

CD4 Monoclonal Mouse Antibody (SK3) - Biotium Choice

CD4 (clone SK3) is a validated mouse monoclonal antibody that recognizes CD4, a membrane glycoprotein of T lymphocytes. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.



Product Description

CD4 (clone SK3) is a mouse monoclonal antibody that recognizes the extracellular domain of CD4, a membrane glycoprotein of T lymphocytes. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance. The antibody is available conjugated to CF® Dyes and Astral™ tandem dyes. They are supplied in PBS, 0.1% BSA, 0.05% azide.

CD4 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. The majority of peripheral T-cell lymphomas are derived from the T-helper/regulatory cell subset so that most mature T-cell neoplasms are CD4/CD8. Anti-CD4 is used in the immunohistochemical staining of lymphoproliferative disorders to evaluate tumors with CD4 aberrant expression.

As an isotype control, we recommend [Isotype Control, Monoclonal Mouse IgG1 Kappa \(IGG1/1331\)](#) for this antibody. [View our full catalog of Biotium Choice antibodies](#), optimized and validated for the best results in flow cytometry.

Product attributes

Antibody number	#P001
Antibody reactivity (target)	CD4
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SK3
Isotype	IgG1, kappa
Molecular weight	55 kDa
Synonyms	L3T4; Leu3; Ly-4; Lymphocyte antigen CD4; p55; T cell antigen 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	Human peripheral blood T cells
Antibody target cellular localization	Plasma membrane
Verified antibody applications	Flow (surface) (verified)
Species reactivity	Human
Antibody application notes	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
Positive control	Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/0.1% BSA/0.05% azide
Antibody research areas	Immunology
Volume per assay	5 uL/test
Cell/tissue expression	Lymphocytes, Myeloid cells
Tumor expression	Leukemia/lymphoma
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (<i>Bos taurus</i>), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

CD4 Monoclonal Mouse Antibody (SK3) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF@405S	411/431 nm	100 ug/mL	25 tests (125 uL)	P001-405S-125	CF@405S Features
			100 tests (500 uL)	P001-405S-500	
CF@405M	416/452 nm	100 ug/mL	25 tests (125 uL)	P001-405M-125	CF@405M Features
			100 tests (500 uL)	P001-405M-500	
CF@405L	413/547 nm	200 ug/mL	25 tests (125 uL)	P001-405L-125	CF@405L Features
			100 tests (500 uL)	P001-405L-500	
CF@488A	490/516 nm	100 ug/mL	25 tests (125 uL)	P001-488A-125	CF@488A Features
			100 tests (500 uL)	P001-488A-500	
CF@503R	503/532 nm	50 ug/mL	25 tests (125 uL)	P001-503R-125	CF@503R Features
			100 tests (500 uL)	P001-503R-500	
CF@514	516/549 nm	100 ug/mL	25 tests (125 uL)	P001-514-125	CF@514 Features
			100 tests (500 uL)	P001-514-500	
CF@ 568	562/584 nm	50 ug/mL	25 tests (125 uL)	P001-568-125	CF@568 Features
			100 tests (500 uL)	P001-568-500	
CF@ 575	575/601 nm	25 ug/mL	25 tests (125 uL)	P001-575-125	
			100 tests (500 uL)	P001-575-500	
RPE-Astral™616	496, 546, 566/617 nm	100 ug/mL	25 tests (125 uL)	P001-R616-125	
			100 tests (500 uL)	P001-R616-500	
CF@647Plus	652/668 nm	25 ug/mL	25 tests (125 uL)	P001-647P-125	
			100 tests (500 uL)	P001-647P-500	
CF@660C	667/685 nm	50 ug/mL	25 tests (125 uL)	P001-660C-125	CF@660C Features
			100 tests (500 uL)	P001-660C-500	
CF@700	696/721 nm	25 ug/mL	25 tests (125 uL)	P001-700-125	CF@700 Features
			100 tests (500 uL)	P001-700-500	
APC-Astral™813	633, 638/813 nm	50 ug/mL	25 tests (125 uL)	P001-A813-125	
			100 tests (500 uL)	P001-A813-500	
R-PE	496, 546, 565/576 nm	25 ug/mL	25 tests (125 uL)	P001-RPE-125	
			100 tests (500 uL)	P001-RPE-500	
APC	651/660 nm	50 ug/mL	25 tests (125 uL)	P001-APC-125	
			100 tests (500 uL)	P001-APC-500	

This datasheet was generated on November 21, 2024 at 02:32:47 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/cd4-monoclonal-mouse-antibodyh-sk3/>