

iFluor™ 860 Anti-human CD279 Antibody
J110Catalog number: 127920R0, 127920R1
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
Isotype	Mouse igg1, κ
Immunogen	CD279 (PD1)
Clone	J110
Conjugate	iFluor™ 860

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 860 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 860
Excitation Wavelength	853 nm
Emission Wavelength	878 nm

Applications

J110 is an anti-human monoclonal antibody that targets the CD279 antigen. CD279 (sometimes called Programmed Death-1 or PD-1) is a 50 - 55 kD member of the Ig superfamily that is found on the surface of cells such as T cells and B cells. CD279 is associated with a variety of biologically interesting macromolecules/ligands, in particular, PDL1. CD279 is a relatively rare antibody target, with fewer than 1000 publications in the last

decade. Even still, CD279 is vital to cancer biomarkers and immunology 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 iFluor™ 860 (ex/em = 853/878 nm).