

## Recombinant Histone H2A.Z-H2B dimers

**Cat. No.** C23010012

**Source:** E. coli

**Lot #:** 001

**Size:** 50 µg/ 50 µl

**Concentration:** 1 µg/µl

**Specificity:** Human

**Purity:** Purified using FPLC, >98% purity as determined by SDS-PAGE

**Storage buffer:** 20 mM Tris-Cl pH 7.9, 1 M NaCl, 1mM EDTA, 0.5 mM PMSF and 1 mM DTT.

**Storage:** Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

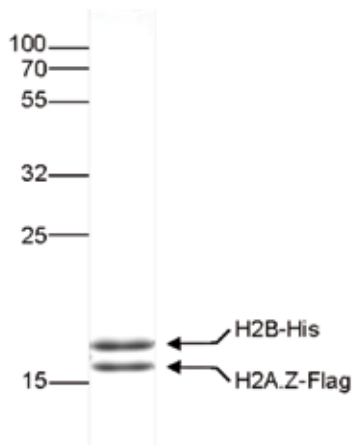
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Dimer of full length N-terminal Flag tagged recombinant histone H2A.Z and N-terminal His tagged recombinant histone H2B, produced in E. coli.

### Protein description

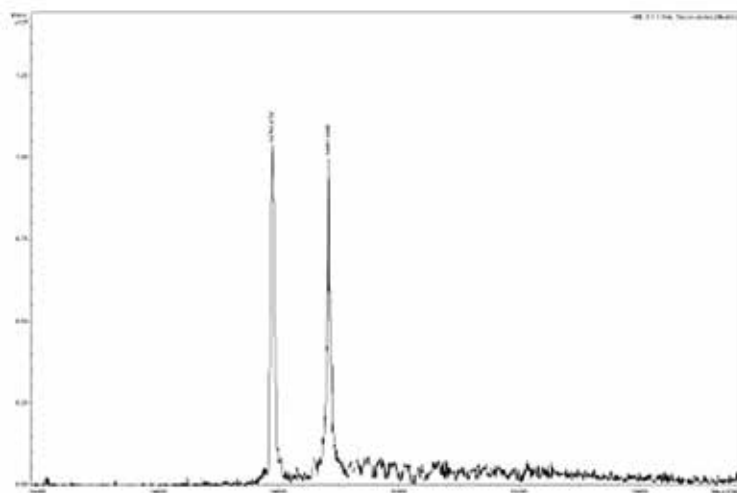
Histones are the main constituents of the protein part of chromosomes of eukaryotic cells. They are rich in the amino acids arginine and lysine and have been greatly conserved during evolution. Histones pack the DNA into tight masses of chromatin. Two core histones of each class H2A, H2B, H3 and H4 assemble and are wrapped by 146 base pairs of DNA to form one octameric nucleosome. Histone tails undergo numerous post-translational modifications, which either directly or indirectly alter chromatin structure to facilitate transcriptional activation or repression or other nuclear processes. The histone variant H2A.Z replaces conventional H2A in a subset of nucleosomes. H2A.Z is involved in transcriptional regulation, antisilencing, silencing, and genome stability. It functions as a key regulator of chromatin function and plays an essential role during mammalian development.

### Quality control



**Figure 1.**

SDS page of the Recombinant Histone H2A.X-H2B dimers. The position of the proteins of interest is indicated on the right; the marker (in kDa) is shown on the left.



**Figure 2.**

ESI-TOF analysis of the Recombinant Histone H2A.X-H2B dimers.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** September 7, 2015