

**XFD680 Anti-human CD19 Antibody *HIB19,
XFD680 Same Structure to Alexa Fluor™
680***Catalog number: 10192190, 10192191
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	HIB19
Conjugate	AF680

Biological Properties

Appearance	liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF680 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF680
Excitation Wavelength	681 nm
Emission Wavelength	704 nm

Applications

HIB19 is an anti-human monoclonal antibody that recognizes the CD19 antigen. CD19 (also known as CVID3) is a 95 kD glycoprotein that is located on the surface of cells like stem cells, dendritic cells and B cells. In certain organisms, CD19 promotes release of sequestered calcium ion into cytosol, plays a role in the upregulation of phosphatidylinositol 3-kinase activity and acts to positively regulate protein kinase B signaling. In addition, it is a member of critical cellular pathways, for example, the antigen receptor-mediated signaling pathway and B cell receptor signaling

pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as lyn. 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, specifically in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 680 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).