

Instruction Manual

Catalog Number	PK-AB718-4015
Synonyms	RPA Interacting Protein Antibody: Replication protein A interacting protein, RPAIN, RIP, hRIP, RPA IP
Description	Replication protein A (RPA) is a single-stranded-DNA binding protein involved in numerous eukaryotic DNA processes including replication, repair and recombination. RPA interacting protein (RPA IP) has been identified as an adapter protein that is involved in RPA nuclear import instead of the prototypical importin proteins that normally mediate nuclear import. Multiple isoforms of RPA IP are known to exist, with the longest isoform localized to the cytoplasm. Isoform 2 is sumoylated and is located in the PML nuclear body within the nucleus. It has been suggested that this isoform mediates the localization of the RPA complex into the PML nuclear body, thereby participating in RPA function in DNA metabolism.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	RPA Interacting Protein antibody was raised in rabbits against a 17 amino acid peptide from near the carboxy terminus of human RPA Interacting Protein.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 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	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. RPA IP antibody can be used for detection of RPA Interacting Protein by Western blot at 0.5 - 1 µg/mL.
Images	Available upon request.
References	Zou Y, Liu Y, Wu X, et al. Functions of human replication protein A (RPA): from DNA replication to DNA damage and stress responses. <i>J. Cell Physiol.</i> 2006; 208:267-73. 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. <i>EMBO J.</i> 1999; 18:4348-58. Park J, Seo T, Kim H, et al. Sumoylation of the novel protein hRIPbeta is involved in replication protein A deposition in PML nuclear bodies. <i>Mol. Cell Biol.</i> 2005; 25:8202-14.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4015P; RPA Interacting Protein Peptide Cat.No. PK-AB718-1205; Jurkat Cell Lysate

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