

## Instruction Manual

|                                |   |
|--------------------------------|---|
| <b>Catalog Number</b>          | PK-AB718-3943P  |
| <b>Quantity</b>                | 50 µg   |
| <b>Source</b>                  | 14 amino acids near the center terminus of human IRF7   |
| <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 IRF7 peptide at -20°C, stable for one year.   |
| <b>Application</b>             | IRF7 peptide is used for blocking activity of IRF7 antibody.  |
| <b>References</b>              | Fitzgerald KA, Rowe DC, Barnes BJ, et al. LPS-TLR4 signaling to IRF-3/7 and NF-kappaB involves the toll adaptors TRAM and TRIF. J. Exp. Med. 2003; 198:1043-55.   |

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