

## Histone H3.3, Recombinant Human

<b>Catalog No</b>	15-0312	<b>Species</b>	Human
<b>Lot No</b>	16067001	<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>	15 kDa

### DESCRIPTION

Recombinant human histone H3.3 (H3F3A, H3.3A, H3F3, accession P84243) expressed in *E. coli* and purified by FPLC. Histone H3 is one of the four 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). H3.3 is a histone variant, a non-allelic replacement histone found in regions of high chromatin turnover outside of S-phase (e.g. at actively transcribed genes). H3.3-H4 associates *in vivo* with the HIRA chaperone complex.

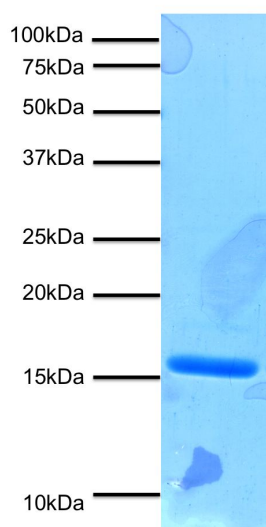
### TECHNICAL INFORMATION

<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 H3.3 is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

### VALIDATION DATA



**FIGURE 1 Protein gel data.** Histone H3.3, 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.