

Histone H2A, Recombinant Human

Catalog No	15-0301	Species	Human
Lot No	25140003-01	Source	<i>E. coli</i>
Pack Size	100 µg	Epitope Tag	None
Concentration	N/A	MW	14 kDa

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).

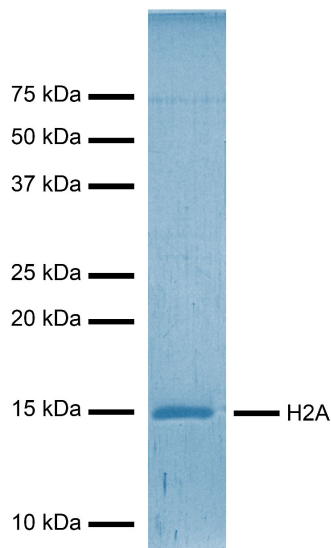
TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	100 µg of lyophilized powder.

APPLICATION NOTES

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

VALIDATION DATA



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.