

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-4331 |
| Synonyms | BRCC36 Antibody: BRCC36, BRCA1, BRCA2-containing complex subunit 3, BRCC3 |
| Description | Together with the breast and ovarian predisposition proteins BRCA1 and BRCA2 and RAD51 and BRCC45, BRCC36 forms a holoenzyme complex that possesses a ubiquitin E3 ligase activity. Aberrant levels of BRCC36 were detected in sporadic breast tumors and depletion of either BRCC36 or BRCC45 by siRNA resulted in increased sensitivity to ionizing radiation and defects in the G2/M checkpoint, indicating that BRCC36 acts to enhance cellular survival following DNA damage. Recent experiments have shown that BRCC36 is essential for ionizing radiation-induced BRCA1 phosphorylation and nuclear foci formation, suggesting that BRCC36 may be an important target in the treatment of radiation-resistant breast tumors. At least two isoforms of BRCC36 are known to exist. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | BRCC36 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human BRCC36. |
| 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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. BRCC36 antibody can be used for detection of BRCC36 by Western blot at 0.5 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL. |
| Images | Available upon request. |
| References | Dong Y, Hakimi MA, Chen X, et al. Regulation of BRCC, a holoenzyme complex containing BRCA1 and BRCA2, by a signalsome-like subunit and its role in DNA repair. <i>Mol. Cell</i> 2003; 12:1087-99. Chen X, Arciero CA, Wang C, et al. BRCC36 is essential for ionizing radiation-induced BRCA1 phosphorylation and nuclear foci formation. <i>Cancer Res.</i> 2006; 66:5039-46. |
| Images | Available upon request. |
| Related Products | Cat.No. PK-AB718-4331P; BRCC36 Peptide Cat.No. PK-AB718-1219; MCF7 Cell Lysate |

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