

APC/Cy7 Anti-mouse CD197 Antibody
4B12Catalog number: 119701C0, 119701C1, 119701C2
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD197 (CCR7, BLR2, EBI1, CMKBR7)
Clone	4B12
Conjugate	APC/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/Cy7
Excitation Wavelength	754 nm
Emission Wavelength	779 nm

Applications

4B12 is an anti-mouse monoclonal antibody that is specific for the CD197 antigen. CD197 (sometimes called BLR2, CCR7, EBI1 or CMKBR7) is a glycoprotein that is expressed on the surface of cells like B cells, T cells and dendritic cells. CD197 is associated with a variety of biologically interesting macromolecules/ligands, for example, CCL19/ECL and CCL21. CD197 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD197 is essential for immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC/Cy7 (ex/em = 754/779 nm).