

**XFD488 Anti-human CD279 Antibody \*2E7,  
XFD488 Same Structure to Alexa Fluor™  
488\***Catalog number: 12791150, 12791151  
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              | 2E7         |
| Conjugate          | AF488       |

**Biological Properties**

|             |                                                                                                      |
|-------------|------------------------------------------------------------------------------------------------------|
| Appearance  | Red liquid                                                                                           |
| Preparation | Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                          |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | AF488  |
| Excitation Wavelength | 499 nm |
| Emission Wavelength   | 520 nm |

**Applications**

2E7 is an anti-human monoclonal antibody that targets the CD279 antigen. CD279 (sometimes referred to as Programmed Death-1 or PD-1) is a 50 - 55 kD glycoprotein that is found on the surface of cells like B cells and T 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 has been widely used in cancer biomarkers, immunology and inhibitory molecules 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 488 (Alexa Fluor® is the trademark of ThermoFisher). It is compatible with the 488 nm laser and 530/40 nm bandpass filter (for example, as in the BD FACSJazz™).