

**PE/XFD610 Anti-human CD8 Antibody**  
**\*OKT-8, XFD610 Same Structure to Alexa**  
**Fluor™ 610\***

Catalog number: 10082100, 10082101, 10082102  
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
Isotype	Mouse igg2a
Immunogen	CD8a (T8, Leu2)
Clone	OKT-8
Conjugate	PE/AF610

### Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)

### Spectral Properties

Conjugate	PE/AF610
Excitation Wavelength	567 nm
Emission Wavelength	627 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 PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 610 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 620/15 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).