

**mFluor™ Green 620 Anti-human CD8  
Antibody \*OKT-8\***Catalog number: 100820U0, 100820U1  
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	mFluor™ Green 620

**Biological Properties**

---

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACSymphony™ A5).