

FITC Anti-human CD122 Antibody *TU27*Catalog number: 112201G0, 112201G1
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 IgG1 kappa
Immunogen	CD122 (IL2R β)
Clone	TU27
Conjugate	FITC

Biological Properties

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

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

TU27 is an anti-human monoclonal antibody that is specific for the CD122 antigen. CD122 (sometimes referred to as IL2R β) is a 70 - 75 kD glycoprotein that is expressed on the surface of cells such as NK cells, T cells, macrophages and B cells. CD122 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD25, jak1, lck and IL-15. CD122 is a fairly uncommon antibody target, with a little more than 2500 publications in the last decade. Even still, CD122 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of neuroscience. This antibody was purified through affinity chromatography and conjugated to FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/40 nm bandpass filter (for example, as in the BD FACSJazz™).