

**XFD555 Anti-human CD56 Antibody *B-A19,
XFD555 Same Structure to Alexa Fluor™
555***Catalog number: 10560160, 10560161
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	CD56 (Leu-19, NKH1, NCAM1)
Clone	B-A19
Conjugate	AF555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF555
Excitation Wavelength	553 nm
Emission Wavelength	568 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 XFD555 (ex/em = 553/568 nm). XFD555 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 555 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSMelody™).