

Instruction Manual

Catalog Number	PK-AB718-4585P
Quantity	50 µg
Source	13 amino acids near the amino terminus of human RGPD5
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 RGPD5 peptide at -20°C, stable for one year.
Application	RGPD5 peptide is used for blocking the activity of RGPD5 antibody.
References	Nothwang HG, Rensing C, Kubler M, et al. Identification of a novel Ran binding protein 2 related gene (RANBP2L1) and detection of a gene cluster on chromosome 2q11-q12. Genomics 1998; 47:383-92.

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