

PE/iFluor™ 647 Anti-human CD143 Antibody
5-369Catalog number: 114301P0, 114301P1, 114301P2
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	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 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 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 BD FACSJazz™).