

**PE/Cy7 Anti-human CD56 Antibody \*B-A19\***Catalog number: 105601N0, 105601N1, 105601N2  
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	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	PE/Cy7

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 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 PE/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 762/35 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).