

TRITC Anti-human CD72 Antibody *3F3*Catalog number: 107201J0, 107201J1
Unit size: 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	TRITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 586/15 nm bandpass filter (for example, as in the BD FACSCelesta™).