

**PE Anti-human CD43 Antibody \*HI165\***Catalog number: 104301L0, 104301L1, 104301L2  
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	CD43 (Leukocyte Sialoglycoprotein, Leukosialin, Galactoglycoprotein, SPN)
Clone	HI165
Conjugate	PE

**Biological Properties**

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

**Spectral Properties**

Conjugate	PE
Excitation Wavelength	566 nm
Emission Wavelength	574 nm

**Applications**

The HI165 monoclonal antibody reacts with human CD43, a 95 - 135 kD transmembrane protein commonly located on the surface of plasma cells, thymocytes, neutrophils, myelomas and T cells. In some organisms, CD43 plays a role in the downregulation of cell adhesion, promotes tumor necrosis factor biosynthetic process and acts to negatively regulate T cell proliferation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as EZR. CD43 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD43 has been widely used in 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 (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).