

XFD647 Anti-human CD235 Antibody *HI264*

Catalog Number: 12350180, 12350181

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	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 IgG2a
Immunogen	CD235a (GYPA, Sialoglycoprotein alpha, PAS-2)
Clone	HI264
Conjugate	AF647

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
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	AF647
Excitation Wavelength	650 nm
Emission Wavelength	671 nm

Applications

HI264 is an anti-human monoclonal antibody that forms an immune complex with the CD235a antigen. CD235a (sometimes called MN sialoglycoprotein or MNS blood group) is a 10 kD single-pass type I membrane protein that is expressed on the surface of cells like erythrocytes. CD235 is associated with a variety of biologically interesting macromolecules/ligands, namely, Plasmodium falciparum erythrocyte binding antigen EBA-175, CD170 and Influenza virus. CD235 is a relatively rare antibody target, with fewer than 6000 publications in the last decade. Even still, CD235a is essential for 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 XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).