

PacBlue Anti-human CD41 Antibody *HIP2*Catalog number: 104111J0, 104111J1
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 IgG3
Immunogen	CD41 (GPIIb, ITGA2B)
Clone	HIP2
Conjugate	PacBlue

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 nm

Applications

The HIP2 monoclonal antibody binds to human CD41, a 22 kD glycoprotein frequently expressed on the surface of megakaryocytes and platelets. In certain organisms, CD41 is a promoter of leukocyte migration. Also, it acts in important cellular pathways, for example, the integrin-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands such as von Willebrand factor (vWF), Fibrinogen and Fibronectin. CD41 is a fairly uncommon antibody target, with a little more than 4000

publications in the last decade. Even still, CD41 is vital to cell adhesion, immunology and cell biology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).