

CD5 Monoclonal Mouse Antibody (B-B8)

Product Description

This antibody recognizes a 67 kDa transmembrane protein which is identified as CD5. The CD5 antigen is found on 95% of thymocytes and 72% of peripheral blood lymphocytes. In lymph nodes, the main reactivity is observed in T cell areas. Anti-CD5 is a pan T-cell marker that also reacts with a range of neoplastic B-cells, e.g. chronic lymphocytic leukemia/small lymphocytic lymphoma (CLL/SLL), mantle cell lymphoma, and a subset (~10%) of diffuse large B-cell lymphoma. CD5 aberrant expression is useful in making a diagnosis of mature T-cell neoplasms. Note that this MAb is not suitable for frozen tissues. Catalog number key for antibody number 0995, Anti-CD5 (B-B8)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#0995		
Antibody reactivity (target)	CD5		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	B-B8		
Isotype	IgG1, kappa		
Molecular weight	67 kDa		
Synonyms	CD5 antigen (p56 62), LEU1, Ly12, LyA, Lymphocyte antigen T1/Leu-1, Lymphocyte glycoprotein T1/Leu1, T-cell surface glycoprotein CD5		
Human gene symbol	CD5		
Entrez gene ID	921		
SwissProt	P06127		
Unigene	58685		
Immunogen	T cells activated with PHA		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Expected antibody applications	Flow (surface) (expected), Functional studies (published for clone), IF (expected)		
Antibody application notes	Not suitable for frozen tissues, Flow cytometry: 0.5-1 ug/million cells, Immunofluorescence: 0.5-1 ug/mL, Optimal dilution for a specific application should be determined by user		
Positive control	293T, Ramos, or MOLT-4 Cells. Tonsil.		
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/m in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL PBS without azide		
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended		
Cell/tissue expression	Lymphocytes, T-cells		
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.		
Tumor expression	Leukemia/lymphoma		
Antibody research areas	Immunology		

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
BNC74	CF®740	742/767	633-685	775/50	CF®740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. Inf Immunity (1996) 64(5): 1595-1599. (cell depletion)
- 2. Br J Haematol (2013) 163: 182-193. (antibody internalization)

This datasheet was generated on November 30, 2025 at 02:08:50 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd5-monoclonal-mouse-antibody-b-b8/