

MLH3 Antibody [#C13087]

Catalog Number:	C13087
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UHC1
Other Names:	DNA mismatch repair protein Mlh3, MutL protein homolog 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 MLH3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MLH3 antibody detects endogenous levels of total MLH3 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Lipkin S.M., Nat. Genet. 24:27-35(2000). Kondo E., Nucleic Acids Res. 29:1695-1702(2001). Lipkin S.M., Hum. Mutat. 17:389-396(2001).