

Mononucleosome, (H3.1ΔN32), 5' Cy5 167x601 DNA, Recombinant Human

Catalog No. 16-2026

Lot No. 19323001-15

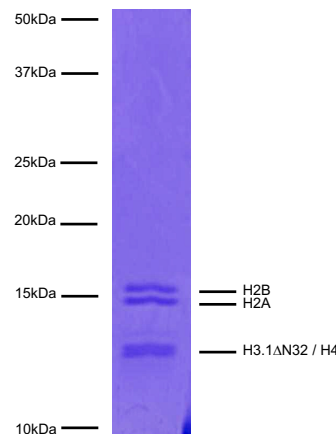
Pack Size 50 µg



EpiCypher®

Product Description:

Mononucleosomes assembled from recombinant 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) with the amino acid sequence of H3 beginning with glycine 33 (amino acids 1-32 are deleted) wrapped in a 167bp DNA sequence containing a 147bp Lowary and Widom 601 positioning sequence. The 601 sequence is flanked by a 10 bp sequence. The 601 DNA contains a 5' Cy5. The nucleosome is the basic subunit of chromatin.



Formulation:

Purified recombinant mononucleosomes (50 µg total mass, 24.9 µg protein and 25.1 µg DNA) in 45.5 µl 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT and 20% glycerol. Concentration of nucleosomes is 5.30 µM. Nucleosome molecular weight = 207,693 Da.

Protein Gel Data: Coomassie stained PAGE gel of proteins in Mononucleosome, (H3.1ΔN32), 5' Cy5 167x601 DNA, Recombinant Human (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.

Storage and Stability:

Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

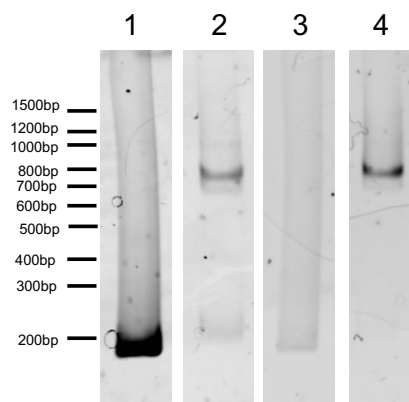
Application Notes:

DNA sequence

5'Cy5-

CGCGGCCGCCCTGGAGAATCCCGGTCTGCAGGCCGCTCAATT
GGTCGTAGACAGCTCTAGCACCGCTTAAACGCACGTACGCGC
TGCCCCCGGTTTTAACCGCCAAGGGGATTACTCCCTAGTCT
CCAGGCACGTGTCAGATATATACATCCTGTGCCGGTCGCG3'

3'GCGCCGGCGGGACCTCTTAGGGCCAGACGTCCGGCGAGTT
AACCAGCATCTGTCGAGATCGTGGCGAATTTGCGTGCATGCG
CGACAGGGGGCGAAAATTGGCGTTCCCCTAATGAGGGAT
CAGAGGTCGTGCACAGTCTATATATGTAGGACACGGCCAGC
GC5'



DNA Gel Data: Mononucleosomes, (H3.1ΔN32), 5' Cy5 167x601 DNA, Recombinant Human resolved via native PAGE and either stained with ethidium bromide and imaged (Lanes 1 and 2) or imaged at 630 nm light (Lanes 3 and 4) **Lane 1:** Free DNA (100 ng). **Lane 2:** Intact nucleosomes (400 ng). **Lane 3:** Free DNA (100 ng). **Lane 4:** Intact nucleosomes (400 ng).

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