

Nucleosome, Recombinant Human, H3K9bu dNuc, Biotinylated



EpiCypher®

Catalog No 16-0371
Lot No 21178003-63
Pack Size 50 µg

Product Description:

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4; accession numbers: H2A-P04908; H2B-O60814; H3.2-Q71DI3*; H4-P62805) wrapped by 147 base pairs of 601 positioning sequence DNA. Histone H3 (created by a proprietary synthetic method) contains butyryl-lysine at position 9. The nucleosome is the basic subunit of chromatin. The 147 bp 601 sequence, identified by Lowary and Widom [1], has high affinity for histone octamers and is useful for nucleosome assembly. The DNA contains a 5' biotin-TEG group.

*H3K9bu also has a Cys to Ala substitution at position 110.

Formulation:

H3K9bu dNuc (27.4 µg protein weight, 50 µg DNA + protein) in 50.4 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molarity = 4.95 µM. MW = 199,903.5 Da

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:

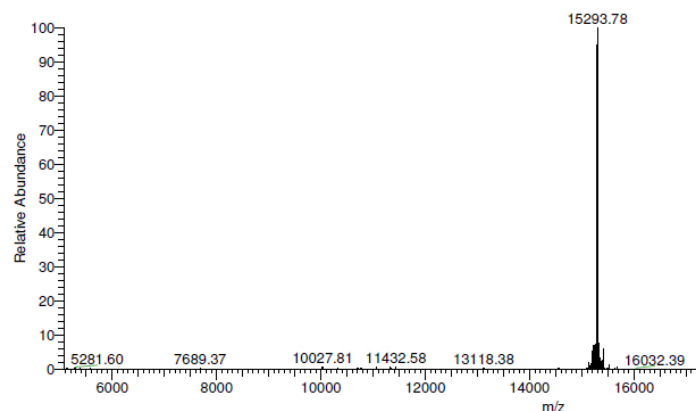
H3K9bu dNuc is highly purified and suitable for a variety of applications, including use as a substrate in enzymatic assays or for effector protein binding experiments.

References:

[1] Lowary PT and Widom J (1998) *J Mol Biol* 276:19-42.

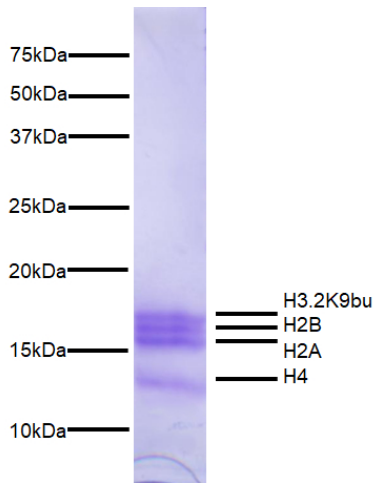


Western Blot Data: Western Analysis of H3K9bu dNuc. **Top Panel:** Unmodified H3 nucleosomes (Catalog No. 16-0006; Lane 1) and H3K9bu nucleosomes (Lane 2) were probed with an anti-H3K9bu antibody and analyzed via ECL readout. Only the H3K9bu sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified nucleosomes (Lane 1) and H3K9bu nucleosomes (Lane 2).

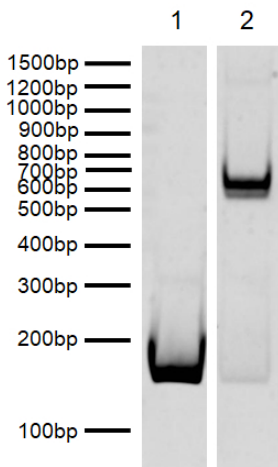


Mass Spec Data: Synthetic H3K9bu histone analyzed by high resolution mass spectrometry. Expected mass = 15,293.8 Da. Determined mass = 15,293.78 Da.

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



Protein Gel Data: Coomassie stained PAGE gel of proteins in H3K9bu dNuc (1 µg) to demonstrate the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3K9bu and H4) are indicated.



DNA Gel Data: H3K9bu dNuc resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (Catalog No 18-0005; 100 ng). **Lane 2:** Intact nucleosomes (400 ng).

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