

CCDC98 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse CCDC98 (NT)

Instruction Manual

Catalog Number	PK-AB718-4321
Synonyms	CCDC98 Antibody; Coiled-coil domain 98, Abraxas 1, ABRA1
Description	CCDC98, also known as Abraxas 1, was identified through protein binding studies using the breast and ovarian predisposition protein BRCA1 as the binding target. CCDC98 recruits RAP80, a ubiquitin-binding protein, to BRCA1, allowing the formation of BRCA1 foci in response to DNA damage caused by ionizing radiation. Both CCDC98 and RAP80 are required for DNA damage resistance, G2-M checkpoint control, and DNA repair. Cells depleted of either CCDC98 or RAP80 exhibited increased sensitivity to ionizing radiation, although not as much as in BRCA1-depleted cells, suggesting that CCDC98 and RAP80 control only part of the DNA damage response role of BRCA1. At least two isoforms of CCDC98 are known to exist.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	CCDC98 antibody was raised in rabbits against a 15 amino acid peptide from near the amino terminus of human CCDC98.
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, IHC Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. CCDC98 antibody can be used for detection of CCDC98 by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 5 µg/mL.
Images	Available upon request.
References	Wang B, Matsuoka S, Balliff BA, et al. Abraxas and RAP80 form a BRCA1 protein complex required for the DNA damage response. <i>Science</i> 2007; 316:1194-1198. Kim H, Huang J, and Chen J. CCDC98 is a BRCA1-BRCT domain-binding protein involved in the DNA damage response. <i>Nat. Struct. Mol. Biol.</i> 2007; 14:710-5. Liu Z, Wu J, and Yu X. CCDC98 targets BRCA1 to DNA damage sites. <i>Nat. Struct. Mol. Biol.</i> 2007; 14:716-20.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4321P; CCDC98 Peptide Cat.No. PK-AB718-1311 - Human Breast Tissue Lysate

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