

PE/XFD700 Anti-human CD19 Antibody
***HIB19, XFD700 Same Structure to Alexa**
Fluor™ 700*

Catalog number: 101921P0, 101921P1, 101921P2
Unit size: 25 tests, 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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/AF700
Excitation Wavelength	566 nm
Emission Wavelength	721 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 700 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 720/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).