

**PE/XFD700 Anti-human CD140a Antibody \*16A1\***

Catalog Number: 114001O0,  
114001O1, 114001O2  
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	Lot specific (please consult certificate of analysis for given lot)
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/AF700

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

**Spectral Properties**

---

Conjugate	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 720/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).