

PE/XFD610 Anti-human CD41 Antibody
***HIP2, XFD610 Same Structure to Alexa**
Fluor™ 610*Catalog number: 10411100, 10411101, 10411102
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 IgG3
Immunogen	CD41 (GPIIb, ITGA2B)
Clone	HIP2
Conjugate	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/AF610
Excitation Wavelength	567 nm
Emission Wavelength	627 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 PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 610 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).