

**XFD555 Anti-human CD147 Antibody \*MEM-M6/6\***

Catalog Number: 11470150, 11470151

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          | CD147 (Basigin, EMMPRIN) |
| Clone              | MEM-M6/6                 |
| Conjugate          | AF555                    |

**Biological Properties**

---

|                       |                                                                                                                                                                                                                                                                                                                                                                                                                                             |
|-----------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Appearance            | Red liquid                                                                                                                                                                                                                                                                                                                                                                                                                                  |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF555 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><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             | AF555  |
| Excitation Wavelength | 553 nm |
| Emission Wavelength   | 568 nm |

## Applications

---

The MEM-M6/6 monoclonal antibody binds to human CD147, a 55 - 65 kD member of the Ig superfamily often expressed on the surface of endothelial cells and leukocytes. In certain organisms, CD147 acts to positively regulate interleukin-6 production. Moreover, it is a component of key cellular pathways, for example, the cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands. CD147 is a fairly uncommon antibody target, with a little more than 3200 publications in the last decade. Even still, CD147 has been widely used in cell adhesion and immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD555 (ex/em = 553/568 nm). XFD555 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 555 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).