

Mononucleosomes, (H3.1ΔN32), Recombinant Human, 5' Cy5

Catalog No. 16-0026
Lot No. 19277001-08
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), with the amino acid sequence of H3 beginning with glycine 33 (amino acids 1-32 are deleted), wrapped by 147 base pairs of 601 positioning sequence DNA with a 5' Cy5 fluor. The nucleosome is the basic subunit of chromatin.

Formulation:

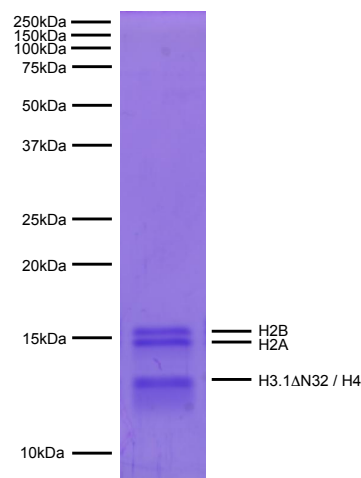
Purified recombinant mononucleosomes (50 µg total mass, 26.4 µg protein and 23.6 µg DNA) in 53.1 µL 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT and 20% glycerol. Concentration of nucleosomes is 4.87 µM. Nucleosome molecular weight = 193,270.96 Da.

Storage and Stability:

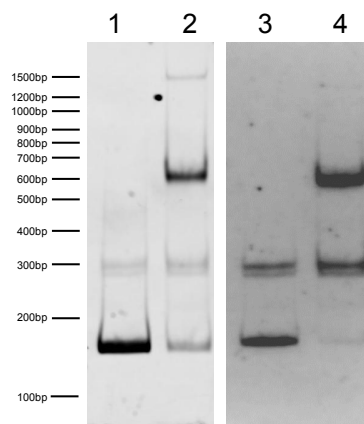
Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Mononucleosomes, (H3.1ΔN32), Human Recombinant, 5' Cy5 are highly purified and are suitable for use as substrates in enzyme screening assays or for nucleosome binding experiments. The absence of post-translational histone modifications makes this product ideal for conducting enzyme activity and screening assays.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Mononucleosomes, (H3.1ΔN32), Recombinant Human, 5' Cy5 (1 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3.1ΔN32, and H4) are indicated. H3.1ΔN32 and H4 co-migrate.



DNA Gel Data: Mononucleosomes, (H3.1ΔN32), Recombinant Human, 5' Cy5 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.