

iFluor™ A7 Anti-human CD56 Antibody *B-A19*Catalog number: 105600S0, 105600S1
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™ A7 |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------|------------|
| Conjugate | iFluor™ A7 |
|-----------|------------|

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™ A7.