

XFD750 Anti-human CD19 Antibody
***SJ25C1, XFD750 Same Structure to Alexa**
Fluor™ 750*

Catalog number: 101911B0, 101911B1
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	AF750

Biological Properties

Appearance	liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF750
Excitation Wavelength	752 nm
Emission Wavelength	776 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 XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher).