

## Recombinant Histone H2B (Human)

**Catalog No:** 31492, 31892

**Expressed In:** *E. coli*

**Quantity:** 100, 1000 µg

**Concentration:** 2 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant Histone H2B (Human) was expressed in *E. coli* and is supplied in 25 mM Tris pH 8.0, 150 mM NaCl, 5% Glycerol. Please refer to product insert upon arrival for lot-specific concentration.

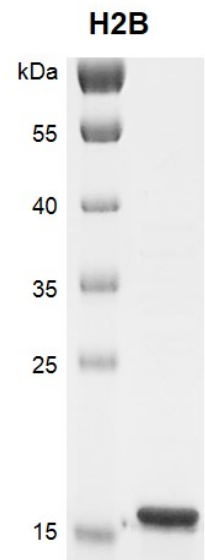
**Background: Histone H2B** is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points.

**Protein Details:** Recombinant Histone H2B (Human) (assessment number: NP\_003509.1) was expressed in *E. coli* cells. This protein does not contain any tag with a molecular weight of 13.9 kDa. The recombinant protein is >85% pure by SDS-PAGE.

**Application Notes:** Recombinant Histone H2B (Human) is suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**Recombinant Histone H2B (Human) gel.**  
Recombinant Histone H2B run on an 15% SDS-PAGE gel and stained with Coomassie blue.