

iFluor™ 405 Anti-human CD147 Antibody
MEM-M6/6Catalog number: 11470020, 11470021
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	CD147 (Basigin, EMMPRIN)
Clone	MEM-M6/6
Conjugate	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 445/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).