

PerCP Anti-human CD19 Antibody *HIB19*

Catalog number: 101921T0, 101921T1, 101921T2

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
Immunogen	CD19 (B4)
Clone	HIB19
Conjugate	PerCP

Biological Properties

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

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 nm

Applications

HIB19 is an anti-human monoclonal antibody that recognizes the CD19 antigen. CD19 (also known as CVID3) is a 95 kD glycoprotein that is located on the surface of cells like stem cells, dendritic cells and B cells. In certain organisms, CD19 promotes release of sequestered calcium ion into cytosol, plays a role in the upregulation of phosphatidylinositol 3-kinase activity and acts to positively regulate protein kinase B signaling. In addition, it is a member of critical cellular pathways, for example, the antigen receptor-mediated signaling pathway and B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as lyn. CD19 is a very popular antibody target, with over 36000 publications in the last decade. CD19 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).