

Recombinant Histone H2A/H2B dimer

Catalog No: 81167, 81867

Lot No: 18218001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 1 µg/µl

Source: Human

Buffer Contents: Recombinant Histone H2A/H2B Dimer is supplied in 10 mM Tris-HCl pH 7.4, 2 M NaCl, 1 mM EDTA, and 5 mM b-mercaptoethanol.

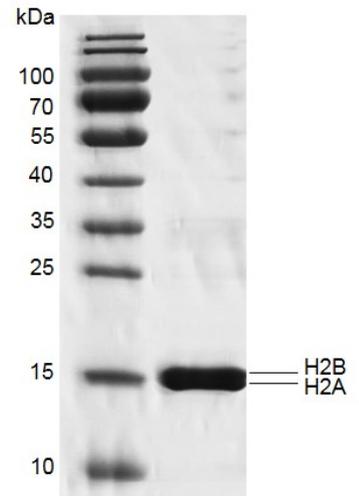
Background: Histone H2A and H2B are the core components of the nucleosome. H2A and H2B assemble H2A/H2B dimer, and two H2A/H2B dimers combine with H3/K4 tetramer to form histone octamer. The histone octamer is wrapped around by 146 bp DNA to assemble a nucleosome, which is the basic structural unit of chromatin *in vivo*.

Protein Details: Recombinant histone H2A/H2B Dimer consists of full length human histones H2A (accession number NP_003503.1) and full length human histone H2B (accession number NP_003509.1). Histone H2A and H2B were expressed in *E. coli* cells. Recombinant Histone H2A/H2B Dimer was assembled by dialysis and purified by gel filtration. The molecular weight of the Histone H2A/H2B Dimer is 27.9 kDa.

Application Notes: Recombinant Histone H2A/H2B Dimer is suitable for use as substrates in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

Histone H2A/H2B Dimer



Recombinant Histone H2A/H2B Dimer
13% SDS-PAGE Coomassie staining

MW:

H2A: 14.1 kDa

H2B: 13.8 kDa

Purity: ≥ 90%