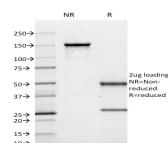


CD10 Monoclonal Mouse Antibody (MME/1620)



Product Description

CD10 (also known as Common Acute Lymphocytic Leukemia Antigen (CALLA)) is a type II transmembrane protein. It is a cell surface enzyme with neutral metalloendopeptidase activity, which inactivates a variety of biologically active peptides. CD10 is expressed on the cells of lymphoblastic, Burkitt's, and follicular germinal center lymphomas, and on cells from patients with chronic myelocytic leukemia (CML). It is also expressed on the surface of normal early lymphoid progenitor cells, immature B cells within adult bone marrow and germinal center B cells within lymphoid tissue.

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 1620, CD10 Monoclonal Mouse Antibody (MME/1620)

Product attributes

Antibody number	#1620
Antibody reactivity (target)	CD10
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MME/1620
Isotype	IgG1, kappa
Molecular weight	100 kDa
Synonyms	CD10; Nephrylsin; Atriopeptidase; Common acute lymphocytic leukemia antigen (CALLA); Enkephalinase; gp100; Membrane metalloendopeptidase; Neutral endopeptidase
Human gene symbol	MME
Entrez gene ID	4311
SwissProt	P08473
Unigene	307734
Immunogen	Recombinant full-length human CD10 protein (MME)
Verified antibody applications	Flow (verified), IF (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Antibody application notes	Flow cytometry: 1-2 ug/million cells; Western blot 0.5-1 ug/mL; Immunofluorescence: 1-2 ug/mL; Optimal dilution for a specific application should be determined. Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody
Positive control	Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney (IHC).
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	Immunology, Proteases
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.
Cell/tissue expression	B-cells, Epithelial cells
Tumor expression	Leukemia/lymphoma

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 26, 2024 at 07:53:05 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cd10-monoclonal-mouse-antibody-mme-1620/>