

XFD488 Anti-human CD112 Antibody *R2.525*

Catalog Number: 11120140, 11120141

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	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 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. <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 AF488

Excitation Wavelength 499 nm

Emission Wavelength 520 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant Analyzer 16).