

Histone H3.3K27M, Recombinant Human

Catalog No. 15-0323
Lot No. 19283001-00
Pack Size 100 µg



EpiCypher®

Type Histone **Expressed In** *E. coli*
Mol. Wgt. 15 kDa **Epitope Tag** None

Product Description:

Recombinant human histone H3.3 (H3F3A, H3.3A, H3F3, accession P84243), containing a lysine to methionine substitution at position 27, 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.

Formulation:

100 µg of lyophilized powder.

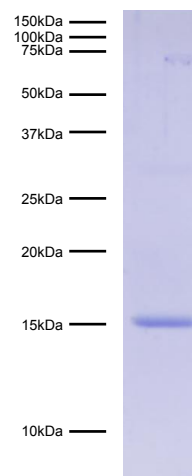
Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Recombinant histone H3.3K27M is suitable for enzyme assays and nucleosome reconstitution. Reconstitute with distilled water prior to usage.

References:



Protein Gel Data: Histone H3.3K27M, 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.

This product is for *in vitro* research use only and is not intended for use in humans or animals.

EpiCypher, Inc • Phone: 855-374-2461 • Fax: 855-420-6111 • www.epicypher.com