

PacOrange Anti-human CD25 Antibody
7G7B6Catalog number: 102511L0, 102511L1
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	PacOrange

Biological Properties

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

Spectral Properties

Conjugate	PacOrange
Excitation Wavelength	400 nm
Emission Wavelength	551 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 PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm laser and 537/65 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).