

**TRITC Anti-human CD25 Antibody \*7G7B6\***Catalog number: 102511J0, 102511J1  
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 igg2a                              |
| Immunogen          | CD25 (IL-2R $\alpha$ , p55, TAC antigen) |
| Clone              | 7G7B6                                    |
| 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 7G7B6 monoclonal antibody recognizes human CD25, a 55 kD glycoprotein often located on the surface of oligodendrocytes, T cells, B cells and thymocytes. In certain organisms, CD25 negatively regulates defense response to virus, downregulates inflammatory response and positively regulates activated T cell proliferation. Additionally, it is involved with critical cellular pathways, for instance, the interleukin-2-mediated signaling pathway, cell surface receptor signaling pathway and cytokine-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like IL-2. CD25 is a very popular antibody target, with over 40000 publications in the last decade. CD25 is vital to immunology research, often serving as a phenotypic marker for

differentiating cell types in flow cytometric applications. 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 586/14 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).