

iFluor™ 840 Anti-human CD8 Antibody
OKT-8Catalog number: 100820Q0, 100820Q1
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	CD8a (T8, Leu2)
Clone	OKT-8
Conjugate	iFluor™ 840

Biological Properties

Appearance	Dark brown liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 840 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 840
Excitation Wavelength	836 nm
Emission Wavelength	879 nm

Applications

OKT-8 is an anti-human monoclonal antibody that forms an immune complex with the CD8a antigen. CD8a (alternatively called T8 or MAL) is a 32 - 34 kD transmembrane glycoprotein that is located on the surface of cells such as T cells. CD8 is a component of essential 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 important macromolecules/ligands like Lck and MHCI. CD8 is a very popular antibody target, with over 120000 publications in the last decade. CD8a has a variety of applications in immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 840 (ex/em = 836/879 nm).