

**mFluor™ Violet 450 Anti-non-human
primates/ human CD157 Antibody
*SY11B5***Catalog number: 115700Z0, 115700Z1
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD157 (BST-1)
Clone	SY11B5
Conjugate	mFluor™ Violet 450

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 452/45 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).