

**APC/XFD750 Anti-human CD279 Antibody**  
**\*J116, XFD750 Same Structure to Alexa**  
**Fluor™ 750\***Catalog number: 127931E0, 127931E1, 127931E2  
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	APC/AF750

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	APC/AF750
Excitation Wavelength	756 nm
Emission Wavelength	785 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 APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher).