

**iFluor™ 594 Anti-human CD279 Antibody
*3D1***Catalog number: 127900C0, 127900C1
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™ 594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 594
Excitation Wavelength	588 nm
Emission Wavelength	604 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™ 594 (ex/em = 588/604 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in

the Luminex Amnis ImageStream).