

PacBlue Anti-human CD8 Antibody *SK1*Catalog number: 100811J0, 100811J1
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
Immunogen	CD8a (T8, Leu2)
Clone	SK1
Conjugate	PacBlue

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 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 MHCI. 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/45 nm bandpass filter (for example, as in the Beckman Coulter DxFLEx).