

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4521P
<b>Quantity</b>	50 µg
<b>Source</b>	19 amino acids near the carboxy terminus of RVF virus
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
<b>Storage &amp; Stability</b>	Store Rift Valley Fever Virus peptide at -20°C, stable for one year.
<b>Application</b>	Rift Valley Fever Virus peptide is used for blocking the activity of RVF virus antibody.
<b>References</b>	Giorgi C, Accardi L, Nicoletti M, et al. Sequences and coding strategies of the S RNAs of Toscana and Rift Valley fever viruses compared to those of Punta Toro, Sicilian sandfly fever, and Uukuniemi viruses. <i>Virology</i> 1991; 180:738-53.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.