

iFluor™ 514 Anti-human CD19 Antibody
SJ25C1Catalog number: 10191060, 10191061
Unit size: 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	CD19 (B4)
Clone	SJ25C1
Conjugate	iFluor™ 514

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 514 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 514
Excitation Wavelength	528 nm
Emission Wavelength	555 nm

Applications

SJ25C1 is an anti-human monoclonal antibody that forms an immune complex with the CD19 antigen. CD19 (sometimes called T-cell surface antigen Leu-12 or B-lymphocyte surface antigen B4) is a 95 kD transmembrane glycoprotein that is located on the surface of cells such as stem cells, B cells and dendritic cells. CD19 is a member of key cellular pathways, in particular, the antigen receptor-mediated signaling pathway and B cell receptor signaling pathway. Additionally, in many organisms, it upregulates release of sequestered calcium ion into cytosol, promotes protein kinase B signaling and is a promoter of phosphatidylinositol 3-kinase activity. From a research standpoint, it is of biological interest due

to its association with key macromolecules/ligands such as CD225. CD19 is a very popular antibody target, with over 36000 publications in the last decade. CD19 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to iFluor™ 514 (ex/em = 528/555 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).