

Mononucleosomes, Recombinant, 177x601 DNA, Desthiobiotinylated

Catalog No. 16-3003

Lot No. 19038001

Pack Size 50 µg



EpiCypher®

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 177 base pairs of 601 positioning sequence DNA with a 5' desthiobiotin-TEG group. The nucleosome is the basic subunit of chromatin.

Formulation:

Mononucleosomes, Recombinant, 177x601 DNA (50 µg DNA+protein, 24.8 µg protein weight) in 36 µl 10 mM Tris pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM EDTA, 20% glycerol. Molarity = 6.37 µmolar. MW = 218,333.45 Da.

Storage and Stability:

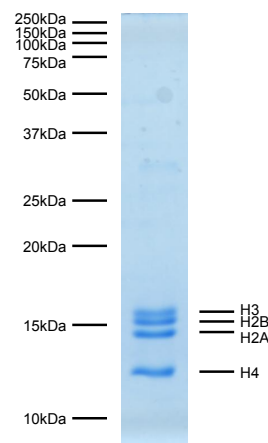
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

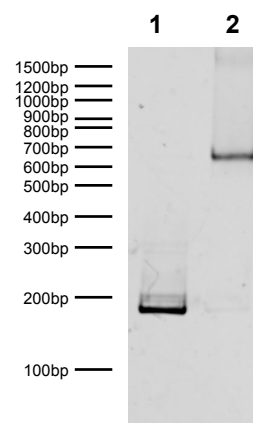
DNA sequence

5'Desthiobiotin-TEG-
CATCAGAATCCCGGTGCCGAGGCCGATCAATTGGTCGTAGAC
AGCTCTAGCACCGCTTAAACGCACGTACGCGCTGTCCCCGC
GTTTTAACCGCCAAGGGGATTACTCCCTAGTCTCCAGGCACG
TGTCAGATATATACATCGATGATGATGGATAGATGGATGATG
GATGGATGG-3'

5'CCATCCATCCATCATCCATCTATCCATCATCATCGATGTATA
TATCTGACACGTGCCTGGAGACTAGGGAGTAATCCCCTTGGC
GGTAAAACGCGGGGGACAGCGCGTACGTGCGTTTAAGCGG
TGCTAGAGCTGTCTACGACCAATTGATCGGCCTCGGCACCGG
GATTCTGATG-3'



Protein Gel Data: Coomassie stained PAGE gel of proteins in Mononucleosomes, Recombinant, 177x601, Desthiobiotinylated (1 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4) are indicated.



DNA Gel Data: Mononucleosomes, Recombinant, 177x601, Desthiobiotinylated resolved via native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (100 ng). **Lane 2:** Intact nucleosomes (400 ng).

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