

XFD555 Anti-human CD229 Antibody *HLy9.25*

Catalog Number: 12290160, 12290161

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	Lot specific (please consult certificate of analysis for given lot)
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	AF555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate AF555

Excitation Wavelength 553 nm

Emission Wavelength 568 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 XFD555 (ex/em = 553/568 nm). XFD555 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 555 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 583/24 nm bandpass filter (for example, as in the Luminex Amnis CellStream).