

**XFD594 Anti-human CD112 Antibody \*R2.525\***

Catalog Number: 11120160, 11120161

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      | Lot specific (please consult certificate of analysis for given lot)         |
| 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          | AF594                               |

**Biological Properties**

---

|                       |                                                                                                                                                                                                                                                                                                                                                                                                                                             |
|-----------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Appearance            | Purple liquid                                                                                                                                                                                                                                                                                                                                                                                                                               |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions                                                                                                                                                                                                                                                                                                                                        |
| Application           | Flow Cytometry (FACS), Fluorescence Imaging                                                                                                                                                                                                                                                                                                                                                                                                 |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.<br><i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i> |

**Spectral Properties**

---

Conjugate                    AF594

Excitation Wavelength   590 nm

Emission Wavelength    618 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).