

**FITC Anti-human CD143 Antibody \*5-369\***Catalog number: 114301G0, 114301G1  
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 IgG1               |
| Immunogen          | CD143 (ACE, kininase II) |
| Clone              | 5-369                    |
| Conjugate          | FITC                     |

**Biological Properties**

|             |                                                                                                     |
|-------------|-----------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                         |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | FITC   |
| Excitation Wavelength | 491 nm |
| Emission Wavelength   | 516 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant X).