

**iFluor™ 450 Anti-human CD25 Antibody
*7G7B6***Catalog number: 10251040, 10251041
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 α , p55, TAC antigen)
Clone	7G7B6
Conjugate	iFluor™ 450

Biological Properties

Appearance	Brown liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 450
Excitation Wavelength	451 nm
Emission Wavelength	502 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 iFluor™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACS Aria™ III).