

## BRCA1 (Ab-1524) Antibody [#B7031]

<b>Catalog Number:</b>	B7031
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P38398
<b>Other Names:</b>	BRCA1; Breast cancer type 1 susceptibility protein; familial breast/ovarian cancer gene 1
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human BRCA1 around the phosphorylation site of serine 1524 (Y-P-S <sup>P</sup> -Q-E).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	BRCA1 (Ab-1524) antibody detects endogenous levels of total BRCA1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Miki Y., Science 266:66-71(1994). Smith T.M., Genome Res. 6:1029-1049(1996). Wilson C.A., Oncogene 14:1-16(1997).