

**XFD700 Anti-human CD5 Antibody *HISM2,
XFD700 Same Structure to Alexa Fluor™
700***Catalog number: 100501A0, 100501A1
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	CD5 (Leu1, Ly-1, T1, Tp67)
Clone	HISM2
Conjugate	AF700

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF700
Excitation Wavelength	696 nm
Emission Wavelength	719 nm

Applications

HISM2 is an anti-human monoclonal antibody that forms an immune complex with the CD5 antigen. CD5 (sometimes referred to as LEU1, Tp67 or Lyt-1) is a 67 kD single-pass type I membrane protein that is located on the surface of cells like B cells and T cells. CD5 is a component of vital

cellular pathways, for instance, the apoptotic signaling pathway. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like ZAP-70, TCR and BCR. CD5 is a moderately popular antibody target, with over 11000 publications in the last decade. CD5 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to XFD700 (ex/em = 696/719 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).