

**APC/Cy7 Anti-human CD108 Antibody**  
**\*MEM-150\***Catalog number: 110801C0, 110801C1, 110801C2  
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 IgM
Immunogen	CD108 (JM8 blood group antigen, semaphorin 7A)
Clone	MEM-150
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**

The MEM-150 monoclonal antibody binds with human CD108, a 80 kD transmembrane glycoprotein often found on the surface of thymus and T cells. CD108 is associated with a variety of biologically interesting macromolecules/ligands, for example, CD232 and tyrosine kinases. CD108 is a relatively rare antibody target, with fewer than 50 publications in the last decade. Even still, CD108 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to APC/Cy7 (ex/em = 754/779 nm).