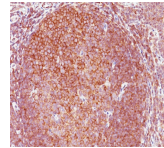


Bcl10 Monoclonal Mouse Antibody (151)



Product Description

B-cell lymphoma/leukemia 10 (BCL10) is a signaling protein that promotes apoptosis, pro-caspase-9 maturation, and activation of NF-kappa-B via NIK and IKK. BCL10 contains an N-terminal caspase recruitment domain (CARD), which is found in a number of apoptotic regulatory molecules. BCL10 is involved in the adaptive immune response, and may be an adapter protein between upstream TNFR1-TRADD-RIP complex and the downstream NIK-IKK-IKAP complex. A BCL10 translocation is recurrent in low-grade mucosa-associated lymphoid tissue (MALT lymphoma). 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 1852, BCL-10 Monoclonal Mouse Antibody (151)**

Product attributes

Antibody number	#1852
Antibody reactivity (target)	BCL10
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	151
Isotype	IgG1, kappa
Molecular weight	33 kDa
Synonyms	BCL10; BCL-10; B-cell CLL/lymphoma 10; B-cell leukemia/lymphoma 10
Human gene symbol	BCL10
Entrez gene ID	8915
SwissProt	O95999
Unigene	193516
Immunogen	Human BCL10 recombinant protein (epitope aa122-168)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoplasmic, Nucleus
Species reactivity	Human
Positive control	HepG2 cells. 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, Immunology, Signal transduction

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
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, LI-COR Bioscience.