

**Purified Rabbit Anti-GFP Antibody**  
**\*Polyclonal\***Catalog number: V102110  
Unit size: 0.1 mg**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	, 0.2% (w/v) BSA

**Antibody Properties**

---

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	GFP

**Biological Properties**

---

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

**Applications**

---

Green fluorescent protein is a protein with a molecular weight of 31 kDa. In *aequorea victoria*, green fluorescent protein has been found to be involved in organismal processes, for example, protein-chromophore linkage, generation of precursor metabolites and energy, and bioluminescence.