

PE/XFD700 Anti-non-human primates/ human CD177 Antibody *MEM-166*

Catalog Number: 117701O0,
117701O1, 117701O2
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	Lot specific (please consult certificate of analysis for given lot)
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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 695/40 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).