

APC/XFD750 Anti-human CD41 Antibody

HIP8, XFD750 Same Structure to Alexa Fluor™ 750

Catalog number: 104101E0, 104101E1, 104101E2
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 IgG1
Immunogen	CD41 (GPIIb, ITGA2B)
Clone	HIP8
Conjugate	APC/AF750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/AF750
Excitation Wavelength	756 nm
Emission Wavelength	785 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 APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher).