

APC/XFD750 Anti-human CD235 Antibody *HIR2*

Catalog Number: 123511E0,
123511E1, 123511E2
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 IgG2b
Immunogen	CD235a (GYPA, Sialoglycoprotein alpha, PAS-2)
Clone	HIR2
Conjugate	APC/AF750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 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	APC/AF750
-----------	-----------

Excitation Wavelength 651 nm

Emission Wavelength 785 nm

Applications

HIR2 is an anti-human monoclonal antibody that targets the CD235a antigen. CD235a (also known as Glycophorin A) is a 10 kD single-pass type I membrane protein that is found on the surface of cells such as stem cells and erythrocytes. CD235 is associated with a variety of biologically interesting macromolecules/ligands, in particular, Plasmodium falciparum erythrocyte binding antigen EBA-175. CD235 is a relatively rare antibody target, with less than 10 publications in the last decade. Even still, CD235a has been widely used in cell adhesion and immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC/AF750 (ex/em = 660/785 nm). AF750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).