

Histone H3.1, Recombinant Human

Catalog No	15-0303	Species	Human
Lot No	24254004-01	Source	<i>E. coli</i>
Pack Size	100 µg	Epitope Tag	None
Concentration	N/A	MW	15 kDa

DESCRIPTION

Recombinant human histone H3.1 (HIST1H3A, accession number P68431) expressed in *E. coli* and purified by FPLC. Histone H3.1 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 (two each of H2A, H2B, H3, and H4). Histone H3.1 is the primary histone H3 variant incorporated into chromatin during cell division and DNA replication.

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 H3.1 is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to use.

VALIDATION DATA

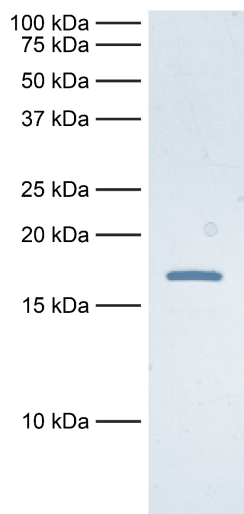


FIGURE 1 Protein gel data. Histone H3.1, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards are indicated.