

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3799P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the center of human IL-27
<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 IL-27 peptide at -20°C, stable for one year.
<b>Application</b>	IL-27 peptide is used for blocking activity of IL-27 antibody.
<b>References</b>	Pfanz S, Timans JC, Cheung J et al. IL-27, a heterodimeric cytokine composed of EB13 and p28 protein, induces proliferation of naïve CD4(+) T cells. <i>Immunity</i> 2002; 16:779-90

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