

MLH1 Antibody [#C13086]

Catalog Number:	C13086
Concentration:	1mg/ml
Swiss-Prot No.:	P40692
Other Names:	DNA mismatch repair protein Mlh1, MutL protein homolog 1, MLH1, COCA2
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 MLH1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	MLH1 antibody detects endogenous levels of total MLH1 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Bronner C.E., Nature 368:258-261(1994). Papadopoulos N., Science 263:1625-1629(1994). Kolodner R.D., Cancer Res. 55:242-248(1995).