

**iFluor™ 820 Anti-cow/ human CD46
Antibody *MEM-258***Catalog number: 104600P0, 104600P1
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	Cow, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD46 (Membrane Cofactor Protein)
Clone	MEM-258
Conjugate	iFluor™ 820

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 820 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 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 iFluor™ 820 (ex/em = 822/850 nm).