

**mFluor™ Red 700 Anti-human CD56  
Antibody \*My31\***Catalog number: 105620V0, 105620V1  
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, κ              |
| Immunogen          | CD56 (Leu-19, NKH1, NCAM1) |
| Clone              | My31                       |
| Conjugate          | mFluor™ Red 700            |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 700 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                    |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | mFluor™ Red 700 |
| Excitation Wavelength | 680 nm          |
| Emission Wavelength   | 695 nm          |

**Applications**

My31 is an anti-human monoclonal antibody that is specific for the CD56 antigen. CD56 (sometimes referred to as NCAM1, NKH1 or Leu-19) is a single-pass type I membrane protein that is found on the surface of cells like NK cells and T cells. CD56 is a component of key cellular pathways, namely, the interferon-gamma-mediated signaling pathway and regulation of semaphorin-plexin signaling pathway. Also, it has been closely

linked to key biological processes like axon guidance, particularly commissural neuron axon guidance. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as NCAM-1. CD56 is a moderately popular antibody target, with over 18000 publications in the last decade. CD56 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of . This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).