

iFluor™ 820 Anti-human CD235 Antibody
HIR2Catalog number: 123510P0, 123510P1
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	iFluor™ 820

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 820 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 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 iFluor™ 820 (ex/em = 822/850 nm).