

**iFluor™ 450 Anti-human/ non-human
primates CD89 Antibody *A59***Catalog number: 10890040, 10890041
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™ 450 |

Biological Properties

| | |
|-------------|--|
| Appearance | Brown liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 450 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 450 |
| Excitation Wavelength | 451 nm |
| Emission Wavelength | 502 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™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSARIA™ III).