

CD19 Monoclonal Mouse Antibody (SJ25C1) - Biotium Choice



CD19 (clone SJ25C1) is a validated mouse monoclonal antibody that recognizes CD19, a transmembrane glycoprotein in B 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

CD19 (clone SJ25C1) is a mouse monoclonal antibody that recognizes the transmembrane glycoprotein CD19. 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.

CD19 is a transmembrane glycoprotein that contains two extracellular immunoglobulin-like domains. CD19 is present in both benign and malignant B-cells and is considered to be the most reliable surface marker of this lineage over a wide range of maturational stages. In normal lymphoid tissue, CD19 is observed in germinal centers, in mantle zone cells, and in scattered cells of the inter-follicular areas. Anti-CD19 exhibits an overall immunoreactivity pattern similar to those of the antibodies against CD20 and CD22. However, in contrast to CD20, expression of CD19 is continuous throughout B-cell development and through terminal differentiation of B-cells into plasma cells. Anti-CD19 positivity is seen in the vast majority of B-cell neoplasms commonly at a lower intensity than normal B-cell counterparts. Plasma cell neoplasms are nearly always negative, as are T-cell neoplasms.

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	#P010
Antibody reactivity (target)	CD19
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Flow Antibodies, Primary
Host species	Mouse
Clonality	Monoclonal
Clone	SJ25C1
Isotype	IgG1, kappa
Molecular weight	95 kDa
Synonyms	B-Lymphocytes Surface Antigen B4; T-cell Surface Antigen Leu-12; CVID3; B4
Human gene symbol	CD19
Entrez gene ID	930
SwissProt	P15391
Unigene	652262
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)
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
Cell/tissue expression	B-cells, Lymphocytes
Volume per assay	5 uL/test
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.

CD19 Monoclonal Mouse Antibody (SJ25C1) - Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF@405M	416/452 nm	100 ug/mL	25 tests (125 uL)	P010-405M-125	CF@405M Features
			100 tests (500 uL)	P010-405M-500	
CF@488A	490/516 nm	100 ug/mL	25 tests (125 uL)	P010-488A-125	CF@488A Features
			100 tests (500 uL)	P010-488A-500	
CF@575	575/601 nm	200 ug/mL	25 tests (125 uL)	P010-575-125	
			100 tests (500 uL)	P010-575-500	
RPE-Astral™616	496, 546, 566/617 nm	50 ug/mL	25 tests (125 uL)	P010-R616-125	
			100 tests (500 uL)	P010-R616-500	
CF@647Plus	652/668 nm	25 ug/mL	25 tests (125 uL)	P010-647P-125	
			100 tests (500 uL)	P010-647P-500	
CF@700	696/721 nm	100 ug/mL	25 tests (125 uL)	P010-700-125	CF@700 Features
			100 tests (500 uL)	P010-700-500	
APC-Astral™813	633, 638/813 nm	50 ug/mL	25 tests (125 uL)	P010-A813-125	
			100 tests (500 uL)	P010-A813-500	
R-PE	496, 546, 565/576 nm	50 ug/mL	25 tests (125 uL)	P010-RPE-125	
			100 tests (500 uL)	P010-RPE-500	

This datasheet was generated on December 21, 2024 at 07:21:51 PM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/cd19-monoclonal-mouse-antibody-sj25c1/>