

## Recombinant Histone H1T

**Catalog No: 81127, 81827**

**Lot No:** 14018001

**Expressed In:** *E. coli*

**Quantity:** 100, 1000 µg

**Concentration:** 0.6 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant Histone H1T is supplied in 25 mM Tris-HCl pH 8.0, 300 mM NaCl, 10% glycerol.

**Background:** Histone H1T, also called as HIST1H1T (Histone Cluster 1 H1 Family Member T) or Testicular H1 Histone, is a replacement subtype, and replication-independent histone, which is a member of the histone H1 family.

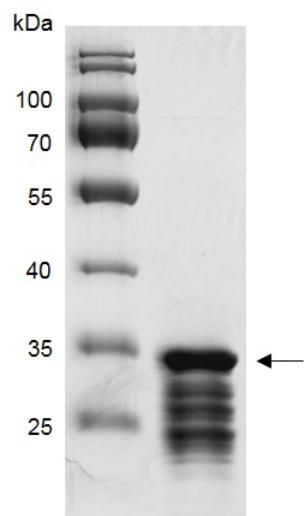
Histones are basic nuclear proteins that are responsible for the nucleosome structure of the chromosomal fiber in eukaryotes. Nucleosomes consist of approximately 146 bp of DNA wrapped around a histone octamer composed of pairs of each of the four core histones (H2A, H2B, H3, and H4). The chromatin fiber is further compacted through the interaction of a linker- histone H1, with the DNA between the nucleosomes to form higher order chromatin structures. Histone H1T forms less compacted chromatin compared to other H1 histone subtypes. Formation of more relaxed chromatin may be required to promote chromatin architecture required for proper chromosome regulation during meiosis, such as homologous recombination.

**Protein Details:** Full length Histone H1T (accession number AAA35944.1) was expressed in *E. coli* cells with no tag. The molecular weight of Histone H1T is 22 kDa.

**Application Notes:** Recombinant Histone H1T 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 H1T**



**Recombinant Histone H1T protein gel**

12.5% SDS-PAGE Coomassie staining

MW: 22 kDa

Purity: >65%