

**XFD350 Anti-human CD5 Antibody *L17F12,
XFD350 Same Structure to Alexa Fluor™
350***Catalog number: 10051140, 10051141
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	CD5 (Leu1, Ly-1, T1, Tp67)
Clone	L17F12
Conjugate	AF350

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF350
Excitation Wavelength	343 nm
Emission Wavelength	441 nm

Applications

L17F12 is an anti-human monoclonal antibody that is specific for the CD5 antigen. CD5 (also known as Lyt-1 or Ly-1) is a 67 kD single-pass type I membrane protein that is found on the surface of cells such as B cells and T cells. CD5 is a component of important cellular pathways, namely, the apoptotic signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as gp35-37, TCR and CD21. CD5 is a moderately popular antibody target, with over 11000 publications in the last

decade. CD5 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 350 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).