CD6 Monoclonal Mouse Antibody (SPV-L14)

Product Description

CD6 is a type I transmembrane glycoprotein that contains a 24-amino acid signal sequence, three extracellular scavenger receptor cysteine-rich (SRCR) domains, a membrane-spanning domain and a 44-amino acid cytoplasmic domain. The CD6 glycoprotein is tyrosine phosphorylated during TCR-mediated T cell activation. CD6 shows significant homology to CD5. CD6 is present on mature thymocytes, peripheral T cells and a subset of B cells. Antibodies to CD6 are used to deplete T cells from bone marrow transplants to prevent graft versus host disease. Catalog number key for antibody number 0402, Anti-CD6 (SPV-L14)

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

Product attributes

Product attributes				
Antibody number	#0402			
Antibody reactivity (target)	CD6			
Antibody type	Primary			
Host species	Mouse			
Clonality	Monoclonal			
Clone	SPV-L14			
Isotype	IgG1, kappa			
Molecular weight	90-130 kDa			
Synonyms	T12; TP120			
Human gene symbol	CD6			
Entrez gene ID	923			
SwissProt	P30203			
Unigene	744366			
Immunogen	CD8+ cytotoxic T-cell clone			
Antibody target cellular localization	Plasma membrane			
Verified antibody applications	Flow (verified), IF (verified)			
Species reactivity	Human, Mouse			
Expected antibody applications	Flow, surface (published for clone), Functional studies (published for clone), IF (published for clone)			
Antibody application notes	Flow cytometry: 1-2 ug/mL, 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, Optimal dilution for a specifiapplication should be determined by user			
Positive control	CCRF-CEM, Jurkat cells, 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 i PBS without azide			
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended			
Cell/tissue expression	B-cells, T-cells			
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.			
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.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

- 1. PNAS USA (1987) 84: 3861-3865. (Flow, surface; IF)
- 2. J Inv Dermatol (1994) 102(6): 833-837. (Flow, surface) 3. J Immunol (2006) 177:1152-1159. (functional studies)
- 4. Nat Immunol (2014) 15(4): 384-392. (functional studies)