

# Nucleosome, Recombinant Human, H3K4,K27me3 dNuc, Biotinylated

**Catalog No** 16-0403  
**Lot No** 22122001-05  
**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-P20671; 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 trimethylated-lysines at positions 4 and 27. 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.

\* H3K4,K27me3 has a Cys to Ala substitution at position 110.

## Formulation:

H3K4,K27me3 dNuc (27.2 µg protein weight, 50 µg DNA + protein) in 50.0 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molarity = 5.0 µM. MW = 199,935.2 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:

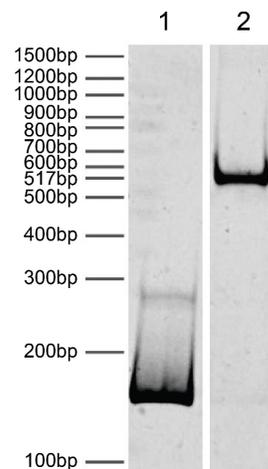
H3K4,K27me3 dNuc is highly purified and suitable for a variety of applications, including use as a substrate in enzymatic assays or for effector protein binding experiments.

## References:

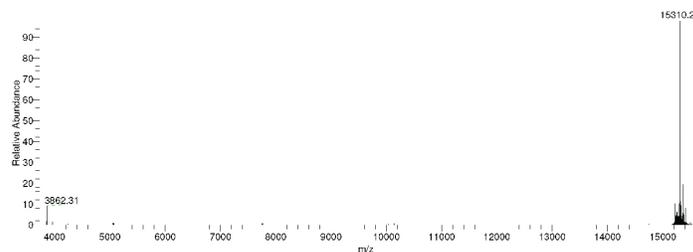
[1] Lowary & Widom *J Mol. Biol.* (1998) PMID: 9514715



# EpiCypher®

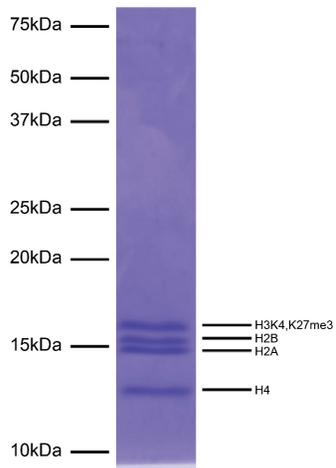


**Figure 1: DNA gel data.** H3K4,K27me3 dNuc resolved via native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (EpiCypher 18-0005; 100 ng). **Lane 2:** Intact H3K4,K27me3 nucleosomes (400 ng).



**Figure 2: Mass spec data.** Semi-synthetic H3K4,K27me3 histone analyzed by high resolution mass spectrometry. Expected mass = 15,308.8 Da. Determined mass = 15,310.20 Da.

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



**Figure 3: Protein gel data.** Coomassie stained PAGE gel of proteins in H3K4,K27me3 dNuc (1  $\mu$ g) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3K4,K27me3, and H4) are indicated.

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