

Instruction Manual

Catalog Number	PK-AB718-3125P
Quantity	50 µg
Source	Amino acids near the carboxy terminus of human IRAK-4
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 IRAK4 peptide at -20°C, stable for one year.
Application	IRAK4 peptide is used for blocking the activity of IRAK-4 antibody.
References	<p>Li S, Strelow A, Fontana EJ, Wesche H. IRAK-4: a novel member of the IRAK family with the properties of an IRAK-kinase. Proc Natl Acad Sci USA. 2002;99(8):5567-72.</p> <p>Xie J and Guo Q. AATF protects neural cells against oxidative damage induced by amyloid b-peptide. Neurobiol. of Dis. 2004; 16:150-7.</p> <p>Xie J and Guo Q. AATF protects neural cells against oxidative damage induced by amyloid b-peptide. Neurobiol. of Dis. 2004; 16:150-7.</p>

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