

PE/Cy5 Anti-human CD172 Antibody *B4B6*Catalog number: 117211L0, 117211L1, 117211L2
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
Immunogen	CD172b (SIRPb)
Clone	B4B6
Conjugate	PE/Cy5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy5
Excitation Wavelength	565 nm
Emission Wavelength	666 nm

Applications

The B4B6 monoclonal antibody binds to human CD172b, a 43 kD transmembrane protein often located on the surface of dendritic cells, myeloid cells and monocytes. CD172 is associated with a variety of biologically interesting macromolecules/ligands, in particular, TYROBP. CD172 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD172b has a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 661/20 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).