

iFluor™ 405 Anti-human CD56 Antibody *B-A19*Catalog number: 10560020, 10560021
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	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 445/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).