

**iFluor™ 568 Anti-human CD19 Antibody**  
**\*SJ25C1\***Catalog number: 101910B0, 101910B1  
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™ 568

**Biological Properties**

---

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	iFluor™ 568
Excitation Wavelength	568 nm
Emission Wavelength	587 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™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).