

Tailless Recombinant Nucleosome with Linker DNA, Biotinylated

Catalog No	16-2027	Species	Human
Lot No	23347001-01	Source	<i>E. coli</i> & synthetic DNA
Pack Size	50 µg	Tag	Biotinylated
Concentration	4.6 µM	MW	215,300 Da

DESCRIPTION

Recombinant mononucleosomes consist of 199 base pairs of DNA wrapped around an octamer core of histone proteins (two each of H2A, H2B, H3.1, and H4) to form a nucleosome, the basic repeating unit of chromatin. The 5' biotin-TEG DNA consists of a core 147 bp 601 nucleosome assembly sequence [1] flanked by 26 bp linker sequences as underlined below. After assembly, histone tails were enzymatically removed. This product is ideal for use as a negative control in binding assays

TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	0.99 mg/mL mononucleosome in 50.5 µL 10 mM Tris-HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. (21.4 µg protein, 50 µg DNA + protein).

APPLICATION NOTES

Tailless Recombinant Nucleosome with Linker DNA is highly purified and suitable for use as a negative control or substrate in enzyme screening assays and nucleosome binding experiments. The biotinylated DNA enables affinity binding applications.

DNA SEQUENCE

5'-Bio-TEG

GGACCCTATACGCGGCCGCCGAATTCCTGGAGAATCCCGGTCTGCAGGCCGCTCAATTGGTCGTAGACAGCTCTAGCACCG
CTTAAACGCACGTACGCGCTGTCCCCGCGTTTTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACGTGTCAGATATA
TACATCCTGTGGATCCGCCGGTTCGCGAACAGCGACC3'

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) H3.1 - P68431 (alt. names: H3, H3/a, H3/b, H3/c, H3/d) H4 - P62805
-------------------	---

REFERENCES

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

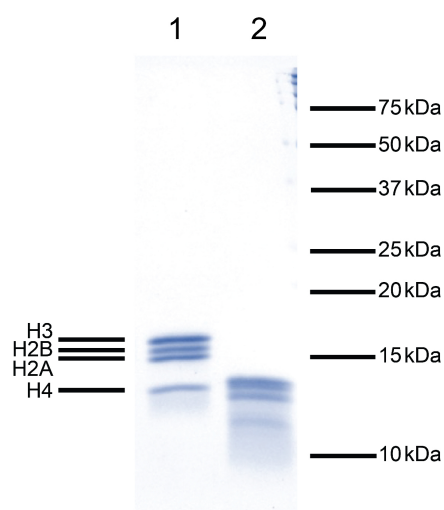


FIGURE 1 Protein gel data. Coomassie stained SDS-PAGE gel of proteins in Tailless Nucleosome (1 μ g) demonstrates the purity of histones in the preparation. **Lane 1:** Histone proteins in the nucleosome before enzymatic removal of tails. Sizes of molecular weight markers and positions of the intact core histones (H2A, H2B, H3, and H4) are indicated. **Lane 2:** Histone proteins in the nucleosome after tail removal via trypsin-digestion. Heterogeneity of the trypsin-treated histone proteins can be observed.

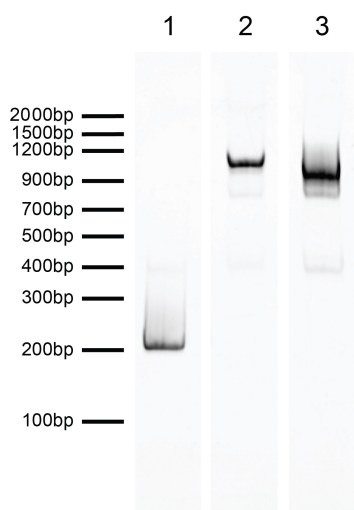


FIGURE 2 DNA gel data. Tailless Nucleosomes resolved via native PAGE and stained with ethidium bromide to visualize DNA. All lanes are from the same gel. **Lane 1:** Free DNA (EpiCypher 18-2044; 100 ng). **Lane 2:** Intact nucleosomes (400 ng). **Lane 3:** Nucleosomes after enzymatic removal of histone tails.