

## Nucleosome, Recombinant Human, H3K4me1 dNuc, Biotinylated

<b>Catalog No</b>	16-0321	<b>Species</b>	Human
<b>Lot No</b>	17031002	<b>Source</b>	<i>E. coli</i> & synthetic DNA
<b>Pack Size</b>	50 µg	<b>Tag</b>	Biotinylated
<b>Concentration</b>	8.04 µM	<b>MW</b>	199,790 Da

### DESCRIPTION

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3\*, and H4) wrapped by 147 base pairs of 601 positioning sequence DNA. Histone H3 (created by a proprietary semi-synthetic method) contains monomethyl-lysine at position 4. 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. \*Histone H3.2 has a Cys to Ala substitution at position 110.

### TECHNICAL INFORMATION

<b>Storage</b>	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
<b>Formulation</b>	1.61 mg/mL mononucleosome in 31 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol (27.3 µg protein, 50 µg DNA + protein).

### APPLICATION NOTES

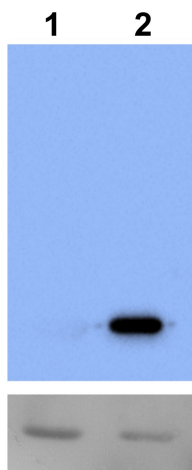
H3K4me1 dNuc is highly purified and suitable for a variety of applications, including use as a substrate in enzyme assays, high-throughput screening and inhibitor testing, chromatin binding studies, protein-protein interaction assays, structural studies, and in effector protein binding experiments. For a corresponding unmodified control, we recommend EpiCypher 16-0006.

### GENE & PROTEIN INFORMATION

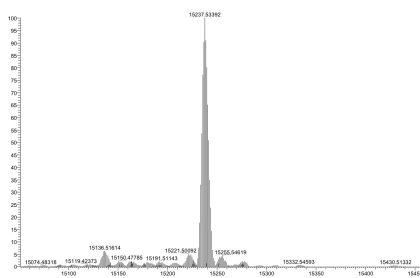
<b>UniProt ID</b>	H2A - P04908 (alt. names: H2A type 1-B/E, H2A.2, H2A/a, H2A/m) H2B - O60814 (alt. names: H2B K, HIRA-interacting protein 1) H3.2 - Q71DI3 H4 - P62805
-------------------	--

### REFERENCES

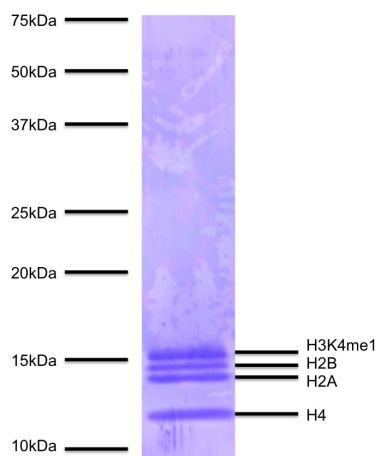
[1] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715



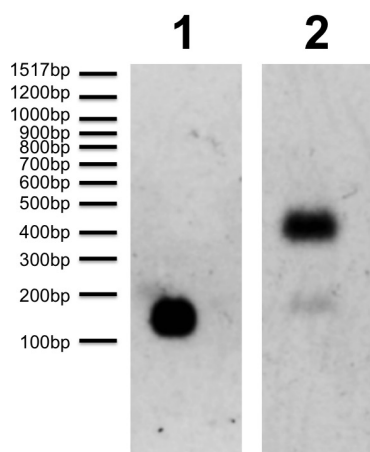
**FIGURE 1 Western blot data.** Western Analysis of H3K4me1 dNuc. **Top Panel:** Unmodified nucleosomes (EpiCypher 16-0006; Lane 1) and H3K4me1 nucleosomes (Lane 2) were probed with an anti-H3K4me1 antibody and analyzed via ECL readout. Only the H3K4me1 sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified octamer (Lane 1) and H3K4me1 octamer (Lane 2).



**FIGURE 2 Mass spec data.** Synthetic H3K4me1 histone analyzed by high resolution mass spectrometry. Expected mass = 15,238 Da. Determined mass = 15,237.5 Da.



**FIGURE 3 Protein gel data.** Coomassie stained PAGE gel of proteins in H3K4me1 dNuc (1 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3K4me1, and H4) are indicated.



**FIGURE 4 DNA gel data.** H3K4me1 dNuc resolved via agarose gel and stained with ethidium bromide to visualize DNA. Both lanes are from the same gel. **Lane 1:** Free DNA (EpiCypher 18-0005; 200 ng). **Lane 2:** Intact H3K4me1 nucleosomes (400 ng).