

---

## Histone H2BK5me1 (H2BK5 Monomethyl) Polyclonal Antibody

(Catalog #A68344)

---

### Background

Variant histone specifically required to direct the transformation of dissociating nucleosomes to protamine in male germ cells (By similarity). Entirely replaces classical histone H2B prior nucleosome to protamine transition and probably acts as a nucleosome dissociating factor that creates a more dynamic chromatin, facilitating the large-scale exchange of histones (By similarity). Core component of nucleosome (By similarity). Nucleosomes wrap and compact DNA into chromatin, limiting DNA accessibility to the cellular machineries which require DNA as a template (By similarity). Histones thereby play a central role in transcription regulation, DNA repair, DNA replication and chromosomal stability (By similarity). DNA accessibility is regulated via a complex set of post-translational modifications of histones, also called histone code, and nucleosome remodeling (By similarity). Also found in fat cells, its function and the presence of post-translational modifications specific to such cells are still unclear.

### Description

Histone H2BK5me1 (H2BK5 Monomethyl) Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

PBS with 0.02% sodium azide, 50% glycerol, pH 7.4.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q96A08/P33778/P62807

### Purification

Affinity Purified

### Immunogen

Synthetic Peptide

### Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

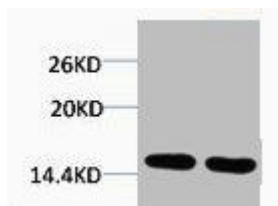
### Alternative Names

HIST1H2BA; TSH2B; Histone H2B type 1-A; Histone H2B, testis; Testis-specific histone H2B; HIST1H2BB; H2BFF; Histone H2B type 1-B; Histone H2B.1; Histone H2B.f; H2B/f; HIST1H2BC; H2BFL; HIST1H2BE; H2BFH; HIST1H2BF; H2BFG; HIST1H2BG; H2BFA; HIST1H2BI; H2BFK

### Application

WB, ELISA; Recommended dilution: WB:1:500-1:1000, ELISA:1:20000

---



Western blot analysis of 1) Hela, 2) 3T3, diluted at 1:2000 using Histone H2BK5me1 (H2BK5 Monomethyl) Polyclonal Antibody.