

**XFD647 Anti-human CD25 Antibody *HI25a,
XFD647 Same Structure to Alexa Fluor™
647***Catalog number: 10250180, 10250181
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
Immunogen	CD25 (IL-2R α , p55, TAC antigen)
Clone	HI25a
Conjugate	AF647

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF647
Excitation Wavelength	650 nm
Emission Wavelength	671 nm

Applications

HI25a is an anti-human monoclonal antibody that is specific for the CD25 antigen. CD25 (sometimes called IL2RA) is a 55 kD transmembrane protein that is found on the surface of cells such as T cells, macrophages, B cells and NK cells. In certain organisms, CD25 is involved in the

positive regulation of activated T cell proliferation, is a negative regulator of T cell proliferation and plays a role in the upregulation of T cell differentiation. Additionally, it is a component of important cellular pathways, for example, the cytokine-mediated signaling pathway, interleukin-2-mediated signaling pathway and Notch signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like IL-2. CD25 is a very popular antibody target, with over 40000 publications in the last decade. CD25 has a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 647 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).