

PacBlue Anti-human CD57 Antibody
HI57aCatalog number: 105701J0, 105701J1
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 IgM
Immunogen	CD57 (HNK-1, Leu-7)
Clone	HI57a
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

The HI57a monoclonal antibody binds with human CD57, a 100 - 115 kD transmembrane protein often expressed on the surface of T cells, natural killer cells, granulocytes and platelets. CD57 is associated with a variety of biologically interesting macromolecules/ligands, for example, P-Selectin and Laminin. CD57 is a fairly uncommon antibody target, with a little more than 3000 publications in the last decade. Even still, CD57

has a variety of applications in costimulatory molecules 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 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 FACSJazz™).