 0.5-1 μ g/mL. Optimal dilution for a specific application should be determined by user
Positive control	Jurkat, K562, HL-60, or HeLa Cells. Tonsil or follicular lymphomas.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 $^{\circ}$ C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 $^{\circ}$ C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Supplied As	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, BSA-free: 1 mg/mL in PBS without azide, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide
Antibody research areas	Apoptosis, Cancer