

Mononucleosomes, Recombinant, Hemi-methylated



EpiCypher™

Catalog No. 16-2103
Lot No. 18058001
Pack Size 50 µg

Product Description:

Mononucleosomes assembled from recombinant histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4; accession numbers: H2A-P06897; H2B-P02281; H3-Q92133; H4-P62799) wrapped by 187 base pairs of DNA containing the 601 positioning sequence DNA. The 187 bp DNA sequence contains a 147 base-pair 601 nucleosome positioning sequence. The 601 sequence is flanked by a hemi-methylated 20 bp sequence as shown in application notes.

Formulation:

Mononucleosomes, Recombinant, 187x601 DNA (50 µg DNA + protein, 24.3 µg protein weight) in 52 µl 10 mM Tris pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM EDTA, 20% glycerol. Molarity = 4.32 µmolar. MW = 224,169 Da.

Storage and Stability:

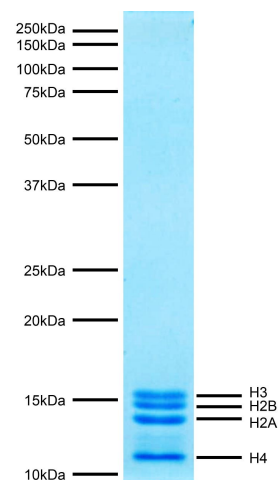
Stable for six months at -20°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

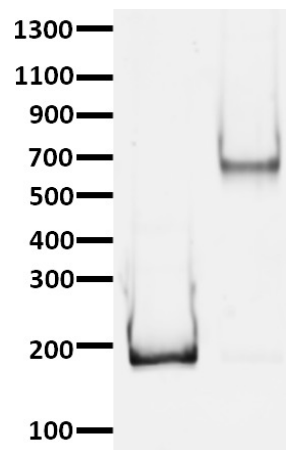
DNA sequence with methylation sites in **RED**

5'GGACCCTATAC**CGCGGCCG**CCCTGGAGAATCCCGGTCTGCA
GGCCGCTCAATTGGTCGTAGACAGCTCTACGTGGCGAATTTG
CGTGTCATGCGCCTGTCCCCGCGTTTTAACCGCCAAGGGGAT
TACTCCCTAGTCTCCAGGCACGTGTCAGATATATACATCCTGT
GCCGGTCGCGAACAGCGACC3'

3'**CCTGGGATATGCGCCGGCGG**GACCTCTTAGGGCCAGACGT
CCGGCGAGTTAACAGCATCTGTGCGAGATGCACCGCTTAAAC
GCACGTACGCGGACAGGGGGCGCAAATTGGCGGTTCCCT
AATGAGGGATCAGAGGTCCGTGCACAGTCTATATATGTAGG
ACACGGCCAGCGCTTGTGCTGG5'



Protein Gel Data: Coomassie stained PAGE gel of proteins in Mononucleosomes, Recombinant, 187x601 DNA (0.75 µg) to demonstrate 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, 187x601 DNA run on a native PAGE gel and stained with ethidium bromide to visualize DNA. Lane 1: Free DNA. Lane 2: Intact nucleosomes (200 ng).

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