

iFluor™ 750 Anti-human CD1 Antibody
OKT-6Catalog number: 100110L0, 100110L1
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 igg1
Immunogen	CD1a (R4, T6)
Clone	OKT-6
Conjugate	iFluor™ 750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 750
Excitation Wavelength	757 nm
Emission Wavelength	779 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 iFluor™ 750 (ex/em = 757/779 nm).