

BRCA-2 antibody (pAb)

Rabbit Anti-Human/Mouse BRCA-2

Instruction Manual

Catalog Number	PK-AB577-3675
Description	BRCA-1 and -2 are breast cancer susceptibility proteins that play a role in DNA damage, repair, cell cycle progression, transcription and apoptosis, etc. The BRCA-1 gene is expressed in numerous tissues, including breast and ovary, and encodes a predicted protein of 1863 amino acids. BRCA2 has been shown to be required for localization of Rad51 to sites of DSBs (double stranded breaks) in DNA. Cells lacking BRCA-1 and BRCA-2 cannot repair DSBs through the Rad51-dependent process.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide mapping to the N-terminus of human BRCA2.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) affinity purified rabbit anti-BRCA2 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml). The optimal concentrations should be determined individually. Other applications have not yet been determined. HeLa cell lysate can be used as a positive control.

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