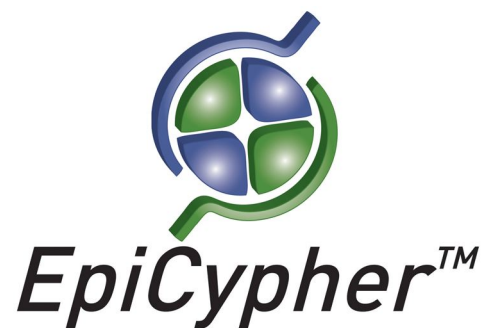


Histone H4, Recombinant Human

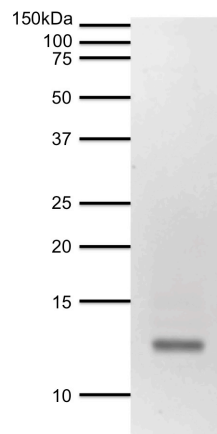
Catalog No. 15-0304
Lot No. 16223001
Pack Size 100 µg



Type Histone **Expressed In** *E. coli*
Mol. Wgt. 11 kDa **Epitope Tag** None

Product 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 unit subunit of chromatin, consisting of 147 base pairs of DNA wrapped around an octamer of core histone proteins (H2A, H2B, H3 and H4).



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.

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 multiple freeze/thaws.

Application Notes:

Recombinant histone H4 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.