

PacBlue Anti-human CD1 Antibody *L161*Catalog number: 100131I0, 100131I1
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	PacBlue

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 nm

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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSCelesta™).