

## RAD51L1 Antibody [#C13106]

<b>Catalog Number:</b>	C13106
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	O15315
<b>Other Names:</b>	DNA repair protein RAD51 homolog 2, R51H2, RAD51-like protein 1, Rad51B, RAD51L1, RAD51B, REC2
<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 peptide derived from internal of human RAD51L1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	RAD51L1 antibody detects endogenous levels of total RAD51L1 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Rice M.C., Proc. Natl. Acad. Sci. U.S.A. 94:7417-7422(1997). Albala J.S., Genomics 46:476-479(1997). Cartwright R., Nucleic Acids Res. 26:1653-1659(1998).