

**PE/iFluor™ 647 Anti-mouse/ human CD59
Antibody *MEM-43/5***Catalog number: 105901S0, 105901S1, 105901S2
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, MIRL, P-18) |
| Clone | MEM-43/5 |
| Conjugate | PE/iFluor™ 647 |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 647 under optimal conditions |
| Application | Flow Cytometry (FACS) |

Spectral Properties

| | |
|-----------------------|----------------|
| Conjugate | PE/iFluor™ 647 |
| Excitation Wavelength | 569 nm |
| Emission Wavelength | 666 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-α and Ick. 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/iFluor™ 647 (ex/em = 569/666 nm). It is

compatible with the 561 nm laser and 697/58 nm bandpass filter (for example, as in the BD FACSMelody™).