

**PerCP Anti-non-human primates/ human
CD157 Antibody *SY11B5***Catalog number: 115701S0, 115701S1, 115701S2
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD157 (BST-1)
Clone	SY11B5
Conjugate	PerCP

Biological Properties

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

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 nm

Applications

The SY11B5 monoclonal antibody binds with non-human primates/ human CD157, a 40 - 46 kD transmembrane glycoprotein frequently found on the surface of spleens, thymus, endothelial cells and lungs. CD157 is associated with a variety of biologically interesting macromolecules/ligands, for example, NAD and cyclic ADP-ribose. CD157 is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD157 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of innate immunity. This antibody was purified through affinity chromatography and conjugated to PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 692/40 nm bandpass filter (for example, as in the BD FACSJazz™).