

**iFluor™ 560 Anti-human CD279 Antibody**  
**\*3D1\***Catalog number: 127900A0, 127900A1  
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	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	iFluor™ 560

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 560 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 560
Excitation Wavelength	560 nm
Emission Wavelength	571 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 iFluor™ 560 (ex/em = 560/571 nm). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).