

Purified Anti-human CD140a Antibody
16A1Catalog number: 11400000
Unit size: 100 ug**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD140a (PDGFRa)
Clone	16A1

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.