

XFD594 Anti-cow/ human CD46 Antibody *MEM-258*

Catalog Number: 10460170, 10460171

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	Cow, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD46 (Membrane Cofactor Protein)
Clone	MEM-258
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

MEM-258 is an anti-cow/ human monoclonal antibody that targets the CD46 antigen. CD46 (alternatively called Membrane Cofactor Protein) is a 45 - 75 kD transmembrane protein that is expressed on the surface of cells like platelets, macrophages, granulocytes and B cells. CD46 is associated with a variety of biologically interesting macromolecules/ligands, namely, CD29, SCR2/3/4 and CD9. CD46 is a fairly uncommon antibody target, with a little more than 3600 publications in the last decade. Even still, CD46 is vital to cell biology, complement and immunology research, frequently 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).