

PE/Texas Red® Anti-human CD56 Antibody
My31Catalog number: 105621R0, 105621R1, 105621R2
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, κ
Immunogen	CD56 (Leu-19, NKH1, NCAM1)
Clone	My31
Conjugate	PE/Texas Red®

Biological Properties

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

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 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 PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 611/31 nm bandpass filter (for example, as in the Luminex Amnis CellStream).