

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2471P
<b>Quantity</b>	50 µg
<b>Source</b>	13 amino amino terminus of human IL-21 Receptor
<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-21 Receptor peptide at -20°C, stable for one year.
<b>Application</b>	IL-21 receptor peptide is used for blocking the activity of IL-21 receptor antibody.
<b>References</b>	Arrish-Novak J, Dillon SR, Nelson A, et al. Interleukin 21 and its receptor are involved in NK cell expansion and regulation of lymphocyte function. <i>Nature</i> . 2000;408(6808):57-63. Yu J, Zhang L, Hwang PM, et al. PUMA induces the rapid apoptosis of colorectal cancer cells. <i>Mol. Cell</i> 2001; 7:673-82. Eferl R, Ricci R, Kenner L, et al. Liver tumor development. c-Jun antagonizes the proapoptotic activity of p53. <i>Cell</i> 2003; 112:181-92.

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