

Recombinant Histone H3K27ac

Catalog No: 31290

Expressed In: Synthetic

Quantity: 25 µg

Source: Human

Buffer Contents: 25 µg supplied as lyophilized powder. Recombinant histones can be resuspended in water or any suitable buffer. We recommend a starting concentration of 1 mg/ml. To fully solubilize the histone we suggest resuspension in the buffer of choice at room temperature for 20-30 minutes with occasional pipetting. Addition of salt or Tris to the resuspension buffer may enhance histone solubility.

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 protein, present at the interface between the nucleosome core and DNA entry/exit points.

Protein Details: Recombinant Histone H3K27ac proteins are synthetic modified histones containing acetylation at K27. The Recombinant Histone H3K27ac proteins correspond to the native histone sequence and do not contain any amino acid substitutions or residue analogs. Protein concentration was determined using Bradford assay. The recombinant histone is >95% pure by SDS-PAGE and confirmed by high-resolution ESI-TOF mass spectrometry. The molecular weight of the recombinant histone is 15,313 Daltons. To see a representative mass spectrum, please visit our website at www.activemotif.com and view the product details page.

Application Notes: Recombinant histones are suitable for use as positive controls in the analysis of histone post-translational modifications, as substrates for histone modification enzymes, or to generate chromatin *in vitro*.

References:

This product was used in the following publications:

Nat. Immunol. (2016). 17(6):695-703. PMID: 27111144. (HDAC Assay)

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.

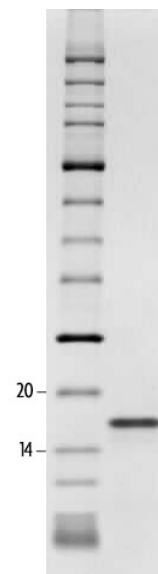


Figure 1: SDS-PAGE gel analysis of Recombinant Histone H3K27ac protein.