

iFluor™ A7 Anti-human CD8 Antibody *SK1*

Catalog number: 100810S0, 100810S1

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	iFluor™ A7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ A7
-----------	------------

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 iFluor™ A7.