

Mononucleosomes (H2AZ.1), Human Recombinant, Non-biotinylated

Catalog No. 16-1014

Lot No. 20195003-40

Pack Size 50 µg



EpiCypher®

Product Description:

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2AZ.1, H2B, H3 and H4. Accession numbers: H2AZ.1-POC0S5; H2B-O60814; H3.1-P68431; H4-P62805) wrapped by 147 base pairs of 601 positioning sequence DNA. The nucleosome is the basic subunit of chromatin. H2AZ.1 and H2AZ.2 are variants of histone H2A and implicated in diverse cellular functions ranging from transcriptional regulation, chromosome transmission and DNA damage repair. H2AZ has a dedicated deposition machinery (the SWR-C ATPase).

Formulation:

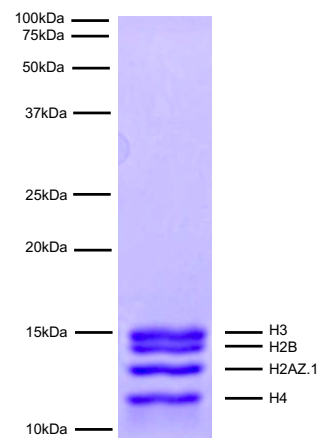
Purified recombinant mononucleosomes (27.4 µg protein weight, 50.0 µg DNA+protein) in 43.8 µl in 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT, & 20% glycerol. Concentration of nucleosomes is 5.04 µM. Nucleosome molecular weight = 198,425.5 Da.

Storage and Stability:

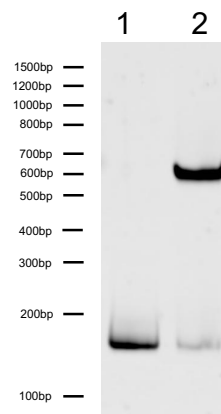
Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Mononucleosomes (H2AZ.1), Human Recombinant Biotinylated 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 them ideal for conducting enzyme activity and screening assays.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Nucleosome (H2AZ.1), Human Recombinant Biotinylated (1 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2AZ.1, H2B, H3 and H4) are indicated.



DNA Gel Data: Nucleosome (H2AZ.1), Human Recombinant Biotinylated resolved via native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** DNA extracted from nucleosomes (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.