

Recombinant Histone H1

Catalog No: 81126, 81826

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Source: Human

Buffer Contents: Recombinant Histone H1 in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, and 5% glycerol.

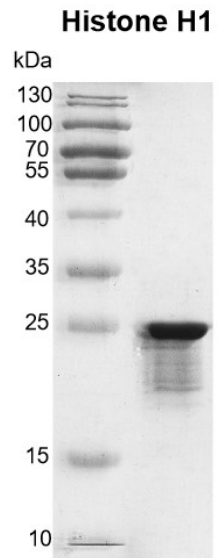
Background: Histone H1, also called as H1F0 (H1 Histone Family Member 0), is a replacement subtype, and replication independent histone, which is a member of the histone H1 family. It involves in basal repression of gene expression. H1 can act as a regulator of individual gene transcription through chromatin remodeling, nucleosome spacing and DNA methylation. Diseases associated with H1F0 include Retinal Cancer and Deafness, Autosomal Dominant.

Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker- histone H1, with the DNA between the nucleosomes to form higher order chromatin structures. Histone H1 present at the interface between the nucleosome core and DNA entry/exit points.

Protein Details: Full length Histone H1 (accession number NP_005309.1) was expressed in *E. coli* cells with no tag. The molecular weight of Histone H1 is 21 kDa.

Application Notes: This product was manufactured as described in Protein Details. Where possible, Active Motif has developed functional or activity assays for recombinant proteins. Additional characterization such as enzyme kinetic activity assays, inhibitor screening or other biological activity assays may not have been performed for every product. All available data for a given product is shown on the lot-specific Technical Data Sheet.

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.



Recombinant Histone H1 protein gel
12.5% SDS-PAGE Coomassie staining
MW: 21 kDa
Purity: >85%