

iFluor™ 568 Anti-human CD229 Antibody
HLy9.25Catalog number: 122900B0, 122900B1
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD229 (Ly9)
Clone	HLy9.25
Conjugate	iFluor™ 568

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 568
Excitation Wavelength	568 nm
Emission Wavelength	587 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 iFluor™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).