

PE/XFD700 Anti-human CD41 Antibody *HIP2*

Catalog Number: 104111P0,
104111P1, 104111P2
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	Lot specific (please consult certificate of analysis for given lot)
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/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 697/58 nm bandpass filter (for example, as in the BD FACSMelody™).