

Instruction Manual

Catalog Number	PK-AB718-3795P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human IL-23
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-23 peptide at -20°C, stable for one year.
Application	IL-23 peptide is used for blocking activity of IL-23 antibody.
References	<p>Oppmann B, Lesley R, Blom B, et al. Novel p19 protein engages IL-12p40 to form a cytokine, IL-23, with biological activities similar as well as distinct from IL-12. <i>Immunity</i> 2000; 13:715-25.</p> <p>Marsters SA, Ayres TM, Skubatch M, et al. Herpesvirus entry mediator, a member of the tumor necrosis factor receptor (TNFR) family, interacts with members of the TNFR-associated factor family and activates the transcription factors NF-kappaB and AP-1. <i>J. Biol. Chem.</i> 1997; 272:14029-32.</p> <p>Mauri DN, Ebner R, Montgomery RI, et al. LIGHT, a new member of the TNF superfamily, and lymphotoxin alpha are ligands for herpesvirus entry mediator. <i>Immunity</i> 1998; 8:21-30.</p>

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