

**PE/iFluor™ 594 Anti-mouse/ human CD59
Antibody *MEM-43/5***Catalog number: 105901Y0, 105901Y1, 105901Y2
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™ 594

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 594 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 594
Excitation Wavelength	566 nm
Emission Wavelength	606 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™ 594 (ex/em = 566/606 nm). It is

compatible with the 561 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).