

## Histone H2A Antibody [#B8247]

|                           |  |
|---------------------------|--|
| <b>Catalog Number:</b>    | B8247  |
| <b>Concentration:</b>     | 1mg/ml   |
| <b>Swiss-Prot No.:</b>    | P0C0S8/P04908/P20671/Q96KK5/Q6F113/Q7L7L0  |
| <b>Other Names:</b>       | H2A.1b; H2A/c; H2A/d; H2A/i; H2A/n; H2A/p; H2A1; H2AC; H2AFC; H2AFD ANDHIST1H2AL; H2AFI; H2AFN ANDHIST1H2AG; H2AFP; HIST1H2AI; HIST1H2AJ; HIST1H2AM; Histone H2A.c/d/i/n/p |
| <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 Histone H2A.   |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.  |
| <b>Specificity:</b>       | Histone H2A antibody detects endogenous levels of total Histone H2A protein.   |
| <b>Reactivity:</b>        | Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)  |
| <b>Applications:</b>      | IHC: 1:50~1:100 ELISA: 1:5000  |
| <b>References:</b>        | Albig W., Hum. Genet. 101:284-294(1997).<br>Albig W., Biol. Chem. 380:7-18(1999).<br>Dobner T., DNA Seq. 1:409-413(1991).  |