

**PE/AF700 Anti-human CD112 Antibody**  
**\*R2.525\***Catalog number: 11120100, 11120101, 11120102  
Unit size: 25 tests, 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 kappa
Immunogen	CD112 (HVEB, Nectin-2, PVRL2, PRR2)
Clone	R2.525
Conjugate	PE/AF700

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

Conjugate	PE/AF700
Excitation Wavelength	566 nm
Emission Wavelength	721 nm

**Applications**

R2.525 is an anti-human monoclonal antibody that forms an immune complex with the CD112 antigen. CD112 (also known as PVRL2, Nectin-2 or HVEB) is a glycoprotein that is located on the surface of cells such as macrophages and epithelial cells. CD112 is associated with a variety of biologically interesting macromolecules/ligands, in particular, PRR3, CD112 and afadin. CD112 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD112 is essential for immunology 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 PE/AF700 (ex/em = 566/721 nm). It is compatible with the 561 nm laser and 720/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell

Analyzer).