

Instruction Manual

Catalog Number	PK-AB718-3091P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human AID
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 AID peptide at -20°C, stable for one year.
Application	AID peptide is used for blocking the activity of AID antibody.
References	<p>Muramatsu M, Sankaranand VS, Anant S, et al. Specific expression of activation-induced cytidine deaminase (AID), a novel member of the RNA-editing deaminase family in germinal center B cells. <i>J. Biol. Chem.</i> 1999; 274:18470-6.</p> <p>Paavola,P., Horelli-Kuitunen,N., Palotie,A. and Peltonen,L. Characterization of a novel gene, PNUTL2, on human chromosome 17q22-q23 and its exclusion as the Meckel syndrome gene. <i>Genomics</i> 1999;55 (1):122-125</p> <p>Zieger,B., Tran,H., Hainmann, I., Wunderle,D., Zgaga-Griesz,A., Blaser,S. and Ware,J. Characterization and expression analysis of two human septin genes, PNUTL1 and PNUTL2. <i>Gene</i> 2000;261(2):197-203</p>

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