

Instruction Manual

Catalog Number	PK-AB718-3199P
Quantity	50 µg
Source	16 amino acids near the center of human ILPIP
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 ILPIP peptide at -20°C, stable for one year.
Application	ILPIP peptide is used for blocking the activity of the ILPIP antibody.
References	<p>Sanna MG, Correia JS, Luo Y, et al. (2002) ILPIP, a novel anti-apoptotic protein that enhances XIAP-mediated activation of JNK and protection of apoptosis. <i>J. Biol. Chem.</i> 277:30454-62.</p> <p>Wajant H and Scheurich P. Analogies between Drosophila and mammalian TRAF pathways. <i>Prog. Mol. Subcell. Biol.</i> 2004; 34:47-72.</p> <p>Kobayashi T, Walsh PT, Walsh MC, et al. TRAF6 is a critical factor for dendritic cell maturation and development. <i>Immunity</i> 2003; 19:353-63.</p> <p>O'Neill LAJ, Fitzgerald FA, and Bowie AG. The Toll-IL-1 receptor adaptor family grows to five members. <i>Trends in Imm.</i> 2003; 24:286-9.</p>

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