

## XFD750 Anti-human CD279 Antibody \*J116\*

Catalog Number: 127931B0,

127931B1

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, $\kappa$ |
| Immunogen          | CD279 (PD1)          |
| Clone              | J116                 |
| Conjugate          | AF750                |

### Biological Properties

---

|                       |  |
|-----------------------|--|
| Appearance            | Dark blue liquid   |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions   |
| Application           | Flow Cytometry (FACS), Fluorescence Imaging  |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 $\mu$ L/million cells in 100 $\mu$ 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 | AF750 |
|-----------|-------|

Excitation Wavelength 752 nm

Emission Wavelength 776 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 XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).