

## VP35 Polyclonal Antibody

(Catalog # A67074)

---

### Background

Acts as a polymerase cofactor in the RNA polymerase transcription and replication complex. Prevents establishment of cellular antiviral state by blocking virus-induced phosphorylation and activation of interferon regulatory factor 3 (IRF3), a transcription factor critical for the induction of interferons alpha and beta. This blockage is produced through the interaction with and inhibition host IKBKE and TBK1 producing a strong inhibition of the phosphorylation and activation of IRF3. Also inhibits the antiviral effect mediated by the interferon-induced, double-stranded RNA-activated protein kinase EIF2AK2/PKR.

### Description

VP35 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Reston ebolavirus

### Isotype

IgG

### Uniprot ID

Q91DE0

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Reston ebolavirus Polymerase cofactor VP35 protein (1-329AA)

### 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

Polymerase cofactor VP35, VP35

### Application

ELISA

---