

Purified Anti-human CD45 Antibody
HI151Catalog number: 10452000
Unit size: 100 ug**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD45 (Leukocyte Common Antigen (LCA), T200, PTPRC)
Clone	HI151

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

Applications

HI151 is an anti-human monoclonal antibody that targets the CD45 antigen. CD45 (also known as Leukocyte Common Antigen (LCA) or PTPRC) is a 180 - 240 kD transmembrane protein that is expressed on the surface of cells like B cells, macrophages, granulocytes and NK cells. CD45 has been associated with essential biological processes like dephosphorylation, specifically protein dephosphorylation. Also, in some organisms, it plays a role in the upregulation of stem cell proliferation, is an inhibitor of cytokine-mediated signaling pathway and enhances hematopoietic stem cell migration. CD45 acts in vital cellular pathways, for example, the B cell receptor signaling pathway, positive regulation of antigen receptor-mediated signaling pathway and negative regulation of cytokine-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like Src kinases. CD45 is a very popular antibody target, with over 50000 publications in the last decade. CD45 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of cell biology and neuroscience. This antibody was purified through affinity chromatography.