

XFD750 Anti-human CD72 Antibody *3F3*

Catalog Number: 107201B0,

107201B1

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	Lot specific (please consult certificate of analysis for given lot)
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 IgG2b
Immunogen	CD72 (Ly-19.2, Ly-32.2, Lyb-2)
Clone	3F3
Conjugate	AF750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF750
-----------	-------

Excitation Wavelength 752 nm

Emission Wavelength 776 nm

Applications

The 3F3 monoclonal antibody binds with human CD72, a 39 - 43 kD transmembrane glycoprotein often located on the surface of B cells, dendritic cells, liver kupffer cells and macrophages. CD72 is associated with a variety of biologically interesting macromolecules/ligands, for example, CD5. CD72 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD72 is commonly 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 XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).