

**APC Anti-human CD143 Antibody \*5-369\***Catalog number: 114301B0, 114301B1, 114301B2  
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 IgG1
Immunogen	CD143 (ACE, kininase II)
Clone	5-369
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**

---

The 5-369 monoclonal antibody reacts with human CD143, a 170 kD transmembrane protein typically expressed on the surface of dendritic cells, epithelial cells, endothelial cells and macrophages. CD143 is associated with a variety of biologically interesting macromolecules/ligands, for instance, ANG-1. CD143 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD143 is vital to immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).