

H3K27me3 Recombinant Nucleosome, Biotinylated

Catalog No	16-0317-20	Species	Human
Lot No	25295006-01	Source	<i>E. coli</i> & synthetic DNA
Pack Size	20 µg	Tag	Biotinylated
Concentration	4.38 µM	MW	199,732 Da

DESCRIPTION

H3K27me3 (histone H3 lysine 27 trimethylation) Recombinant Nucleosome, Biotinylated consists of 147 base pairs of DNA wrapped around an octamer core of histone proteins (two each of H2A, H2B, H3.2, and H4) to form a nucleosome, the basic repeating unit 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. H3K27me3 nucleosome contains trimethylated lysine at position 27 and a Cys to Ala substitution at position 110 on histone H3.2. The DNA contains a 5' biotin-TEG group.

TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	0.875 mg/mL mononucleosome in 22.9 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol (10.9 µg protein, 20 µg DNA + protein).

APPLICATION NOTES

H3K27me3 nucleosome 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

UniProt ID	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

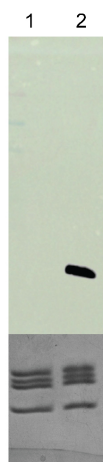


FIGURE 1 Western blot data. Western Analysis of H3K27me3 nucleosome. **Top Panel:** Unmodified (EpiCypher 16-0006; Lane 1) and H3K27me3 (Lane 2) nucleosomes were probed with an anti-H3K27me3 antibody and analyzed via enhanced chemiluminescence (ECL). Only the H3K27me3 sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified (Lane 1) and H3K27me3 (Lane 2) nucleosomes.

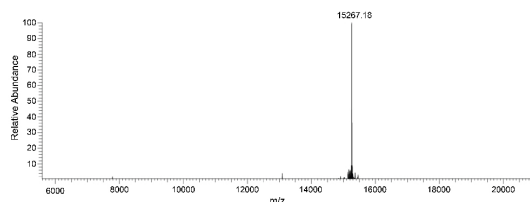


FIGURE 2 Mass spec data. Synthetic H3K27me3 histone analyzed by high resolution mass spectrometry. Expected mass = 15,266 Da. Determined mass = 15,267.18 Da.

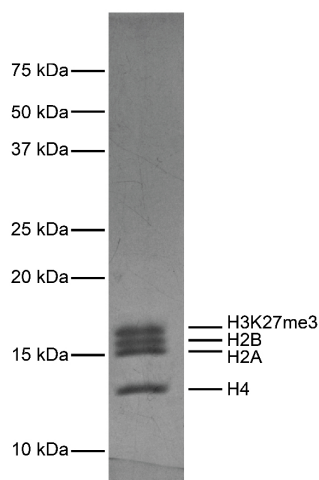


FIGURE 3 Protein gel data. Coomassie stained SDS-PAGE gel of proteins in H3K27me3 nucleosome (2 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3K27me3, and H4) are indicated.

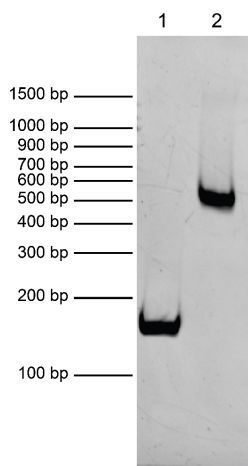


FIGURE 4 DNA gel data. H3K27me3 nucleosome resolved via native PAGE gel and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (EpiCypher 18-0005; 100 ng). **Lane 2:** Intact H3K27me3 nucleosomes (400 ng).