

iFluor™ 660 Anti-human CD25 Antibody
HI25aCatalog number: 102500G0, 102500G1
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	iFluor™ 660

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 660
Excitation Wavelength	663 nm
Emission Wavelength	678 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 iFluor™ 660 (ex/em = 663/678 nm). It is compatible with the 638 nm laser and 660/20 nm bandpass filter (for example, as in the Beckman Coulter Navios EX).