

Instruction Manual

Catalog Number	PK-AB718-2131P
Quantity	50 µg
Source	21 amino acids near the carboxy terminus of human IL-1RAcP
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-1RAcP peptide at -20°C, stable for one year.
Application	IL-1RAcP peptide is used for blocking the activity of IL-1RAcP antibody.
References	<p>Greenfeder SA, Nunes P, Kwee L, Labow M, Chizzonite RA, Ju G. Molecular cloning and characterization of a second subunit of the interleukin 1 receptor complex. <i>J Biol Chem</i> 1995;270:13757-65.</p> <p>Zandi E, Rothwarf DM, Delhase M, Hayakawa M, Karin M. The IkappaB kinase complex (IKK) contains two kinase subunits, IKKalpha and IKKbeta, necessary for IkappaB phosphorylation and NF-kappaB activation. <i>Cell</i> 1997;91:243-52</p> <p>Woronicz JD, Gao X, Cao Z, Rothe M, Goeddel DY. IkappaB kinase-alpha: NF-kappaB activation and complex formation with IkappaB kinase-alpha and NIK. <i>Science</i> 1997;278:866-9</p>

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