

ARF-BP1 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat ARF-BP1

Instruction Manual

Catalog Number	PK-AB718-4213
Synonyms	ARF-BP1 Antibody; ARF binding protein 1; HECT, UBA and WWE domain containing 1; HUWE1; Mcl-1 ubiquitin ligase E3; Mule
Description	The ARF tumor suppressor is a critical regulator of p53 stability. In addition to p53, ARF1 binds to other proteins such as MDM2 and ARF-BP1, a large protein containing HECT, UBA and WWE motifs. ARF-BP1 directly binds and ubiquitinates p53; this activity is inhibited by ARF, indicating that ARF-BP1 is a critical mediator of the p53-dependent and p53-independent tumor suppressor functions of ARF. ARF-BP1 can also catalyze the polyubiquitination of Mcl-1, an anti-apoptotic Bcl-2 family member involved in DNA damage-induced apoptosis. Elimination of ARF-BP1 expression by RNA interference stabilized Mcl-1 protein, resulting in an attenuation of apoptosis induced by DNA-damage agents.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	ARF-BP1 antibody was raised in rabbits against a 19 amino acid peptide from near the carboxy terminus of human ARF-BP1.
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, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. ARF-BP1 antibody can be used for detection of ARF-BP1 by Western blot at 1 µg/mL. Antibody can also be used for immunocytochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Gallagher SJ, Kefford RF and Rizos H. The ARF tumour suppressor. <i>Int. J. Biochem. Cell Biol.</i> 2006; 38:1637-41. Pomerantz J, Schreiber-Agus N, Leigeois NJ, et al. The Ink4a tumor suppressor gene product, p19ARF interacts with MDM2 and neutralizes MDM2's inhibition of p53. <i>Cell</i> 1998; 92:725-34. Chen D, Kon N, Li M, et al. ARF-BP1/Mule is a critical mediator of the ARF tumor suppressor. <i>Cell</i> 2005; 121:1071-83. Zhong Q, Gao W, Du F, et al. Mule/ARF-BP1, a BH3-only E3 ubiquitin ligase, catalyzes the polyubiquitination of Mcl-1 and regulates apoptosis. <i>Cell</i> 2005; 121:1085-95.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4213P; ARF-BP1 Peptide Cat.No. PK-AB718-1224; Daudi Cell Lysate

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