

**APC/Cy7 Anti-human CD279 Antibody
*2E7***Catalog number: 127911D0, 127911D1, 127911D2
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	2E7
Conjugate	APC/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/Cy7
Excitation Wavelength	754 nm
Emission Wavelength	779 nm

Applications

2E7 is an anti-human monoclonal antibody that targets the CD279 antigen. CD279 (sometimes referred to as Programmed Death-1 or PD-1) is a 50 - 55 kD glycoprotein that is found on the surface of cells like B cells and T 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 has been widely used in cancer biomarkers, immunology and inhibitory molecules research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC/Cy7 (ex/em = 754/779 nm).