

**iFluor™ 680 Anti-non-human primates/
human CD157 Antibody *SY11B5***Catalog number: 11570010, 11570011
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	iFluor™ 680

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 680 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 680
Excitation Wavelength	684 nm
Emission Wavelength	701 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 iFluor™ 680 (ex/em = 684/701 nm).