

Histone H2A, Recombinant Human

Catalog No 15-0301
Lot No 21280002-61
Pack Size 100 µg

Type Histone **Expressed In** *E. coli*

Mol. Wgt. 14 kDa **Epitope Tag** None

Product Description:

Recombinant human histone H2A (HIST1H2AB, H2AC4, accession P04908) expressed in *E. coli* and purified by FPLC. Histone H2A is one of the four core histone proteins that are present in the nucleosome, the basic repeating unit subunit of chromatin, consisting of 147 base pairs of DNA wrapped around an octamer of core histone proteins (H2A, H2B, H3 and H4).

Formulation:

100 µg of lyophilized powder.

Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.

Application Notes:

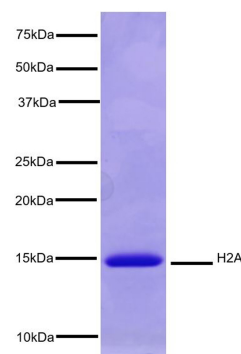
Recombinant histone H2A is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

References:

This product is for *in vitro* research use only and is not intended for use in humans or animals.



EpiCypher®



Protein Gel Data: Histone H2A, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.