

## Recombinant Histone H4, His-Tag (Human)

**Catalog No:** 31493, 31893

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

**Quantity:** 100, 1000 µg

**Concentration:** 0.75 µg/µl

**Source:** Human

**Buffer Contents:** Full length recombinant Histone H4, His-Tag (Human) is supplied in 25 mM Tris pH 7.4, 150 mM NaCl, 5% Glycerol, and 0.04% Triton X-100.

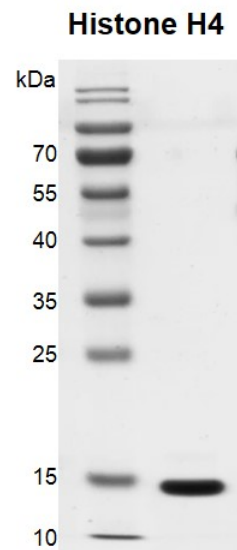
**Background: Histone H4** is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker histone, present at the interface between the nucleosome core and DNA entry/exit points.

**Protein Details:** Recombinant Histone H4, His-Tag (Human) (accession number: NP\_003539.1) was expressed in *E. coli* cells and contains a C-terminal 6×His-Tag with a molecular weight of 12.19 kDa. The recombinant histone H4 is >95% pure by SDS-PAGE.

**Application Notes:** Recombinant Histone H4, His-Tag (Human) is suitable for use as substrate for histone modification enzymes, or to generate chromatin in vitro.

**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 guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



**Recombinant Histone H4, His-Tag (Human) protein gel.**

Recombinant Histone H4, His-Tag run on a 12% SDS-PAGE gel and stained with Coomassie blue.

Purity: >95%