

Instruction Manual

Catalog Number	PK-AB718-3017P
Quantity	50 µg
Source	12 amino amino terminus of human ILP-2
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 ILP-2 peptide at -20°C, stable for one year.
Application	ILP-2 peptide is used for blocking the activity of ILP-2 antibody.
References	<p>Richter BW, Mir SS, Eiben LJ, Lewis J, Reffey SB, Frattini A, Tian L, Frank S, Youle RJ, Nelson DL, Notarangelo LD, Vezzoni P, Fearnhead HO, Duckett CS. Molecular cloning of ILP-2, a novel member of the inhibitor of apoptosis protein family. <i>Mol Cell Biol.</i> 2001;21(13):4292-301.</p> <p>Hugot JP, Chamaillard M, et al. Association of NOD2 leucine-rich repeat variants with susceptibility to Crohn's disease. <i>Nature</i> 2001; 411:599-603.</p> <p>Ogura Y, Bonen DK, Inohara N, et al. A frameshift mutation in NOD2 associated with susceptibility to Crohn's disease. <i>Nature</i> 2001; 411:603-6.</p>

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