

PerCP Anti-mouse/ human CD59 Antibody
MEM-43/5Catalog number: 105901V0, 105901V1, 105901V2
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	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-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 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 693/37 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant X).