

XFD350 Anti-human CD25 Antibody *7G7B6*

Catalog Number: 10251140, 10251141

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	Lot specific (please consult certificate of analysis for given lot)
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	AF350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF350
Excitation Wavelength	343 nm
Emission Wavelength	441 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 XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).