

Purified Anti-human CD20 Antibody *HI47*Catalog number: 10200000
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 IgG3
Immunogen	CD20 (Bp35, B1)
Clone	HI47

Biological Properties

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

Applications

HI47 is an anti-human monoclonal antibody that recognizes the CD20 antigen. CD20 (alternatively called Leukocyte surface antigen Leu-16) is a 33 - 37 kD transmembrane protein that is located on the surface of cells such as B cells and T cells. CD20 is a member of critical cellular pathways, in particular, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands like lyn. CD20 is a very popular antibody target, with over 28000 publications in the last decade. CD20 has been widely used in immunology and costimulatory molecules research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography.