

**PE/Cy5 Anti-human CD31 Antibody**  
**\*MEM-05\***Catalog number: 103101N0, 103101N1, 103101N2  
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	CD31 (PECAM-1, EndoCAM, Platelet endothelial cell adhesion molecule)
Clone	MEM-05
Conjugate	PE/Cy5

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	PE/Cy5
Excitation Wavelength	565 nm
Emission Wavelength	666 nm

**Applications**

---

The MEM-05 monoclonal antibody reacts with human CD31, a 130 - 140 kD transmembrane glycoprotein commonly found on the surface of monocytes and endothelial cells. CD31 has been closely linked to key biological processes like cell-cell adhesion, particularly leukocyte cell-cell adhesion. In addition, in certain organisms, it is an enhancer of peptidyl-tyrosine phosphorylation, plays a role in the upregulation of cell migration and plays a role in the upregulation of protein localization to cell-cell junction. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as CD38. CD31 is a very popular antibody target, with over 32000 publications in the last decade. CD31 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the

study of angiogenesis and immunology. This antibody was purified through affinity chromatography and conjugated to PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 660/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).