

## MSH6 Antibody [#C13091]

|                           |                                                                                                                                                        |
|---------------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Catalog Number:</b>    | C13091                                                                                                                                                 |
| <b>Concentration:</b>     | 1mg/ml                                                                                                                                                 |
| <b>Swiss-Prot No.:</b>    | P52701                                                                                                                                                 |
| <b>Other Names:</b>       | DNA mismatch repair protein Msh6, MutS-alpha 160 kDa subunit, G/T mismatch-binding protein, GTBP, GTMBP, p160, MSH6, GTBP                              |
| <b>Storage/Stability:</b> | Store at -20°C/1 year                                                                                                                                  |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), 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 MSH6.                                                            |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                                  |
| <b>Specificity:</b>       | MSH6 antibody detects endogenous levels of total MSH6 protein.                                                                                         |
| <b>Reactivity:</b>        | Human, Mouse                                                                                                                                           |
| <b>Applications:</b>      | WB: 1:500~1:1000   ELISA: 1:10000                                                                                                                      |
| <b>References:</b>        | Acharya S., Proc. Natl. Acad. Sci. U.S.A. 93:13629-13634(1996).<br>Shiwaku H.O., DNA Res. 4:359-362(1997).<br>Palombo F., Science 268:1912-1914(1995). |