

**Purified Anti-human CD279 Antibody  
\*J110\***Catalog number: 12792000  
Unit size: 100 ug**Product Details**

---

|                    |   |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.         |
| Expiration Date    | 12 months upon receiving                                    |
| Concentration      | 0.5 mg/mL   |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide |

**Antibody Properties**

---

|                    |               |
|--------------------|---------------|
| Species Reactivity | Human         |
| Class              | Primary       |
| Clonality          | Monoclonal    |
| Host               | Mouse         |
| Isotype            | Mouse igg1, κ |
| Immunogen          | CD279 (PD1)   |
| Clone              | J110          |

**Biological Properties**

---

|             |  |
|-------------|--|
| Appearance  | Liquid   |
| Preparation | Antibody purified by affinity chromatography   |
| Application | Flow Cytometry (FACS), ELISA, HC, Western Blot |

**Applications**

---

J110 is an anti-human monoclonal antibody that targets the CD279 antigen. CD279 (sometimes called Programmed Death-1 or PD-1) is a 50 - 55 kD member of the Ig superfamily that is found on the surface of cells such as T cells and B cells. CD279 is associated with a variety of biologically interesting macromolecules/ligands, in particular, PDL1. CD279 is a relatively rare antibody target, with fewer than 1000 publications in the last decade. Even still, CD279 is vital to cancer biomarkers and immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography.