

## Histone H4, Recombinant Human

<b>Catalog No</b>	15-0304	<b>Species</b>	Human
<b>Lot No</b>	24044003-03	<b>Source</b>	<i>E. coli</i>
<b>Pack Size</b>	100 µg	<b>Epitope Tag</b>	None
<b>Concentration</b>	N/A	<b>MW</b>	11 kDa

### DESCRIPTION

Recombinant human histone H4 (HIST1H4A, accession P62805) expressed in *E. coli* and purified by FPLC. Histone H4 is one of the four proteins that are present in the nucleosome, the basic repeating subunit of chromatin, consisting of 147 base pairs of DNA wrapped around an octamer of core histone proteins (H2A, H2B, H3, and H4).

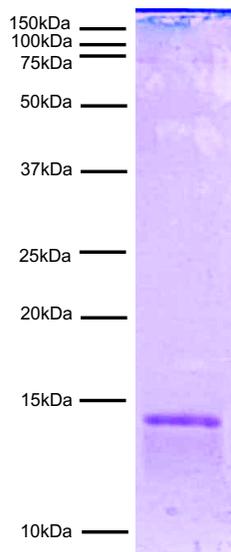
### TECHNICAL INFORMATION

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

### APPLICATION NOTES

Recombinant histone H4 is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

### VALIDATION DATA



**FIGURE 1 Protein gel data.** Histone H4 (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.