

PE/XFD610 Anti-human CD56 Antibody *My31*

Catalog Number: 105621N0,
105621N1, 105621N2
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	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	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L 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	PE/AF610
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 627 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 PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).