

iFluor™ 710 Anti-human CD140a Antibody
16A1Catalog number: 114000K0, 114000K1
Unit size: 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	iFluor™ 710

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 710
Excitation Wavelength	717 nm
Emission Wavelength	739 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 iFluor™ 710 (ex/em = 717/739 nm).