

PE/Cy7 Anti-human CD8 Antibody *HIT8a*

Catalog number: 100801N0, 100801N1, 100801N2

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 | CD8a (T8, Leu2) |
| Clone | HIT8a |
| Conjugate | PE/Cy7 |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions |
| Application | Flow Cytometry (FACS) |

Spectral Properties

| | |
|-----------------------|--------|
| Conjugate | PE/Cy7 |
| Excitation Wavelength | 566 nm |
| Emission Wavelength | 778 nm |

Applications

The HIT8a monoclonal antibody binds to human CD8a, a 32 - 34 kD member of the Ig superfamily often located on the surface of dendritic cells, t cells and thymocytes. CD8 is a component of important cellular pathways, namely, the transmembrane receptor protein tyrosine kinase signaling pathway and cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like Lck and MHCI. CD8 is a very popular antibody target, with over 125000 publications in the last decade. CD8a is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).