

**PE/Cy5 Anti-human CD42b Antibody \*HIP1\***Catalog number: 104201M0, 104201M1, 104201M2  
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 $\alpha$ , Platelet glycoprotein Ib $\alpha$ ) |
| Clone              | HIP1                                                       |
| Conjugate          | PE/Cy5                                                     |

**Biological Properties**

|             |                                                                                                       |
|-------------|-------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                 |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PE/Cy5 |
| Excitation Wavelength | 565 nm |
| Emission Wavelength   | 666 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 PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 661/20 nm

bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).