

**mFluor™ Blue 570 Anti-human CD122  
Antibody \*TU27\***Catalog number: 112200T0, 112200T1  
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 $\beta$ ) |
| Clone              | TU27                  |
| Conjugate          | mFluor™ Blue 570      |

**Biological Properties**

|             |                                                                                                                 |
|-------------|-----------------------------------------------------------------------------------------------------------------|
| Appearance  | Red liquid                                                                                                      |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                     |

**Spectral Properties**

|                       |                  |
|-----------------------|------------------|
| Conjugate             | mFluor™ Blue 570 |
| Excitation Wavelength | 553 nm           |
| Emission Wavelength   | 565 nm           |

**Applications**

TU27 is an anti-human monoclonal antibody that is specific for the CD122 antigen. CD122 (sometimes referred to as IL2R $\beta$ ) 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 mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZES Cell Analyzer).