

APC/Cy7 Anti-human CD319 Antibody
162.1Catalog number: 131901D0, 131901D1, 131901D2
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	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	CD319 (SLAMF7, CRACC)
Clone	162.1
Conjugate	APC/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/Cy7
Excitation Wavelength	754 nm
Emission Wavelength	779 nm

Applications

The 162.1 monoclonal antibody binds to human CD319, a 66 kD single-pass type I membrane protein often found on the surface of B cells, dendritic cells, lymphocytes and tregs. CD319 is associated with a variety of biologically interesting macromolecules/ligands, for example, CD319. CD319 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD319 has been widely used in 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/Cy7 (ex/em = 754/779 nm).