

**PerCP Anti-human CD108 Antibody**  
**\*MEM-150\***Catalog number: 11080150, 11080151, 11080152  
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 (JMH blood group antigen, semaphorin 7A) |
| Clone              | MEM-150                                        |
| Conjugate          | PerCP                                          |

**Biological Properties**

|             |                                                                                                      |
|-------------|------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PerCP  |
| Excitation Wavelength | 477 nm |
| Emission Wavelength   | 678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 692/80 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).