

**PacBlue Anti-rabbit/ ferret/ cow/ human
CD88 Antibody *S5/1***Catalog number: 108801K0, 108801K1
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	PacBlue

Biological Properties

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

Spectral Properties

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).