

Instruction Manual

Catalog Number	PK-AB718-2489P
Quantity	50 µg
Source	15 amino acids near the carboxy terminus of human IL-22 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-22 Receptor peptide at -20°C, stable for one year.
Application	IL-22 receptor peptide is used for blocking the activity of IL-22 receptor antibody.
References	<p>Dumoutier L, Louahed J, Renauld JC. Cloning and characterization of IL-10-related T cell-derived inducible factor (IL-TIF), a novel cytokine structurally related to IL-10 and inducible by IL-9. <i>J Immunol.</i> 2000;164(4):1814-9.</p> <p>Oda E, Ohki R, Murasawa H, et al. Noxa, a BH3-only member of the Bcl-2 family and candidate mediator of p53-induced apoptosis. <i>Science.</i> 2000;288(5468):1053-8.</p> <p>Oda K, Arakawa H, Tanaka T, et al. p53AIP1, a potential mediator of p53-dependent apoptosis, and its regulation by Ser-46-phosphorylated p53. <i>Cell.</i> 2000 Sep 15;102(6):849-62.</p>

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