

Recombinant Histone H3 phospho Ser10

Catalog No: 31272

Source: Human

Quantity: 25 µg

Expressed In: *E. coli*

Contents: Lyophilized from 50 µl sterile water containing 30% 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 HPLC. The purified protein was subsequently ligated to a **phosphorylated serine 10** peptide via a native peptide bond and repurified prior to lyophilization. Protein concentration was determined by spectrophotometry and molar extinction coefficient. Protein was determined to have ≥ 98% purity by SDS-PAGE. The molecular weight is 15,306 Daltons.

Protein Applications:

Each lot has been evaluated by SDS-PAGE and confirmed by mass spec.

Resuspend the lyophilized pellet in sterile water, or other appropriate buffer for downstream applications. We suggest resuspension to a concentration of 1 mg/ml. It will take about 20-30 minutes for the histone to become solubilized in the resuspension buffer. We recommend adding the resuspension buffer at room temperature and periodically pipetting up and down during the 20 minute incubation to completely solubilize the histone.

Recombinant Histone H3 phospho Ser10 is suitable for use as a substrate for *in vitro* enzymatic reactions.

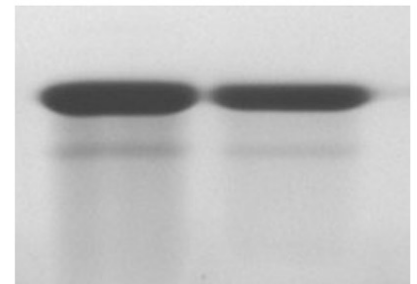
Storage:

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.

Recombinant Histone H3 phospho Ser10 2 µg Recombinant Histone H3 phospho Ser10 1 µg



SDS-PAGE analysis of 2 µg Recombinant Histone H3 phospho Ser10 (lane 1) and 1 µg Recombinant Histone H3 phospho Ser10 (lane 2).