

**mFluor™ Red 780 Anti-human CD279
Antibody *J110***Catalog number: 127920W0, 127920W1
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 | mFluor™ Red 780 |

Biological Properties

| | |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Appearance | Dark blue liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 780 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-----------------|
| Conjugate | mFluor™ Red 780 |
| Excitation Wavelength | 629 nm |
| Emission Wavelength | 767 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 mFluor™ Red 780 (ex/em = 629/767 nm). It is compatible with the 633 nm laser and 783/56 nm bandpass filter (for example, as in the BD FACSVerse™).