

PerCP/Cy5.5 Anti-human CD8 Antibody
SK1Catalog number: 100811U0, 100811U1, 100811U2
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 |
| Immunogen | CD8a (T8, Leu2) |
| Clone | SK1 |
| Conjugate | PerCP/Cy5.5 |

Biological Properties

| | |
|-------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions |
| Application | Flow Cytometry (FACS) |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | PerCP/Cy5.5 |
| Excitation Wavelength | 489 nm |
| Emission Wavelength | 679 nm |

Applications

SK1 is an anti-human monoclonal antibody that is specific for the CD8a antigen. CD8a (sometimes called MAL or Leu2) is a 32 - 34 kD member of the Ig superfamily that is expressed on the surface of cells such as NK cells and T cells. CD8 is a component of important cellular pathways, for example, 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 important macromolecules/ligands like MHC1. CD8 is a very popular antibody target, with over 125000 publications in the last decade. CD8a is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PerCP/Cy5.5 (ex/em = 489/679 nm). It is compatible with the 488 nm laser and 693/37 nm bandpass filter (for example, as in the Miltenyi

Biotec MACSQuant Analyzer 10).