

**mFluor™ Violet 540 Anti-human CD108  
Antibody \*MEM-150\***Catalog number: 11080120, 11080121  
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 IgM
Immunogen	CD108 (JMH blood group antigen, semaphorin 7A)
Clone	MEM-150
Conjugate	mFluor™ Violet 540

**Biological Properties**

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 540 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	mFluor™ Violet 540
Excitation Wavelength	394 nm
Emission Wavelength	537 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 mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 510/50 nm bandpass filter (for example, as in the BD FACSaria™ III).