

**APC/iFluor™ 750 Anti-human CD229  
Antibody \*HLy9.25\***Catalog number: 122901G0, 122901G1, 122901G2  
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 IgG1
Immunogen	CD229 (Ly9)
Clone	HLy9.25
Conjugate	APC/iFluor™ 750

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	APC/iFluor™ 750
Excitation Wavelength	754 nm
Emission Wavelength	793 nm

**Applications**

---

The HLy9.25 monoclonal antibody reacts with human CD229, a 100 - 120 kD transmembrane protein frequently found on the surface of thymocytes, B cells, T cells and natural killer cells. In some organisms, CD229 positively regulates interleukin-17 production, and is associated with a variety of biologically interesting macromolecules/ligands, namely, SAP. CD229 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD229 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to APC/iFluor™

750 (ex/em = 754/793 nm).