

PerCP/Cy5.5 Anti-human CD42b Antibody
HIP1Catalog number: 104201U0, 104201U1, 104201U2
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	CD42b (gplb α , Platelet glycoprotein Ib α)
Clone	HIP1
Conjugate	PerCP/Cy5.5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP/Cy5.5
Excitation Wavelength	489 nm
Emission Wavelength	679 nm

Applications

HIP1 is an anti-human monoclonal antibody that targets the CD42b antigen. CD42b (sometimes called GP1BA or Glycocalicin) is a 145 kD single-pass type I membrane protein that is found on the surface of cells such as stem cells and platelets. CD42b is involved with important cellular pathways, for instance, the cell surface receptor signaling pathway and blood coagulation, intrinsic pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as von Willebrand factor (vWF), Thrombin, c and d. CD42b is a relatively rare antibody target, with fewer than 800 publications in the last decade. Even still, CD42b is vital to cell adhesion and cell biology

research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PerCP/Cy5.5 (ex/em = 489/679 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACSCelesta™).