

# Histone H2A/H2B Dimer, Recombinant Human



## EpiCypher®

**Catalog No** 15-0311  
**Lot No** 22122004-04  
**Pack Size** 50 µg

**Type** Histone  
**Expressed In** *E. coli*  
**Mol. Wgt.** 27.8 kDa  
**Epitope Tag** None

### Product Description:

Histone H2A/H2B dimer made from recombinant histones expressed in *E. coli* (accession numbers: H2A-P04908; H2B-O60814). Histones H2A and H2B were expressed and purified individually by FPLC, and then assembled into dimers that were further purified using gel filtration chromatography.

### Formulation:

50 µg of H2A/H2B dimer (1.18 mg/mL) in 10 mM Tris-HCl pH 7.5, 2 M NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol.

### Storage and Stability:

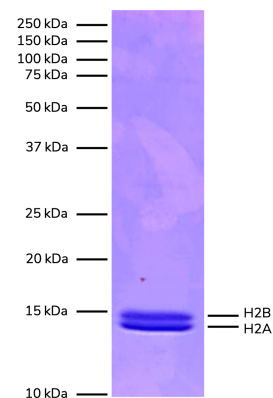
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

### Application Notes:

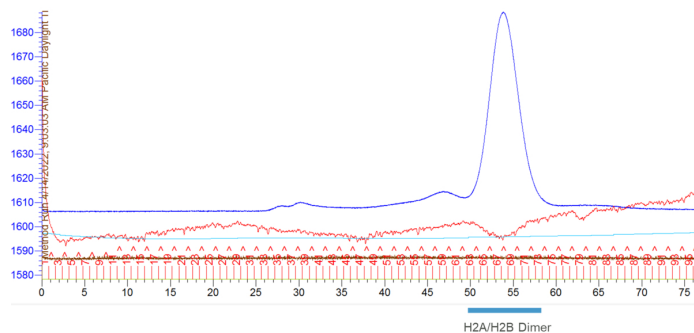
Recombinant histone H2A/H2B dimer is suitable for enzyme assays and nucleosome reconstitution.

### References:

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



**Figure 1: Protein gel data.** Recombinant H2A/H2B dimer (2 µg) was resolved via SDS PAGE and stained with Coomassie blue. The migration and molecular weight of the protein standards and migration of H2A and H2B are indicated.



**Figure 2: Purification data.** Chromatogram from gel filtration purification of histone H2A/H2B dimer. Only fractions corresponding to the intact dimer (blue bar) were collected and pooled.