

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 angioma 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

Antibody number	#P009
Antibody reactivity (target)	CD8A
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SK1
Isotype	IgG1, kappa
Molecular weight	32-34 kDa (alpha chain)
Synonyms	CD8, Leu2, MAL, p32
Human gene symbol	CD8A
Entrez gene ID	925
SwissProt	P01732
Unigene	85258
Immunogen	Human peripheral blood T cells
Antibody target cellular localization	Plasma membrane
Verified antibody	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, T-cells
Tumor expression	Leukemia/lymphoma

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

This datasheet was generated on July 17, 2024 at 09:34:46 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cd8a-monoclonal-mouse-antibodyh-sk1/>