

XFD350 Anti-human CD279 Antibody *J116*

Catalog Number: 12793140, 12793141

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, κ
Immunogen	CD279 (PD1)
Clone	J116
Conjugate	AF350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L 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 AF350

Excitation Wavelength 343 nm

Emission Wavelength 441 nm

Applications

The J116 monoclonal antibody binds to human CD279, a 50 - 55 kD member of the Ig superfamily commonly located on the surface of B cells and thymocytes. CD279 is associated with a variety of biologically interesting macromolecules/ligands, for instance, PDL1. CD279 is a relatively rare antibody target, with fewer than 1000 publications in the last decade. Even still, CD279 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of inhibitory molecules, cancer biomarkers and immunology. This antibody was purified through affinity chromatography and conjugated to XFD350 (ex/em = 343/441 nm). XFD350 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 350 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).