

## Instruction Manual

Catalog Number	PK-AB718-4015P
Quantity	50 µg
Source	19 amino acids near the amino terminus of human RPA interacting protein
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 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 RPA Interacting Protein peptide at -20°C, stable for one year.
Application	RPA interacting protein peptide is used for blocking the activity of RPA interacting protein antibody.
References	Jullien D, Gorlich D, Laemmli UK, et al. Nuclear import of RPA in Xenopus egg extracts requires a novel protein XRIPalpha but not importin alpha . EMBO J. 1999; 18:4348-58.

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