

XFD350 Anti-human CD8 Antibody *OKT-8*

Catalog Number: 10082140, 10082141

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	Lot specific (please consult certificate of analysis for given lot)
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	AF350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF350
Excitation Wavelength	343 nm
Emission Wavelength	441 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 XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).