

Instruction Manual

Catalog Number	PK-AB718-2213P
Quantity	50 µg
Source	20 amino acids near the carboxy terminus of human IRAK2
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 IRAK2 peptide at -20°C, stable for one year.
Application	IRAK2 peptide is used for blocking the activity of IRAK2 antibody.
References	<p>Muzio M, Ni J, Feng P, Dixit VM. IRAK (Pelle) family member IRAK-2 and MyD88 as proximal mediators of IL-1 signaling. <i>Science</i> 1997;278:1612-5.</p> <p>Sheridan JP, Marsters SA, Pitti RM, et al. A. Control of TRAIL-induced apoptosis by a family of signaling and decoy receptors. <i>Science</i> 1997;277:818-21</p> <p>Degli-Esposti MA, Smolak PJ, Walczak H, et al, Smith CA. Cloning and characterization of TRAIL-R3, a novel member of the emerging TRAIL receptor family. <i>J Exp Med</i> 1997;186(7):1165-70</p>

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