

PerCP Anti-human CD172a Antibody
15-414Catalog number: 117201S0, 117201S1, 117201S2
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, SIRPα, BIT, MFR, MYD-1, P84)
Clone	15-414
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

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 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).