

PacOrange Anti-human CD206 Antibody
15-2Catalog number: 120601J0, 120601J1
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	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	PacOrange

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacOrange
Excitation Wavelength	400 nm
Emission Wavelength	551 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

PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm laser and 528/46 nm bandpass filter (for example, as in the Luminex Amnis CellStream).