

PacBlue Anti-human CD19 Antibody
SJ25C1Catalog number: 101911J0, 101911J1
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	PacBlue

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).