

APC Anti-human CD57 Antibody *AHC0061*Catalog number: 105711C0, 105711C1, 105711C2
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	IgM, kappa light chain
Immunogen	CD57 (HNK-1, Leu-7)
Clone	AHC0061
Conjugate	APC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC
Excitation Wavelength	651 nm
Emission Wavelength	660 nm

Applications

AHC0061 is an anti-human monoclonal antibody that is specific for the CD57 antigen. CD57 (sometimes referred to as LEU7, NK-1 or HNK-1) is a 100 - 115 kD glycoprotein that is expressed on the surface of cells such as T cells and NK cells. CD57 is associated with a variety of biologically interesting macromolecules/ligands, for example, laminin, L-Selectin and P-Selectin. CD57 is a fairly uncommon antibody target, with a little more than 3800 publications in the last decade. Even still, CD57 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to APC (ex/em = 651/660 nm). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).