

PE/Cy7 Anti-human CD172a Antibody
15-414Catalog number: 117201M0, 117201M1, 117201M2
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 IgG2a
Immunogen	CD172a (SHPS-1, SIRP α , BIT, MFR, MYD-1, P84)
Clone	15-414
Conjugate	PE/Cy7

Biological Properties

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

Spectral Properties

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

The 15-414 monoclonal antibody binds with human CD172a, a 85 - 90 kD glycoprotein often found on the surface of dendritic cells and neurons. CD172a is associated with a variety of biologically interesting macromolecules/ligands, in particular, CD47 and PTPN11. CD172a is a relatively rare antibody target, with fewer than 700 publications in the last decade. Even still, CD172a has a variety of applications in immunology and cell biology 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/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSAria™ III).