

**TRITC Anti-human CD21 Antibody \*HI21a\***Catalog number: 10210110, 10210111  
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          | CD21 (CR2, C3dR, Epstein-Barr virus receptor) |
| Clone              | HI21a   |
| Conjugate          | TRITC   |

**Biological Properties**

---

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

**Spectral Properties**

---

|                       |        |
|-----------------------|--------|
| Conjugate             | TRITC  |
| Excitation Wavelength | 544 nm |
| Emission Wavelength   | 570 nm |

**Applications**

---

The HI21a monoclonal antibody recognizes human CD21, a 145 kD transmembrane glycoprotein frequently expressed on the surface of thymocytes and dendritic cells. CD21 plays a role in vital cellular pathways, namely, the complement activation, classical pathway. In addition, it has been associated with vital biological processes like immune response, specifically innate immune response. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as CD19, CD23 and C3d. CD21 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD21 is essential for complement, innate immunity and 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 TRITC (ex/em = 544/570 nm). It is compatible with the 532 nm laser and 583/24 nm bandpass filter (for example, as in the Luminex Amnis CellStream).