

Recombinant Histone H3/H4 tetramer

Catalog No: 81169, 81869

Lot No: 19318001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.5 µg/µl

Source: Human

Buffer Contents: Recombinant Histone H3/H4 Tetramer is in 10 mM Tris-HCl pH 7.4, 2 M NaCl, 1 mM EDTA, 10% glycerol, and 5 mM b-mercaptoethanol.

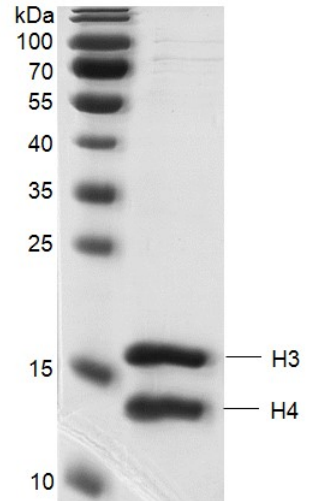
Background: Histone H3 and H4 are the core components of the nucleosome. H3 and H4 assemble H3/H4 tetramer, and H3/K4 tetramer combines with two H2A/H2B dimers to form histone octamer. The histone octamer is wrapped around by 146 bp DNA to assemble a nucleosome, which is the basic structural unit of chromatin *in vivo*.

Protein Details: Recombinant histone H3/K4 Tetramer consists of full length human histones H3 (accession number NP_003520.1) and full length human histone H4 (accession number NP_003539.1). Histone H3 and H4 were expressed in *E. coli* cells. Recombinant histone H3/H4 Tetramer was assembled by dialysis and purified by gel filtration. The molecular weight of the Histone H3/H4 Tetramer is 53 kDa.

Application Notes: Recombinant Histone H3/H4 Tetramer 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.

H3/H4 tetramer



Recombinant Histone H3/H4 Tetramer
13% SDS-PAGE Coomassie staining

MW:

H3: 15.3 kDa

H4: 11.2 kDa

Purity: ≥ 90%