

Instruction Manual

Catalog Number	PK-AB718-1117P
Quantity	50 µg
Source	19 amino acids near the carboxy terminus of human RAIDD
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 RAIDD peptide at -20°C, stable for one year.
Application	RAIDD peptide is used for blocking the activity of RAIDD antibody.
References	Duan H, Dixit VM. RAIDD is a new 'death' adaptor molecule. Nature 1997;385:86-89. Deng, H, et al. Nature 1996;381:661-6 Dragic, T, et al. Nature 1996;381:667-73

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