

PE/iFluor™ 594 Anti-human CD72 Antibody
3F3Catalog number: 107201Y0, 107201Y1, 107201Y2
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 IgG2b
Immunogen	CD72 (Ly-19.2, Ly-32.2, Lyb-2)
Clone	3F3
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 3F3 monoclonal antibody binds with human CD72, a 39 - 43 kD transmembrane glycoprotein often located on the surface of B cells, dendritic cells, liver kupffer cells and macrophages. CD72 is associated with a variety of biologically interesting macromolecules/ligands, for example, CD5. CD72 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD72 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. 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

615/20 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).