

PARC antibody (pAb)

Rabbit Anti-Human/Mouse PARC (p53 Associated Parkin-Like Cytoplasmic Protein)

Instruction Manual

Catalog Number	PK-AB718-3171
Synonyms	PARC Antibody: p53-associated parkin-like cytoplasmic protein, UbcH7-associated protein 1
Description	The continued localization of p53 to the nucleus is essential for its function as a tumor suppressor. PARC, a large, Parkin-like ubiquitin ligase has recently been identified as a cytoplasmic anchor protein in p53-associated protein complexes. In the absence of stress, PARC inactivation results in nuclear localization of p53 and activation of p53-dependent apoptosis, while overexpression of this protein promoted cytoplasmic sequestration of p53. Surprisingly, PARC knockout mice were viable and exhibited no obvious phenotype, suggesting that other proteins, such as the highly related cullin family of E3 ubiquitin ligases, may perform similar functions in the absence of PARC. Additionally, it has been suggested that p53 binding to PARC may serve to control PARC function.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	PARC antibody was raised against a 17 amino acid peptide from near the carboxy terminus of human PARC.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. PARC antibody can be used for the detection of PARC by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunocytochemistry starting at 1 µg/mL. For immunofluorescence start at 2 µg/mL.
Images	Available upon request.
References	Nikolaev AY, Li M, Puskas N, et al. Parc: a cytoplasmic anchor for p53. Cell 2003; 112:29-40. Skaar JR, Arai T, and DeCaprio JA. Dimerization of CUL7 and PARC is not required for all CUL7 functions and mouse development. Mol. Cell. Biol. 2005; 25:5579-89.
Images	Available upon request.
Related Products	Daudi Lysate, Cat. No. PK-AB718-1224 PARC Peptide, Cat. No. PK-AB718-3171P

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