

**PerCP Anti-human CD279 Antibody \*3D1\***Catalog number: 127901S0, 127901S1, 127901S2  
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       |
| Immunogen          | CD279 (PD1) |
| Clone              | 3D1         |
| Conjugate          | PerCP       |

**Biological Properties**

|             |                                                                                                      |
|-------------|------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PerCP  |
| Excitation Wavelength | 477 nm |
| Emission Wavelength   | 678 nm |

**Applications**

The 3D1 monoclonal antibody reacts with human CD279, a 50 - 55 kD member of the Ig superfamily typically located on the surface of t cells and b cells. CD279 is associated with a variety of biologically interesting macromolecules/ligands, namely, PDL1. CD279 is a relatively rare antibody target, with less than 1000 publications in the last decade. Even still, CD279 is vital to cancer biomarkers research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 675/25 nm bandpass filter (for example, as in the BD Accuri™ C6 Plus).