

PerCP Anti-mouse CD197 Antibody *4B12*Catalog number: 119701S0, 119701S1, 119701S2
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	PerCP

Biological Properties

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

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 660/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).