

## CD8a Monoclonal Mouse Antibody (SK1) - Biotium Choice



CD8a (clone SK1) is a validated mouse monoclonal antibody that recognizes CD8a, a transmembrane glycoprotein and co-receptor found in T cells. The antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for optimal performance.

### Product Description

CD8a (clone SK1) is a mouse monoclonal antibody derived from the CD8 alpha chain. 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.

CD8 is a cell surface receptor expressed either as a heterodimer with the CD8 beta chain (CD8 alpha/beta) or as a homodimer (CD8 alpha/alpha). A majority of thymocytes and a subpopulation of mature T cells and NK cells express CD8a. CD8 binds to MHC class 1 and through its association with protein tyrosine kinase p56lck plays a role in T cell development and activation of mature T cells. For mature T-cells, CD4 and CD8 are mutually exclusive, so anti-CD8, generally used in conjunction with anti-CD4. It is a useful marker for distinguishing helper/inducer T-lymphocytes, and most peripheral T-cell lymphomas are CD4 /CD8-. Anaplastic large cell lymphoma is usually CD4 and CD8-, and in T-lymphoblastic lymphoma/leukemia, CD4 and CD8 are often co-expressed. CD8 is also found in littoral cell angiomata of the spleen.

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>	#P009
<b>Antibody reactivity (target)</b>	CD8A
<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>	SK1
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight</b>	32-34 kDa (alpha chain)
<b>Synonyms</b>	CD8, Leu2, MAL, p32
<b>Human gene symbol</b>	CD8A
<b>Entrez gene ID</b>	925
<b>SwissProt</b>	P01732
<b>Unigene</b>	85258
<b>Immunogen</b>	Human peripheral blood T cells
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Verified antibody applications</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, T-cells
<b>Tumor expression</b>	Leukemia/lymphoma
<b>Product origin</b>	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.

## CD8a Monoclonal Mouse Antibody (SK1) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	
<a href="#">CF@405S</a>	411/431 nm	200 ug/mL	25 tests (125 uL)	<a href="#">P009-405S-125</a>	<a href="#">CF@405S Features</a>
			100 tests (500 uL)	<a href="#">P009-405S-500</a>	
<a href="#">CF@405M</a>	416/452 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-405M-125</a>	<a href="#">CF@405M Features</a>
			100 tests (500 uL)	<a href="#">P009-405M-500</a>	
<a href="#">CF@405L</a>	413/547 nm	300 ug/mL	25 tests (125 uL)	<a href="#">P009-405L-125</a>	<a href="#">CF@405L Features</a>
			100 tests (500 uL)	<a href="#">P009-405L-500</a>	
<a href="#">CF@488A</a>	490/516 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-488A-125</a>	<a href="#">CF@488A Features</a>
			100 tests (500 uL)	<a href="#">P009-488A-500</a>	
<a href="#">CF@503R</a>	503/532 nm	50 ug/mL	25 tests (125 uL)	<a href="#">P009-503R-125</a>	<a href="#">CF@503R Features</a>
			100 tests (500 uL)	<a href="#">P009-503R-500</a>	
<a href="#">CF@514</a>	516/549 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-514-125</a>	<a href="#">CF@514 Features</a>
			100 tests (500 uL)	<a href="#">P009-514-500</a>	
<a href="#">CF@568</a>	562/584 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-568-125</a>	<a href="#">CF@568 Features</a>
			100 tests (500 uL)	<a href="#">P009-568-500</a>	
<a href="#">CF@575</a>	575/601 nm	200 ug/mL	25 tests (125 uL)	<a href="#">P009-575-125</a>	
			100 tests (500 uL)	<a href="#">P009-575-500</a>	
<a href="#">RPE-Astral™616</a>	496, 546, 566/617 nm	200 ug/mL	25 tests (125 uL)	<a href="#">P009-R616-125</a>	
			100 tests (500 uL)	<a href="#">P009-568-500</a>	
<a href="#">CF@647Plus</a>	652/668 nm	50 ug/mL	25 tests (125 uL)	<a href="#">P009-647P-125</a>	
			100 tests (500 uL)	<a href="#">P009-647P-500</a>	
<a href="#">CF@660C</a>	667/685 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-660C-125</a>	<a href="#">CF@660C Features</a>
			100 tests (500 uL)	<a href="#">P009-660C-500</a>	
<a href="#">CF@700</a>	696/721 nm	100 ug/mL	25 tests (125 uL)	<a href="#">P009-700-125</a>	<a href="#">CF@700 Features</a>
			100 tests (500 uL)	<a href="#">P009-700-500</a>	
<a href="#">CF@820</a>	822/835 nm	40 ug/mL	25 tests (125 uL)	<a href="#">P009-820-125</a>	<a href="#">CF@820 Features</a>
			100 tests (500 uL)	<a href="#">P009-820-500</a>	
<a href="#">APC-Astral™813</a>	633, 638/813 nm	50 ug/mL	25 tests (125 uL)	<a href="#">P009-A813-125</a>	
			100 tests (500 uL)	<a href="#">P009-A813-500</a>	
<a href="#">R-PE</a>	496, 546, 565/576 nm	60 ug/mL	25 tests (125 uL)	<a href="#">P009-RPE-125</a>	
			100 tests (500 uL)	<a href="#">P009-RPE-500</a>	
<a href="#">APC</a>	651/660 nm	50 ug/mL	25 tests (125 uL)	<a href="#">P009-APC-125</a>	
			100 tests (500 uL)	<a href="#">P009-APC-500</a>	

This datasheet was generated on March 28, 2025 at 01:59:22 PM. Visit product page to check for updated information before use.  
 Product link: <https://biotium.com/product/cd8a-monoclonal-mouse-antibodyh-sk1/>