

## XRCC1 Antibody [#C0395]

<b>Catalog Number:</b>	C0395
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P18887
<b>Other Names:</b>	DNA-repair protein XRCC1, XRC1
<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 XRCC1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	XRCC1 antibody detects endogenous levels of total XRCC1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
<b>References:</b>	Ming-Whei Yu, J Natl Cancer Inst, Oct 2003; 95: 1485 - 1488. Nicolas Lévy, Nucleic Acids Res., Jan 2006; 34: 32 - 41. Claire Seedhouse, Blood, Nov 2002; 100: 3761 - 3766. Eric J. Duell, Cancer Res., Aug 2002; 62: 4630 - 4636.