

MRE11A Antibody [#C13088]

Catalog Number:	C13088
Concentration:	1mg/ml
Swiss-Prot No.:	P49959
Other Names:	Double-strand break repair protein MRE11A, MRE11 homolog 1, MRE11 meiotic recombination 11 homolog A, MRE11A, HNGS1, MRE11
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 MRE11A.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MRE11A antibody detects endogenous levels of total MRE11A protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Petrini J.H.J., Genomics 29:80-86(1995). Paull T.T., Mol. Cell 1:969-979(1998). Pitts S.A., Hum. Mol. Genet. 10:1155-1162(2001).