

**FITC Anti-non-human primates/ human
CD177 Antibody *MEM-166***Catalog number: 117701G0, 117701G1
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD177 (NB1gp, HNA-2a, NB1, Neutrophil-specific antigen 1, PRV1)
Clone	MEM-166
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

MEM-166 is an anti-non-human primates/ human monoclonal antibody that targets the CD177 antigen. CD177 (sometimes referred to as Neutrophil-specific antigen 1, PRV1 or NB1gp) is a 60 kD member of the uPAR family that is expressed on the surface of cells like granulocytes. CD177 plays a role in critical cellular pathways, namely, the regulation of integrin-mediated signaling pathway. In addition, it has been associated with vital biological processes such as leukocyte cell-cell adhesion, particularly cell-cell adhesion via plasma-membrane adhesion molecules. In some organisms, CD177 is an enhancer of superoxide anion generation, and is associated with a variety of biologically interesting

macromolecules/ligands. CD177 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD177 is essential for immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 528/65 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).