

## HRP Rabbit Anti-V5 Antibody \*Polyclonal\*

Catalog Number: V102200

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
Formulation	HEPES – saline with 0.002% Thimerosal and stabilizer

### Antibody Properties

---

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	V5

### Biological Properties

---

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with HRP under optimal conditions
Application	FACS

### Spectral Properties

---

Conjugate	HRP
-----------	-----

### Applications

---

Identifies the 14 amino acid long epitope tag, V5, derived from paramyxovirus P/V proteins (amino acid sequence: GKPIPPELLGLDST).