

**PE/iFluor™ 700 Anti-human CD140a
Antibody *16A1***Catalog number: 114001U0, 114001U1, 114001U2
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	CD140a (PDGFRa)
Clone	16A1
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 nm

Applications

16A1 is an anti-human monoclonal antibody that is specific for the CD140a antigen. CD140a (sometimes called PDGFRa) is a member of the platelet-derived growth factor family that is expressed on the surface of cells such as endothelial cells and platelets. CD140a is associated with a variety of biologically interesting macromolecules/ligands, for instance, PDGF. CD140a is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD140a is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of neuroscience, stem cells and immunology. This antibody was purified through affinity chromatography and

conjugated to PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 720/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).