

**Purified Mouse Anti-Alexa Fluor 647  
Antibody \*Monoclonal\***Catalog number: V102025  
Unit size: 1.0 mg**Product Details**

---

Storage Conditions	Freeze (<-15°C) with minimized light exposure.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	PBS, pH 7.2 and 0.09% sodium azide

**Antibody Properties**

---

Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Alexa Fluor 647

**Biological Properties**

---

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

**Applications**

---

Alexa Fluor 647 has shown evidence of use in a diverse array of applications, specifically in the field of biotechnology and pharmacological research. Enzyme luciferase, ammonia monooxygenase and target proteases are, for instance, just a few of the enzymes that it has helped scientists to better understand. Alexa Fluor 647 has a variety of applications in the study of vital cellular processes such as endocytosis tracking, dismal prognosis and BRI1-AFCS endocytosis, and it has been widely used in the application of key cell-based pathways like FRET pathway and metabolic pathway.