

## HMG17 Antibody [#B1039]

|                           |  |
|---------------------------|--|
| <b>Catalog Number:</b>    | B1039  |
| <b>Concentration:</b>     | 1mg/ml   |
| <b>Swiss-Prot No.:</b>    | P05204   |
| <b>Other Names:</b>       | HMG17, high mobility group protein N2, high-mobility group (nonhistone chromosomal) protein 17, high-mobility group nucleosomal binding domain 2, high-mobility group nucleosome binding domain 2, nonhistone chromosomal protein HMG-17 |
| <b>Storage/Stability:</b> | Store at -20°C/1 year  |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), 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 HMG17.   |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  |
| <b>Specificity:</b>       | HMG17 antibody detects endogenous levels of total HMG17 protein.   |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%);<br>Rat (Identities = 100%, Positives = 100%)   |
| <b>Applications:</b>      | IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000   |
| <b>References:</b>        | Michael Bustin. Nucleic Acids Res., Jun 1991; 19: 3115 - 3121.<br>Julio E. Herrera. Mol. Cell. Biol., May 1999; 19: 3466 - 3473.<br>Robert Hock. J. Cell Biol., Dec 1998; 143: 1427.   |