

BRCC45 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat BRCC45

Instruction Manual

Catalog Number	PK-AB718-4329
Synonyms	BRCC45 Antibody; BRCA1, BRCA2-containing complex subunit 45, Brain and reproductive organ expressed, BRE, TNFRSF1A modulator
Description	BRCC45 was initially suggested to be a housekeeping protein that is highly expressed in brain and reproductive organs. Later experiments indicated BRCC45 forms a complex with the breast and ovarian predisposition proteins BRCA1 and BRCA2 as well as RAD51 and BRCC36. This complex has a ubiquitin E3 ligase activity and is thought to enhance cellular survival following DNA damage. BRCC45 has also been suggested to function as a death receptor-associated anti-apoptotic protein by inhibiting the BID-induced activation of the mitochondrial apoptotic pathway. Higher levels of BRCC45 were detected in the majority of hepatocellular carcinomas, suggesting that BRCC45 may promote tumorigenesis when overexpressed. At least three isoforms of BRCC45 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	BRCC45 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human BRCC45.
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. BRCC45 antibody can be used for detection of BRCC45 by Western blot at 1 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	Li L, Yoo H, Beckker FF, et al. Identification of a brain- and reproductive-organs-specific gene responsive to DNA damage and retinoic expression. <i>Biochim. Biophys. Res. Commun.</i> 1995; 206:764-74. 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. Li Q, Ching AK, Chan BC, et al. A death receptor-associated anti-apoptotic protein, BRE, inhibits mitochondrial apoptotic pathway. <i>J. Biol. Chem.</i> 2004; 279:52106-16. Chan BC, Ching AK, To KF, et al. BRE is an antiapoptotic protein in vivo and overexpressed in human hepatocellular carcinoma. <i>Oncogene</i> 2007; epub.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4329P; BRCC45 Peptide Cat.No. PK-AB718-1201; HeLa Cell Lysate

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