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)

Droduct attributes

Call us: 800-304-5357

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; Neprilysin; 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			
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			
•	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			
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 Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney			
Antibody application notes Positive control	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 Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney (IHC).			
Antibody application notes Positive control Shipping condition	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 Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,			
Antibody application notes Positive control Shipping condition Storage Conditions	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 Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney (IHC). Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C			
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate	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. Ramos cells (FACS/IF). Jurkat or LnCap lysates (WB). Kidney (IHC). 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/0.05% pack purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% BS			
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation	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 (HC). Ramos cells (FACS/IF), Jurkat or LnCap lysates (WB), Kidney (IHC). 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/.05% azide, Purified; 0.2 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when			
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life	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 conjugates than for indirect detection with secondary antibody. Ramos cells (FACS/IF), Jurkat or LnCap lysates (WB). Kidney (IHC). 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 stored as recommended			
Antibody application notes Positive control Shipping condition Storage Conditions Regulatory status Antibody/conjugate formulation Shelf life Antibody research areas	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 conjugates than for indirect detection with secondary antibody (IHC). Romot 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, Purified: 0.2 mg/mL in PBS without azide Guaranteed for at least 24 months from date of receipt when stored as recommended Immunology, Proteases Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in			

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 30, 2025 at 05:10:06 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/cd10-monoclonal-mouse-antibody-mme-1620/