

**APC/iFluor™ 700 Anti-human CD45  
Antibody \*HI30\***Catalog number: 104501F0, 104501F1, 104501F2  
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	CD45 (Leukocyte Common Antigen (LCA), T200, PTPRC)
Clone	HI30
Conjugate	APC/iFluor™ 700

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 nm

**Applications**

---

The HI30 monoclonal antibody recognizes human CD45, a 180 - 240 kD transmembrane glycoprotein typically found on the surface of B cells, neutrophils, hematopoietic cells and dendritic cells. CD45 is a component of vital cellular pathways, in particular, the regulation of receptor signaling pathway via JAK-STAT, B cell receptor signaling pathway and T cell receptor signaling pathway. Also, in certain organisms, it plays a role in the downregulation of cytokine-mediated signaling pathway, represses interleukin-2 biosynthetic process and is an enhancer of humoral immune response mediated by circulating immunoglobulin. CD45 has been thought to be involved with critical biological processes such as

dephosphorylation, especially protein dephosphorylation, and is associated with a variety of biologically interesting macromolecules/ligands, for instance, p56lck. CD45 is a very popular antibody target, with over 50000 publications in the last decade. CD45 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of neuroscience, immunology and neuroscience cell markers. This antibody was purified through affinity chromatography and conjugated to APC/iFluor™ 700 (ex/em = 685/710 nm).