

XFD750 Anti-human CD56 Antibody *My31*

Catalog Number: 105621A0, 105621A1

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 IgG1, κ
Immunogen	CD56 (Leu-19, NKH1, NCAM1)
Clone	My31
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

My31 is an anti-human monoclonal antibody that is specific for the CD56 antigen. CD56 (sometimes referred to as NCAM1, NKH1 or Leu-19) is a single-pass type I membrane protein that is found on the surface of cells like NK cells and T cells. CD56 is a component of key cellular pathways, namely, the interferon-gamma-mediated signaling pathway and regulation of semaphorin-plexin signaling pathway. Also, it has been closely linked to key biological processes like axon guidance, particularly commissural neuron axon guidance. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as NCAM-1. CD56 is a moderately popular antibody target, with over 18000 publications in the last decade. CD56 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of . 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).