

Recombinant Histone H3K36me1 (MLA)

Catalog No: 31622, 31722

Lot No: 08917001

Expressed In: *E. coli*

Quantity: 100, 1000 µg

Concentration: 0.75 µg/µl

Source: Human

Buffer Contents: Recombinant Histone H3K36me1 (MLA) is supplied in 50 mM Tris-HCl, pH 8.0, 150 mM NaCl and 5% glycerol.

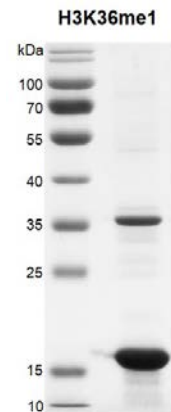
Background: Histone H3 is one of the core components of the nucleosome. The nucleosome is the smallest subunit of chromatin and consists of 146 base pairs of DNA wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3 and H4). Histone H1 is a linker protein, present at the interface between the nucleosome core and DNA entry/exit points.

Recombinant Histone H3K36me1 (MLA) has been generated using the patented Methylated Lysine Analog (MLA) technology. In MLA, methylated histones are generated via a chemical alkylation reaction that substitutes a methylated analog of lysine, aminoethylcysteine, for the existing lysine at the desired residue. Aminoethylcysteine is structurally and chemically similar to lysine, though it contains a sulfide substitution in place of the lysine γ -methylene. The MLA technique provides precise control over the site and degree of methylation. The MLA technology is covered under U.S. Patent No. 8,278,112.

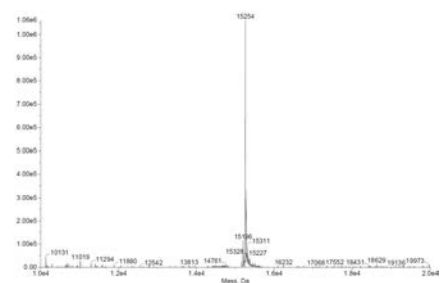
Protein Details: Recombinant human Histone H3 monomethyl Lys36 (H3K36me1) (accession number NP_003520.1) was expressed in *E. coli* cells and contains two substitutions (cysteine to serine at amino acid 96 and cysteine to alanine at amino acid 110). Recombinant methylated histones are specifically methylated via a chemical alkylation reaction that introduces a methyl lysine analog (MLA). The molecular weight of the protein is 15257 Daltons.

Application Notes: Recombinant histones are suitable for use as positive controls in the analysis of histone post-translational modifications, as substrates for histone modification enzymes, or to generate chromatin *in vitro*.

Storage and Guarantee: Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.



Recombinant Histone H3K36me1 (MLA) protein gel.
12% SDS-PAGE Coomassie staining.
Purity: > 78%



ESI-TOF Mass Spec analysis for Recombinant Histone H3K36me1 (MLA).

Expected mass: 15257 Da

Observed mass: 15257 Da