
H2B1 Protein

(Catalog No. E35002)

Description

Histone H2B combines with Histone H2A to form the H2A-H2B heterodimer. Two H2A/H2B heterodimers interact with an H3/H4 tetramer to form the histone octamer. (1,2) Histone H2B is also modified by various enzymes and can act as a substrate for them. These modifications have been shown to be important in gene regulation.

Molecular Weight (kDa)

13.8 kDa

Formulation

20 mM sodium phosphate, 300 mM NaCl, 1 mM EDTA

Source

Expressed in E.coli expression system

Size

50 µg

Storage Temp & Conditions

Store stock solution at -20°C. Avoid multiple freeze/thaw cycles

Purification Notes

SDS-PAGE purified

Application Notes

Enzyme-linked Immunoabsorbent Assay; Western Blot (Recombinant protein); Antibody Production; Protein Array

Ordering Information:

Products	Size	Cat. No.
H2B1 Protein	50 µg	E35002 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.