

# Nucleosome, Recombinant Human, H4K20me1 dNuc, Biotinylated



## EpiCypher®

**Catalog No** 16-0331  
**Lot No** 21056002-05  
**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-P04908; H2B-O60814; H3.1-P68431; H4-P62805) wrapped by 147 base pairs of 601 positioning sequence DNA. Histone H4 (created by a proprietary synthetic method) contains monomethyl-lysine at position 20. 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.

### Formulation:

H4K20me1 dNuc (27 µg protein weight, 50 µg DNA + protein) in 51 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molarity = 4.9 µM. MW = 199,887.86 Da.

### Storage and Stability:

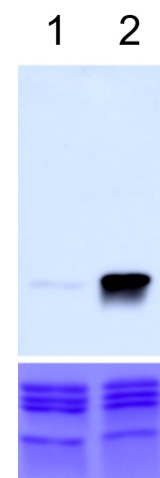
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

### Application Notes:

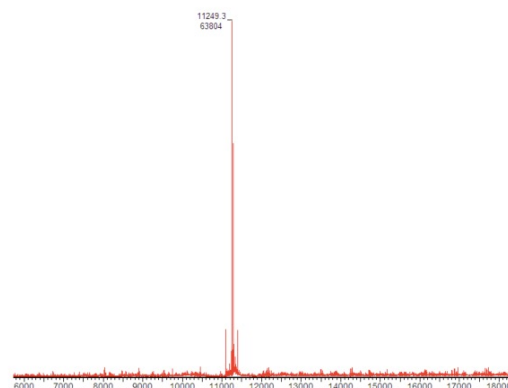
H4K20me1 dNuc is highly purified and suitable for a variety of applications, including use as a substrate in enzymatic assays or for effector protein binding experiments.

### References:

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

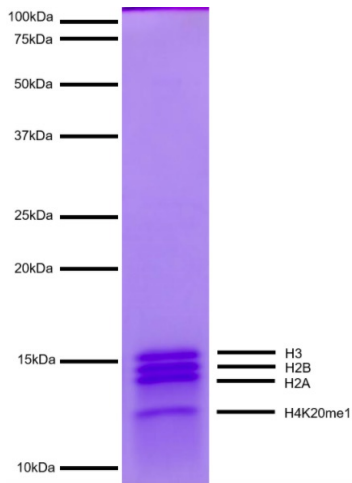


**Western Blot Data:** Western Analysis of H4K20me1 dNuc. **Top Panel:** Unmodified H4 nucleosomes (Catalog No. 16-0006; Lane 1) and H4K20me1 nucleosomes (Lane 2) were probed with an anti-H4K20me1 antibody and analyzed via ECL readout. Only the H4K20me1 sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified (Lane 1) and H4K20me1 nucleosomes (Lane 2).

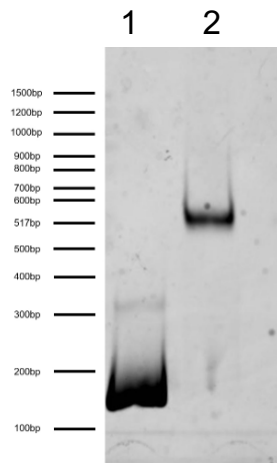


**Mass Spec Data:** Synthetic H4K20me1 histone analyzed by high resolution mass spectrometry. Expected mass = 11,250.1 Da. Determined mass = 11,249.3 Da.

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 H4K20me1 dNuc (2  $\mu$ g) to demonstrate the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4K20me1) are indicated.



**DNA Gel Data:** H4K20me1 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 H4K20me1 nucleosomes (400 ng).

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