

Purified Anti-human CD1 Antibody *L161*

Catalog number: 10013000

Unit size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD1c (R7, M241)
Clone	L161

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.