

XFD680 Anti-human CD56 Antibody *B-A19*

Catalog Number: 10560190, 10560191

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
Immunogen	CD56 (Leu-19, NKH1, NCAM1)
Clone	B-A19
Conjugate	AF680

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF680 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	AF680
Excitation Wavelength	681 nm
Emission Wavelength	704 nm

Applications

B-A19 is an anti-human monoclonal antibody that targets the CD56 antigen. CD56 (alternatively called NKH1 or NCAM-1) is a single-pass type I membrane protein that is expressed on the surface of cells like NK cells and T cells. CD56 has been associated with important biological processes like axon guidance, especially commissural neuron axon guidance. In addition, it plays a role in essential cellular pathways, for instance, the regulation of semaphorin-plexin signaling pathway and interferon-gamma-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as Heparin sulfate. CD56 is a fairly uncommon antibody target, with a little more than 10000 publications in the last decade. Even still, CD56 has a variety of applications in research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 680 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).