

iFluor™ 594 Anti-human CD140b Antibody
18A2Catalog number: 114010C0, 114010C1
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	CD140b (PDGFRb)
Clone	18A2
Conjugate	iFluor™ 594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 594
Excitation Wavelength	588 nm
Emission Wavelength	604 nm

Applications

18A2 is an anti-human monoclonal antibody that recognizes the CD140b antigen. CD140b (also known as PDGFRb) is a transmembrane glycoprotein that is expressed on the surface of cells like endothelial cells. CD140b is associated with a variety of biologically interesting macromolecules/ligands, namely, PDGF. CD140b is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even

still, CD140b is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of neuroscience cell markers. This antibody was purified through affinity chromatography and conjugated to iFluor™ 594 (ex/em = 588/604 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).