

# Histone H2B, Recombinant Human

**Catalog No** 15-0302  
**Lot No** 21015002-25  
**Pack Size** 100 µg

**Type** Histone  
**Expressed In** *E. coli*  
**Mol. Wgt.** 14 kDa  
**Epitope Tag** None

## Product Description:

Recombinant human histone H2B (HIST1H2BK, accession O60814) expressed in *E. coli* and purified by FPLC. Histone H2B is one of the four proteins that are present in the nucleosome, the basic repeating unit subunit of chromatin, consisting of 147 base pairs of DNA wrapped around an octamer of core histone proteins (H2A, H2B, H3 and H4).

## Formulation:

100 µg of lyophilized powder.

## Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.

## Application Notes:

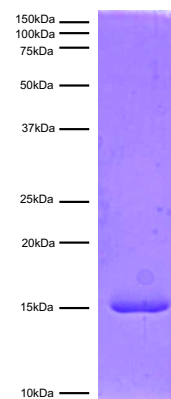
Recombinant histone H2B is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

## References:

This product is for *in vitro* research use only and is not intended for use in humans or animals.



# EpiCypher®



**Protein Gel Data:** Histone H2B, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.