

iFluor™ A7 Anti-human CD1 Antibody
L161Catalog number: 100130S0, 100130S1
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 kappa
Immunogen	CD1c (R7, M241)
Clone	L161
Conjugate	iFluor™ A7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ A7
-----------	------------

Applications

L161 is an anti-human monoclonal antibody that is specific for the CD1c antigen. CD1c (also known as M241 or R7) is a 43 kD glycoprotein that is expressed on the surface of cells such as dendritic cells, macrophages, B cells and T cells. In certain organisms, CD1 acts to positively regulate T cell mediated cytotoxicity, and is associated with a variety of biologically interesting macromolecules/ligands, for example, β -2-microglobulin. CD1 is a moderately popular antibody target, with over 15000 publications in the last decade. CD1c is essential for immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ A7.