CD20 Polyclonal Rabbit Antibody



Product Description

Recognizes a protein of 30-33 kDa, which is identified as CD20. It is a non-Ig differentiation antigen of B-cells and its expression is restricted to normal and neoplastic B-cells, being absent from all other leukocytes and tissues. CD20 is expressed by pre B-cells and persists during all stages of B-cell maturation but is lost upon terminal differentiation into plasma cells. This antibody can be used for immunophenotyping of leukemia and malignant cells, B lymphocyte detection in peripheral blood and B cell localization in tissues. It reacts with the majority of B-cells present in peripheral blood and lymphoid tissues and their derived lymphomas. In lymphoid tissue, germinal center blasts and B-immunoblasts are particularly reactive. It is a reliable antibody for ascribing a B-cell phenotype in known lymphoid tissues. Rarely, CD20-positive T-cell lymphomas have been reported. Reactivity has also been noted with Reed-Sternberg cells in cases of Hodgkin's disease, particularly of lymphocyte predominant type.

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 1498, Anti-CD20|MS4A1

Product attributes

Tumor expression

Call us: 800-304-5357

Product attributes			
Antibody number	#1498		
Antibody reactivity (target)	y reactivity (target) CD20, MS4A1		
Antibody type	rimary		
Host species	Rabbit		
Clonality	Polyclonal		
Isotype	IgG		
Molecular weight	33-37 kDa		
Synonyms	APY; ATOPY; B-lymphocyte antigen CD20; B-lymphocyte cell-surface antigen B1; Bp35; Fc epsilon receptor I beta chain; Fc Fragment of IgE high affinity I receptor for beta polypeptide; FCER1B; High affinity immunoglobulin epsilon receptor subunit beta; IgE Fc receptor subunit beta; IGEF; IGHER; LEU16 Leukocyte surface antigen Leu-16; Ly44; Membrane spanning domains subfamily A member 2; Membrane-spanning 4-domain subfamily A member 1 (MS4A1)		
Human gene symbol	MS4A1		
Entrez gene ID	931		
SwissProt	P11836		
Unigene	712553		
Immunogen	combinant full-length human MS4A1 protein		
Verified antibody applications	IHC (FFPE) (verified)		
Antibody target cellular localization	Plasma membrane		
Species reactivity	Human		
Species reactivity Positive control	Human Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils.		
	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes		
Positive control	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils.		
Positive control Shipping condition	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Positive control Shipping condition Storage Conditions	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) Conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified: 8SA-free: 1 mg/mL in		
Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	Daudi, Raji, and U266, and human lymphocytes. Lymph nodes and tonsils. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C For research use only (RUO) 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 Guaranteed for at least 24 months from date of receipt when		

Leukemia/lymphoma

Email: btinfo@biotium.com

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.

This datasheet was generated on November 2, 2025 at 03:51:39 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd20-polyclonal-rabbit-antibody/