

PE/iFluor™ 700 Anti-human CD206 Antibody
15-2Catalog number: 120601U0, 120601U1, 120601U2
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	CD206 (MMR, MRC1)
Clone	15-2
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 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/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).