

iFluor™ 790 Anti-human/ non-human primates CD89 Antibody *A59*Catalog number: 108900M0, 108900M1
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	Human, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD89 (FCAR)
Clone	A59
Conjugate	iFluor™ 790

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 790 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 790
Excitation Wavelength	787 nm
Emission Wavelength	812 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 iFluor™ 790 (ex/em = 787/812 nm).