

XRCC3 Antibody [#C13130]

Catalog Number:	C13130
Concentration:	1mg/ml
Swiss-Prot No.:	O43542
Other Names:	DNA-repair protein XRCC3, X-ray repair cross-complementing protein 3
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human XRCC3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	XRCC3 antibody detects endogenous levels of total XRCC3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Liu N., Mol. Cell 1:783-793(1998). Miller K.A., J. Biol. Chem. 277:8406-8411(2002). Liu N., Nucleic Acids Res. 30:1009-1015(2002).