

## Recombinant Histone H3 acetyl Lys4 (H3K4ac)

**Catalog No:** 31275

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

**Quantity:** 25 µg µg

**Source:** Human

**Buffer Contents:** 25 µg supplied as lyophilized powder.

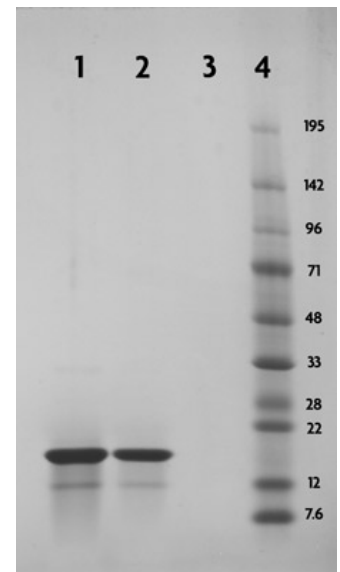
**Background:** Histone H3 is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker histone, present at the interface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

**Protein Details:** Recombinant human Histone H3.2 produced in *E. coli* and purified using HPLC. The purified protein was subsequently ligated to an acetyl lysine 4 peptide via a native peptide bond and repurified prior to lyophilization. The recombinant histone is >98% pure by SDS-PAGE and confirmed by high-resolution ESI-TOF mass spectrophotometry. The molecular weight of the protein is 15,269 Daltons. Resuspend protein in suitable buffer for your assay.

**Application Notes:** Recombinant histones are suitable for use as substrates for *in vitro* enzymatic reactions and chromatin assembly.

**Storage and Guarantee:** Lyophilized proteins can be stored at -20°C or -80°C, preferably desiccated. 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 H3 acetyl Lys4 analyzed by SDS-PAGE gel.

SDS-PAGE gel analysis of 2 µg of Recombinant Histone H3 acetyl Lys4 (Lane 1) and 1 µg Recombinant Histone H3 acetyl Lys4 (Lane 2). Lane 4 contains the molecular weight marker.