

Recombinant Histone H3 acetyl Lys14

Catalog No: 31254

Source: Human

Quantity: 25 µg

Expressed In: *E. coli*

Contents: Lyophilized from 50 µl sterile water containing 20% acetonitrile. 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 FPLC. The purified protein was subsequently ligated to an **acetyl lysine 14** 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. Protein concentration was determined by Nanodrop. The MW of the protein is 15,266 Daltons. Resuspend protein in suitable buffer for your assay.

Protein Applications:

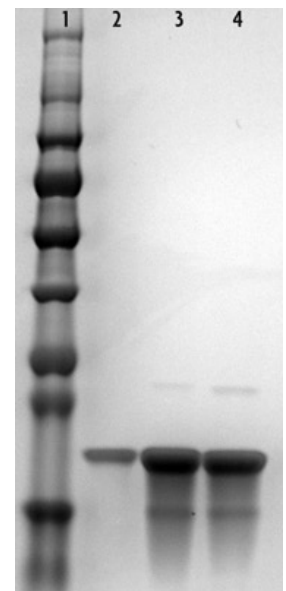
Each lot has been evaluated by SDS-PAGE.

Storage:

Lyophilized protein can be stored at -20°C. Once the protein is resuspended, we recommend making aliquots to avoid multiple freeze-thaw cycles and storing the protein at -80°C.

Guarantee:

For research use only. Product is guaranteed stable for six months from date of receipt when stored properly.



SDS-PAGE gel of Recombinant Histone H3 acetyl Lys14.

Lane 1: Marker

Lane 2: 1.5 µg Recombinant Histone H3

Lane 3: 1.5 µg Recombinant Histone H3 acetyl Lys9

Lane 4: 1.5 µg Recombinant Histone H3 acetyl Lys14