

**HRP Rabbit Anti-T7 Antibody \*Polyclonal\***

Catalog number: V102180

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	T7
Conjugate	HRP

**Biological Properties**

---

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

**Applications**

---

Identifies the 11 amino acid long peptide derived from T7 bacteriophage gene10.