

XFD594 Anti-human CD147 Antibody *MEM-M6/6*

Catalog Number: 11470160, 11470161

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	AF594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF594 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. <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	AF594
Excitation Wavelength	590 nm
Emission Wavelength	618 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).