CD4 Monoclonal Mouse Antibody (C4/206)



Product Description

Recognizes a protein of 55 kDa, identified as CD4. CD4 is a membrane glycoprotein of T lymphocytes that interacts with major histocompatibility complex class II antigens and is also a receptor for the human immunodeficiency virus. This protein is expressed not only in T lymphocytes, but also in B cells, macrophages, and granulocytes. It is also expressed in specific regions of the brain. The protein functions to initiate or augment the early phase of T-cell activation, and may function as an important mediator of indirect neuronal damage in infectious and immune-mediated diseases of the central nervous system. Multiple alternatively spliced transcript variants encoding different isoforms have been identified.

Also view our <u>CD4 Monoclonal Mouse Biotium Choice Antibody (SK3)</u>. The antibody belongs to the Biotium Choice list of select antibodies that have been developed and validated in-house for optimal performance. The antibody has been validated in flow cytometry.

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 0206, Anti-CD4 (C4/206)

Product attributes

Call us: 800-304-5357

Product attributes			
Antibody number	#0206		
Antibody reactivity (target)	CD4		
Antibody type	Primary		
Host species	Mouse		
Clonality	Monoclonal		
Clone	C4/206		
Isotype	IgG1, kappa		
Molecular weight	55 kDa		
Synonyms	L3T4; Leu3; Ly-4; Lymphocyte antigen CD4; p55; T cell antiger T4/LEU3; T cell differentiation antigen L3T4; T-cell surface antigen T4/Leu-3; T-cell surface glycoprotein CD4		
Human gene symbol	CD4		
Entrez gene ID	920		
SwissProt	P01730		
Unigene	631659		
Immunogen	Recombinant human CD4 protein		
Antibody target cellular localization	Plasma membrane		
Verified antibody applications	Flow (surface) (verified), IF (verified)		
Species reactivity	Human		
	orimary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a		
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user		
Antibody application notes Positive control	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a		
	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/ml., Flow Cytometry 0.5-1 ug/mlion cells/0.1 ml., Optimal dilution for a specific application should be determined by user Human peripheral blood mononuclear cells (PBMCs),		
Positive control	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/ml., Flow Cytometry 0.5-1 ug/mlion cells/0.1 ml., Optimal dilution for a specific application should be determined by user Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60. Lymph node or tonsil.		
Positive control Shipping condition	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mlL indirection appetition appetition appetition and the determined by user Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60. Lymph node or tonsil. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light,		
Positive control Shipping condition Storage Conditions	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/ml., Flow Cytometry 0.5-1 ug/mllion cells/0.1 ml., Optimal dilution for a specific application should be determined by user Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60. Lymph node or tonsil. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when		
Positive control Shipping condition Storage Conditions Shelf life	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mlL indirect of the specific application should be determined by user Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60. Lymph node or tonsil. Room temperature Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C Guaranteed for at least 24 months from date of receipt when stored as recommended		
Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/ml., Flow Cytometry 0.5-1 ug/ml. Indiverse of the property of th		
Positive control Shipping condition Storage Conditions Shelf life Regulatory status Antibody/conjugate formulation	primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 0.5-1 ug/mlL, Flow Cytometry 0.5-1 ug/mlL indiversed to the properties of		

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 05:12:19 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-cd4-c4206/