

**iFluor™ 670 Anti-rabbit/ ferret/ cow/
human CD88 Antibody *S5/1***Catalog number: 108800H0, 108800H1
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	Rabbit, ferret, cow, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a kappa
Immunogen	CD88 (C5R1, C5AR)
Clone	S5/1
Conjugate	iFluor™ 670

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 670 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 670
Excitation Wavelength	671 nm
Emission Wavelength	682 nm

Applications

S5/1 is an anti-rabbit/ ferret/ cow/ human monoclonal antibody that forms an immune complex with the CD88 antigen. CD88 (also known as C5R1 or C5AR) is a 43 kD transmembrane glycoprotein that is expressed on the surface of cells like endothelial cells, dendritic cells, granulocytes and macrophages. CD88 is associated with a variety of biologically interesting macromolecules/ligands, for example, anaphylatoxin. CD88 is a relatively rare antibody target, with fewer than 700 publications in the last decade. Even still, CD88 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of innate immunity and cell biology. This antibody was purified

through affinity chromatography and conjugated to iFluor™ 670 (ex/em = 671/682 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).