

iFluor™ 670 Anti-human CD108 Antibody
MEM-150Catalog number: 110800H0, 110800H1
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	iFluor™ 670

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 670 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 670
Excitation Wavelength	671 nm
Emission Wavelength	682 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 iFluor™ 670 (ex/em = 671/682 nm). It is compatible with the 642 nm laser and 702/87 nm bandpass

filter (for example, as in the Luminex Amnis CellStream).