

**XFD488 Anti-human CD19 Antibody \*HI19a,  
XFD488 Same Structure to Alexa Fluor™  
488\***Catalog number: 10190150, 10190151  
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
Isotype	Mouse IgG1
Immunogen	CD19 (B4)
Clone	HI19a
Conjugate	AF488

**Biological Properties**

---

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 nm

**Applications**

---

HI19a is an anti-human monoclonal antibody that targets the CD19 antigen. CD19 (also known as T-cell surface antigen Leu-12 or B-lymphocyte surface antigen B4) is a 95 kD transmembrane glycoprotein that is found on the surface of cells like B cells and stem cells. CD19 plays a role in

critical cellular pathways, in particular, the antigen receptor-mediated signaling pathway and B cell receptor signaling pathway. Furthermore, in certain organisms, it plays a role in the upregulation of phosphatidylinositol 3-kinase activity, positively regulates release of sequestered calcium ion into cytosol and plays a role in the upregulation of protein kinase B signaling. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as CD81 and Leu-13. CD19 is a very popular antibody target, with over 36000 publications in the last decade. CD19 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 488 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 488 nm laser and 525/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).