

**PacOrange Anti-human CD41 Antibody**  
**\*HIP8\***Catalog number: 104101K0, 104101K1  
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	CD41 (GPIIb, ITGA2B)
Clone	HIP8
Conjugate	PacOrange

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	PacOrange
Excitation Wavelength	400 nm
Emission Wavelength	551 nm

**Applications**

HIP8 is an anti-human monoclonal antibody that recognizes the CD41 antigen. CD41 (also known as GP2B) is a 22 kD member of the Integrin family that is found on the surface of cells such as stem cells. CD41 is a component of critical cellular pathways, for instance, the integrin-mediated signaling pathway. Moreover, in some organisms, it enhances leukocyte migration. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as von Willebrand factor (vWF) and Fibrinogen. CD41 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD41 is frequently used in flow cytometry applications as

a phenotypic marker for differentiation of cell types, specifically in the study of cell biology and cell adhesion. This antibody was purified through affinity chromatography and conjugated to PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).