

**XFD488 Anti-human CD56 Antibody \*AHC0060\***

Catalog Number: 10561150, 10561151

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	IgG1, kappa
Immunogen	CD56 (Leu-19, NKH1, NCAM1)
Clone	AHC0060
Conjugate	AF488

**Biological Properties**

---

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 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	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 nm

## Applications

---

AHC0060 is an anti-human monoclonal antibody that is specific for the CD56 antigen. CD56 (sometimes referred to as NCAM-1, Leu-19 or NCAM1) is a member of the Ig superfamily that is expressed on the surface of cells such as T cells and NK cells. CD56 has been associated with important biological processes like axon guidance, specifically commissural neuron axon guidance. Additionally, it is involved with important cellular pathways, for instance, the interferon-gamma-mediated signaling pathway and regulation of semaphorin-plexin signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like NCAM-1 and heparin sulfate. CD56 is a moderately popular antibody target, with over 18000 publications in the last decade. CD56 is essential for research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSCanto™ II).