

**PE/iFluor™ 594 Anti-human CD56 Antibody**  
**\*B-A19\***Catalog number: 105601Y0, 105601Y1, 105601Y2  
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/iFluor™ 594             |

**Biological Properties**

|             |                                                                                                               |
|-------------|---------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 594 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                         |

**Spectral Properties**

|                       |                |
|-----------------------|----------------|
| Conjugate             | PE/iFluor™ 594 |
| Excitation Wavelength | 566 nm         |
| Emission Wavelength   | 606 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/iFluor™ 594 (ex/em = 566/606 nm). It is compatible with the 561 nm laser and 611/31 nm bandpass filter (for example, as in the Luminex Amnis CellStream).