

FITC Anti-human CD41 Antibody *HIP8*Catalog number: 104101H0, 104101H1
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	FITC

Biological Properties

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

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).