

## 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

<b>Antibody number</b>	#P001
<b>Antibody reactivity (target)</b>	CD4
<b>Biotium Choice Antibody</b>	Primary
<b>Antibody type</b>	Biotium Choice Flow Antibodies, Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	SK3
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	55 kDa
<b>Synonyms</b>	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
<b>Human gene symbol</b>	CD4
<b>Entrez gene ID</b>	920
<b>SwissProt</b>	P01730
<b>Unigene</b>	631659
<b>Immunogen</b>	Human peripheral blood T cells
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Verified antibody application</b>	Flow (surface) (verified)
<b>Species reactivity</b>	Human
<b>Antibody application notes</b>	Recommended concentration for flow cytometry: 5 uL per million cells/0.1 mL
<b>Positive control</b>	Human peripheral blood mononuclear cells (PBMCs), CCRF-CEM or HL-60
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	Conjugates: PBS/0.1% BSA/0.05% azide
<b>Antibody research areas</b>	Immunology
<b>Volume per assay</b>	5 uL/test
<b>Cell/tissue expression</b>	Lymphocytes, Myeloid cells
<b>Tumor expression</b>	Leukemia/lymphoma

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

This datasheet was generated on June 30, 2024 at 05:58:36 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/cd4-monoclonal-mouse-antibodyh-sk3/>