

Nucleosome, Recombinant Human, H3K9ac dNuc, Biotinylated

Catalog No. 16-0314

Lot No. 21032002-15

Pack Size 50 µg



EpiCypher®

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 semi-synthetic method) contains acetyl-lysine at position 9. The nucleosome is the basic subunit of chromatin. The 601 sequence, identified by Lowary and Widom, is a 147-base pair sequence that has high affinity for histone octamers and is useful for nucleosome assembly and contains a 5' biotin-TEG group.

* H3K9ac also has a Cys to Ala substitution at position 110.

Formulation:

Nucleosome, Recombinant Human, H3K9ac (27.3 µg protein weight, 50 µg total weight) in 50 µL of 10 mM Tris HCl, pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molarity = 4.99 µmolar. MW = 199,847.66 Da.

Storage and Stability:

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

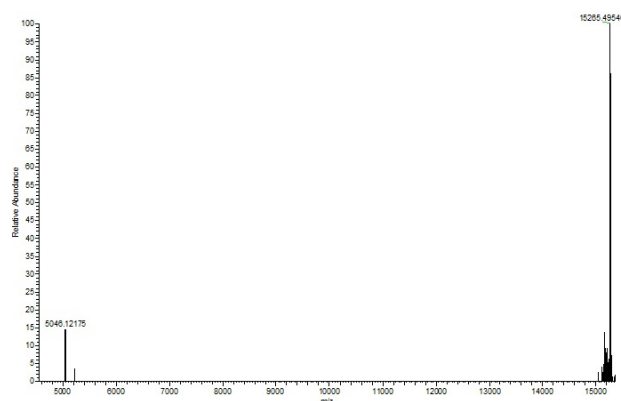
Application Notes:

Nucleosome, Recombinant Human, H3K9ac dNucs are highly purified and are suitable for use as substrates in enzyme screening assays or for effector protein binding experiments. **Nucleosome, Recombinant Human, H3K9ac dNucs from EpiCypher does not contain free DNA which could alter assayed activities.**

References:



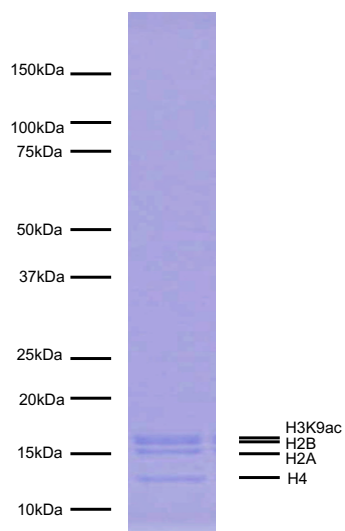
Western Blot Data: Western Analysis of Nucleosome, Recombinant Human, H3K9ac. **Top Panel:** Unmodified H3 (Lane 1) and H3K9ac containing nucleosomes (Lane 2) were probed with an anti-H3K9ac antibody and analyzed via ECL readout. Only the H3K9ac sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified nucleosomes (Lane 1) and H3K9ac nucleosomes (Lane 2).



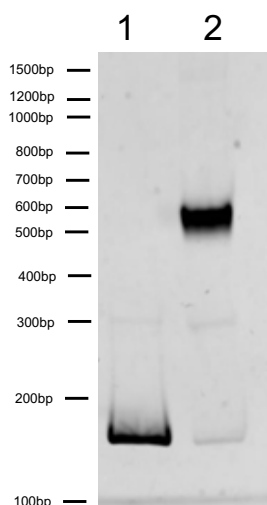
Mass Spec Data: H3K9ac protein analyzed by high resolution mass spectrometry. Expected mass = 15266.8 Da. Determined mass = 15,265.5 Da.

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



Protein Gel Data: Coomassie stained PAGE gel of proteins in Nucleosome, Recombinant Human, H3K9ac (1 µg) to demonstrate the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3K9ac and H4) are indicated.



DNA Gel Data: Nucleosome, Recombinant Human, H3K9ac run on a native PAGE gel and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (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.