

**XFD647 Anti-human CD140b Antibody \*18A2\***

Catalog Number: 11401170, 11401171

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      | 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          | CD140b (PDGFRb) |
| Clone              | 18A2            |
| Conjugate          | AF647           |

**Biological Properties**

---

|                       |                                                                                                                                                                                                                                                                                                                                                                                                                                                 |
|-----------------------|-------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Appearance            | Dark blue liquid                                                                                                                                                                                                                                                                                                                                                                                                                                |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF647 under optimal conditions                                                                                                                                                                                                                                                                                                                                            |
| Application           | Flow Cytometry (FACS), Fluorescence Imaging                                                                                                                                                                                                                                                                                                                                                                                                     |
| 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.<br><br><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 AF647

Excitation Wavelength 650 nm

Emission Wavelength 671 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 XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).