

BRCA-1 antibody (pAb)

Rabbit Anti-Human/Mouse/Rat BRCA-1

Instruction Manual

Catalog Number	PK-AB577-3364
Synonyms	BRCA1, BRCA-1, breast cancer 1, breast and ovarian cancer susceptibility protein 1, RNF53, Ring Finger 53, IRIS, PSCP.
Description	BRCA-1 is a breast cancer susceptibility gene that was localized to chromosome 17q. Mutations within this gene are believed to account for approximately 45% of families with high incidence of breast cancer and at least 80% of families with increased incidence of both early-onset breast cancer and ovarian cancer. The BRCA-1 gene is expressed in numerous tissues, including breast and ovary, and encodes a predicted protein of 1863 amino acids. Like many other genes involved in familial cancer, BRCA-1 appears to encode a tumor suppressor, a protein that acts as a negative regulator of tumor growth.
Quantity	100 µg
Source / Host	Rabbit
Subclass	Rabbit IgG
Immunogen	Synthetic peptide surrounding amino acid 1852 of human BRCA1.
Specific Species React.	human, mouse, rat
Cross-Reactivity	See Applications or Species Reactivity.
Formulation	0.5 mg/ml in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 5 mM EDTA and 0.03% proclin.
Stabilizer	thimerosal
Concentration	0.2 mg/ml
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Applications	Western blotting (1-4 µg/ml), immunoprecipitation (10-20 µg/ml), and immunohistochemistry (10-20 µg/ml). The optimal concentrations should be determined individually. Antibody might be suitable for other applications not tested so far.
Storage	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Shipping Conditions	2-8°C

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