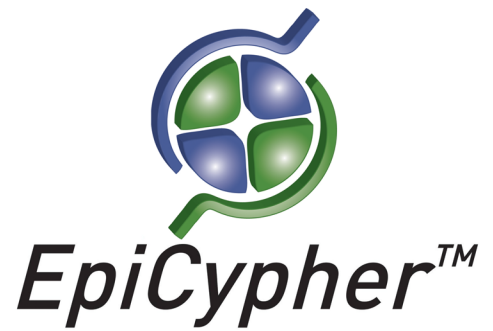


Histone H2A/H2B Dimer, Recombinant Human

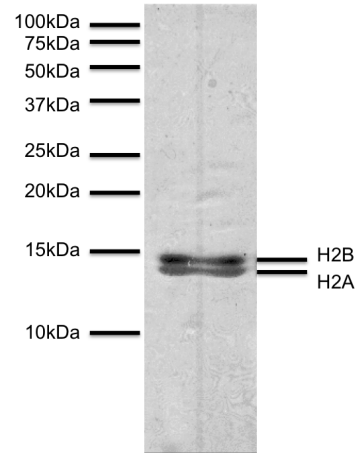
Catalog No. 15-0311
Lot No. 15259001
Pack Size 50 µg



Type Histone **Expressed In** *E. coli*
Mol. Wgt. 29 kDa **Epitope Tag** None

Product Description:

Histone H2A/H2B dimers 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, then assembled into dimers that were further purified using gel filtration chromatography.



Protein Gel Data: Histone H2A/H2B Dimer, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards and migration of H2A and H2B are indicated.

Formulation:

50 µg of H2A/H2B dimer (0.5 mg/ml) in 10mM Tris-HCl pH 7.5, 2M NaCl, 1mM EDTA, 2mM DTT, 20% glycerol.

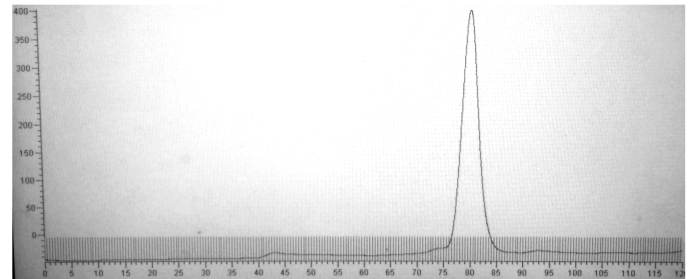
Storage and Stability:

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. Reconstitute with distilled water or suitable buffer prior to usage.

References:



Purification Data: Chromatogram from gel filtration purification of Histone H2A/H2B Dimer, Recombinant Human. Only fractions corresponding to the intact dimer (as indicated, black bar) were collected and pooled.

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