

Recombinant Histone H2B

Catalog No: 31252

Source: *Xenopus laevis*

Quantity: 50 µg

Expressed In: *E. coli*

Contents: Lyophilized from 50 µl sterile water containing 5 mM Beta-mercaptoethanol. Supplied as lyophilized powder.

Background:

Histone H2B is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 147 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points.

Protein Details:

Recombinant *Xenopus laevis* Histone H2B produced in *E. coli* and purified using FPLC. The H2B protein is missing the first three residues (PEP). Protein concentration was determined by spectrophotometry. ≥98% purity by SDS-PAGE. The MW of the protein is 13,494 Daltons. Resuspend protein in suitable buffer for your assay.

The protein sequence is as follows:

```
AKSAPA PKKGSKKAVT KTQKKDGKKR RKS RKESYAI YVYKVLKQVH PDTGISSKAM
SIMNSFVNDV FERIAGEASR LAHYNKRSTI TSREIQTAVR LLLPGELAKH
AVSEGTKAVT KYTSAK
```

Protein Applications:

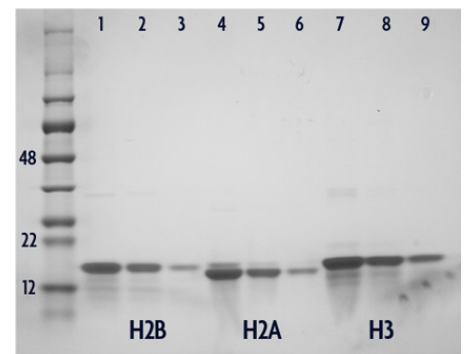
Each lot has been evaluated by SDS-PAGE and dot blot analysis.

Storage:

Lyophilized protein can be stored at -20°C. Once the protein is resuspended, we recommend making aliquots to avoid multiple freeze-thaw cycles and storing the protein at -80°C.

Guarantee:

For research use only. Product is guaranteed stable for six months from date of receipt when stored properly.



SDS-PAGE gel of Recombinant Histone H2B. 5 µg, 2 µg and 1 µg respectively of Recombinant Histone H2B, Catalog No. 31252 (lanes 1-3), Recombinant Histone H2A, Catalog No. 31251 (lanes 4-6) and Recombinant Histone H3, Catalog No. 31207 (lanes 7-9).