

FITC Anti-human CD31 Antibody *MEM-05*Catalog number: 10310110, 10310111
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	CD31 (PECAM-1, EndoCAM, Platelet endothelial cell adhesion molecule)
Clone	MEM-05
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 525/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).