

PE/XFD610 Anti-human CD279 Antibody
***J116, XFD610 Same Structure to Alexa**
Fluor™ 610*Catalog number: 12793100, 12793101, 12793102
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, κ
Immunogen	CD279 (PD1)
Clone	J116
Conjugate	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/AF610
Excitation Wavelength	567 nm
Emission Wavelength	627 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 PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT

Bioquest, and it has the same chemical structure of Alexa Fluor® 610 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 561 nm laser and 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).