

**PE/Cy5 Anti-human CD25 Antibody \*HI25a\***Catalog number: 102501M0, 102501M1, 102501M2  
Unit size: 25 tests, 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 $\alpha$ , p55, TAC antigen) |
| Clone              | HI25a                                    |
| Conjugate          | PE/Cy5                                   |

**Biological Properties**

|             |                                                                                                       |
|-------------|-------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                 |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PE/Cy5 |
| Excitation Wavelength | 565 nm |
| Emission Wavelength   | 666 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 PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 697/58 nm bandpass filter (for example, as in the BD FACSMelody™).