

Instruction Manual

Catalog Number	PK-AB718-1007P
Quantity	50 µg
Source	13 amino acids near the carboxy terminus of human IRAK
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 IRAK peptide at -20°C, stable for one year.
Application	IRAK peptide is used for blocking the activity of IRAK antibody.
References	<p>Cao Z; Henzel WJ; Gao X. IRAK: a kinase associated with the interleukin-1 receptor. <i>Science</i> 1996; 271:1128-31.</p> <p>Jianing Huang, Xiong Gao, Shyun Li, and Zhaodan Cao. Recruitment of IRAK to the interleukin 1 receptor complex requires interleukin-1 receptor accessory protein. <i>Proc Natl Acad Sci USA</i> 1997;94:12829-12832</p> <p>Robinson D, Shibuya K, Mui A, Zonin F, Murphy E, Sana T, Hartley SB, Menon S, Kastelein R, Bazan F, O'Garra A. IGIF does not drive Th1 development but synergizes with IL-12 for interferon-gamma; production and activates IRAK and NF-kappaB. <i>Immunity</i> 1997;7:571-581</p>

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