

**APC/Cy7 Anti-non-human primates/ human  
CD177 Antibody \*MEM-166\***Catalog number: 117701C0, 117701C1, 117701C2  
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	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	APC/Cy7

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

Conjugate	APC/Cy7
Excitation Wavelength	754 nm
Emission Wavelength	779 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 APC/Cy7 (ex/em = 754/779 nm).