

PE/iFluor™ 647 Anti-human CD57 Antibody
HI57aCatalog number: 105701Q0, 105701Q1, 105701Q2
Unit size: 25 tests, 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	PE/iFluor™ 647

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 647
Excitation Wavelength	569 nm
Emission Wavelength	666 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 PE/iFluor™ 647 (ex/em

= 569/666 nm). It is compatible with the 561 nm laser and 670/30 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).