

**mFluor™ Violet 540 Anti-human CD235
Antibody *HI264***Catalog number: 12350120, 12350121
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	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 IgG2a
Immunogen	CD235a (GYPA, Sialoglycoprotein alpha, PAS-2)
Clone	HI264
Conjugate	mFluor™ Violet 540

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 540 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 540
Excitation Wavelength	394 nm
Emission Wavelength	537 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 FAKE6000FAKE 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 mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the BD FACSCanto™).