

**iFluor™ 647 Anti-human CD42b Antibody**  
**\*HIP1\***Catalog number: 104200F0, 104200F1  
Unit size: 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 $\alpha$ , Platelet glycoprotein Ib $\alpha$ )
Clone	HIP1
Conjugate	iFluor™ 647

**Biological Properties**

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 647
Excitation Wavelength	656 nm
Emission Wavelength	670 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 iFluor™ 647 (ex/em = 656/670 nm). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).