

**APC/iFluor™ 750 Anti-human CD122
Antibody *TU27***Catalog number: 112201F0, 112201F1, 112201F2
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 IgG1 kappa
Immunogen	CD122 (IL2R β)
Clone	TU27
Conjugate	APC/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 750
Excitation Wavelength	754 nm
Emission Wavelength	793 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, Ick 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 APC/iFluor™ 750 (ex/em = 754/793 nm).