CD20 / MS4A1 Monoclonal Mouse Antibody (L26)

Product Description

This antibody recognizes a protein of 30-33 kDa, which is identified as CD20. Its epitope is located in the cytoplasmic domain of CD20 and was, therefore, ascribed as CD20cy in the 5th Workshop. CD20 is a non-lg 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 MAb 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. Catalog number key for antibody number 0766, Anti-CD20|MS4A1 (L26)

Call us: 800-304-5357 Email: techsupport@biotium.com

Product attributes

Product attributes				
Antibody number	#0766			
Antibody reactivity (target)	CD20, MS4A1			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	L26			
Isotype	IgG2a, kappa			
Molecular weight	33-37 kDa			
Synonyms	APY; ATOPY; 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; GER; IGER; IGHER; Leukocyte surface antigen Leu-16; Ly44; MS4A1; MS4A2			
Human gene symbol	MS4A1			
Entrez gene ID	931			
SwissProt	P11836			
Unigene	712553			
Immunogen	Human tonsil B cells			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	Flow (verified), IF (verified), IHC (FFPE) (verified), WB (verified)			
Species reactivity	Human			
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistochemistry (formalin-fixed): 0.5-1 ug/mL for 30 minutes at RT, Western blot: 1-2 ug/mL, Flow cytometry: 0.5-1 ug/mlllion cells, Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 minutes followed by cooling at RT for 20 minutes, Optimal dilution for a specific application should be determined by user			
Positive control	Daudi, Raji and U266 cell lines. Human lymphocytes. Human lymph node or 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/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			
Cell/tissue expression	B-cells, Lymphocytes			
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.

This datasheet was generated on October 29, 2025 at 09:37:29 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd20-ms4a1-monoclonal-mouse-antibody-l26/