

---

## **VT A1 Polyclonal Antibody, Biotin Conjugated**

(Catalog #A61735)

---

### **Background**

Involved in the endosomal multivesicular bodies (MVB) pathway. MVBs contain intraluminal vesicles (ILVs) that are generated by invagination and scission from the limiting membrane of the endosome and mostly are delivered to lysosomes enabling degradation of membrane proteins, such as stimulated growth factor receptors, lysosomal enzymes and lipids. Thought to be a cofactor of VPS4A/B, which catalyzes disassembles membrane-associated ESCRT-III assemblies. Involved in the sorting and down-regulation of EGFR. Involved in HIV-1 budding.

### **Description**

VT A1 Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### **Formulation**

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### **Specificity**

Human

### **Isotype**

IgG

### **Uniprot ID**

Q9NP79

### **Purification**

>95%, Protein G purified

### **Immunogen**

Recombinant Human Vacuolar protein sorting-associated protein VT A1 homolog protein (23-307AA)

### **Storage**

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### **Alternative Names**

Vacuolar protein sorting-associated protein VT A1 homolog, VT A1, C6orf55, HSPC228, My012, Dopamine-responsive gene 1 protein, DRG-1, LYST-interacting protein 5, LIP5, SKD1-binding protein 1, SBP1

### **Application**

ELISA; Recommended dilution: ELISA 1:500 - 1:1000

---