

XFD488 Anti-human CD1 Antibody *OKT-6*

Catalog Number: 10011150, 10011151

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 igg1
Immunogen	CD1a (R4, T6)
Clone	OKT-6
Conjugate	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 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	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 nm

Applications

OKT-6 is an anti-human monoclonal antibody that recognizes the CD1a antigen. CD1a (alternatively called T6 or R4) is a 49 kD member of the Ig superfamily that is found on the surface of cells like T cells, B cells, dendritic cells and macrophages. CD1 has been thought to be involved with key biological processes such as immune response, particularly adaptive immune response. Additionally, in some organisms, it acts to positively regulate T cell mediated cytotoxicity. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as β -2-Microglobulin and CD74. CD1 is a moderately popular antibody target, with over 15000 publications in the last decade. CD1a is vital to innate immunity and immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 533/30 nm bandpass filter (for example, as in the BD Accuri™ C6 Plus).