

**mFluor™ Red 780 Anti-human CD235
Antibody *HIR2***Catalog number: 123510W0, 123510W1
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 IgG2b
Immunogen	CD235a (GYPA, Sialoglycoprotein alpha, PAS-2)
Clone	HIR2
Conjugate	mFluor™ Red 780

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 780 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Red 780
Excitation Wavelength	629 nm
Emission Wavelength	767 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 mFluor™ Red 780 (ex/em = 629/767 nm). It is compatible with the 633 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSCanto™ II).