

APC/XFD750 Anti-human/ non-human primates CD89 Antibody *A59*

Catalog Number: 108901E0,
108901E1, 108901E2
Unit Size: 25 tests, 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	Human, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD89 (FCAR)
Clone	A59
Conjugate	APC/AF750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)
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	APC/AF750
-----------	-----------

Excitation Wavelength 651 nm

Emission Wavelength 785 nm

Applications

The A59 monoclonal antibody binds with human/ non-human primates CD89, a 55 - 100 kD single-pass type I membrane protein often expressed on the surface of macrophages, eosinophils and neutrophils. CD89 acts in critical cellular pathways, in particular, the Fc receptor signaling pathway. Furthermore, in certain organisms, it upregulates neutrophil apoptotic process and is a promoter of oxidative stress-induced cell death. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as IgA2 and IgA1. CD89 is a relatively rare antibody target, with fewer than 500 publications in the last decade. Even still, CD89 is essential for immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).