

**iFluor™ 710 Anti-human CD279 Antibody
*3D1***Catalog number: 127900K0, 127900K1
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™ 710

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 710
Excitation Wavelength	717 nm
Emission Wavelength	739 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™ 710 (ex/em = 717/739 nm).