

## PE/XFD610 Anti-human CD206 Antibody \*15-2\*

Catalog Number: 120601N0,  
120601N1, 120601N2  
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	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	CD206 (MMR, MRC1)
Clone	15-2
Conjugate	PE/AF610

### Biological Properties

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)
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	PE/AF610
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 627 nm

## Applications

---

The 15-2 monoclonal antibody recognizes human CD206, a 162 - 175 kD transmembrane protein frequently expressed on the surface of endothelial cells, dendritic cells and macrophages. CD206 is associated with a variety of biologically interesting macromolecules/ligands, namely, plant glycoproteins, viral glycoproteins and lutropin. CD206 is a fairly uncommon antibody target, with a little more than 6900 publications in the last decade. Even still, CD206 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of neuroscience cell markers. This antibody was purified through affinity chromatography and conjugated to PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACSCelesta™).