

XFD488 Anti-rabbit/ ferret/ cow/ human CD88 Antibody

S5/1

Catalog Number: 10880150, 10880151

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	Rabbit, ferret, cow, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a kappa
Immunogen	CD88 (C5R1, C5AR)
Clone	S5/1
Conjugate	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 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	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).