

Instruction Manual

Catalog Number	PK-AB718-2469P
Quantity	50 µg
Source	15 amino acids near the extracellular domain of human IL-21 Receptor
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
Reconstitution	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.
Storage & Stability	Store IL-21 Receptor peptide at -20°C, stable for one year.
Application	IL-21 receptor peptide is used for blocking the activity of IL-21 receptor antibody.
References	<p>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.</p> <p>Kelly K, Manos E, Jensen G, et al. APRIL/TRDL-1, a tumor necrosis factor-like ligand, stimulates cell death. <i>Cancer Res</i>. 2000;60(4):1021-7.</p> <p>Yu G, Boone T, Delaney J, et al. APRIL and TALL-1 and receptors BCMA and TACI: system for regulating humoral immunity. <i>Nat Immunol</i>. 2000;1(3):252-6.</p>

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