

iFluor™ 430 Anti-human CD41 Antibody
HIP2Catalog number: 10411030, 10411031
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	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 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 iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).