

Instruction Manual

Catalog Number	PK-AB718-3377P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human ECSIT
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 ECSIT peptide at -20°C, stable for one year.
Application	ECSIT peptide is used for blocking the activity of the ECSIT antibody.
References	<p>Kopp E, Medzhitov R, Carothers J, et al. ECSIT is an evolutionarily conserved intermediate in the Toll/IL-1 signal transduction pathway. <i>Genes Dev.</i> 1999; 13:2059-71.</p> <p>Vaux DL and Silke J. Mammalian mitochondrial IAP binding proteins. <i>Biochem. Biophys. Res. Comm.</i> 2003; 304:499-504.</p> <p>Suzuki Y, Imai Y, Nakayama H, et al. A serine protease, HtrA2, is released from the mitochondria and interacts with XIAP, inducing cell death. <i>Mol. Cell</i> 2001; 8:613-21.</p> <p>Ahn J-Y, Rong R, Kroll TG, et al. PIKE (Phosphatidylinositol 3-kinase enhancer)-A GTPase stimulates Akt activity and mediates cellular invasion. <i>J. Biol. Chem.</i> 2004; 279:16441-51.</p>

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