

Recombinant Histone H3.2 biotinylated (Human)

Catalog No: 31271

Expressed In: *E. coli*

Quantity: 25 µg

Source: Human

Buffer Contents: Recombinant Histone H3.2 biotinylated (Human) is 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.

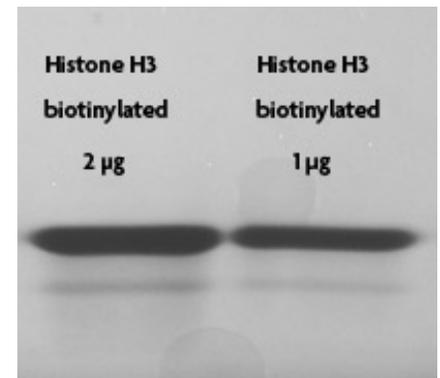
Background: Histone H3 is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker protein, present at the interface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

Protein Details: Truncated recombinant Histone H3.2 biotinylated (Human) is produced in *E. coli* and purified using FPLC. The purified protein is subsequently ligated to a peptide containing the histone tail with a N-terminal biotin via a native peptide bond. The full-length protein is then repurified prior to lyophilization. Protein concentration was determined using the molar extinction coefficient for Histone H3 and absorbance at 280nm. Protein was determined to have ≥ 98% purity by SDS-PAGE. The molecular weight is 15,565 Daltons.

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*.

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 H3.2 biotinylated (Human) tested by SDS-PAGE gel.

SDS-PAGE analysis of 1.5 µg Recombinant Histone H3 biotinylated (lane 1) and 1 µg Recombinant Histone H3 biotinylated (lane 2). Molecular weight marker is in lane 4. Lane 3 is empty.