

**PE Anti-mouse/ human CD59 Antibody**  
**\*MEM-43/5\***Catalog number: 105901M0, 105901M1, 105901M2  
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 | Mouse, human                                     |
| Class              | Primary                                          |
| Clonality          | Monoclonal                                       |
| Host               | Mouse                                            |
| Isotype            | Mouse IgG2b                                      |
| Immunogen          | CD59 (Protectin, H19, 1F-5Ag, MACIF, M1RL, P-18) |
| Clone              | MEM-43/5                                         |
| Conjugate          | PE                                               |

**Biological Properties**

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

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PE     |
| Excitation Wavelength | 566 nm |
| Emission Wavelength   | 574 nm |

**Applications**

The MEM-43/5 monoclonal antibody binds to mouse/ human CD59, a 19 - 25 kD member of the Ly6 superfamily frequently found on the surface of all cell types. CD59 is associated with a variety of biologically interesting macromolecules/ligands, in particular, fyn, c8- $\alpha$  and lck. CD59 is a fairly uncommon antibody target, with a little more than 3900 publications in the last decade. Even still, CD59 is vital to costimulatory molecules and neuroinflammation research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSMelody™).