

# Mononucleosomes (H3.3K27M), Recombinant Human, Biotinylated



EpiCypher®

**Catalog No** 16-0323  
**Lot No** 21214004-07  
**Pack Size** 100 µg

## Product Description:

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3.3 and H4; accession numbers: H2A-P04908; H2B-O60814; H3.3-P84243; H4-P62805) wrapped by 147 base pairs of 601 positioning sequence DNA. Histone H3.3 is a non-allelic replacement found in regions of high chromatin turnover outside of S-phase (e.g. at actively transcribed genes). The substitution of M for K at position 27 is a dominant negative inhibitor of the EZH2 methyltransferase complex PRC2 and is associated with pediatric brain cancer (DIPG) [1]. The nucleosome is the basic subunit of chromatin. The 147 bp 601 sequence, identified by Lowary and Widom [2], has high affinity for histone octamers and is useful for nucleosome assembly. The DNA contains a 5' biotin-TEG group.

## Formulation:

H3.3K27M Mononucleosomes (54.6 µg protein weight, 100.0 µg DNA + protein) in 90.9 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molarity = 5.5 µM. MW = 199,844 Da.

## Storage and Stability:

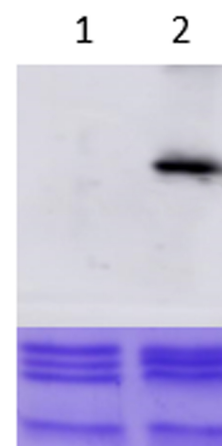
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

## Application Notes:

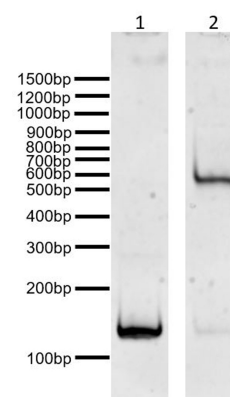
H3.3K27M Mononucleosomes is highly purified and suitable for a variety of applications, including use as a substrate in enzymatic assays or for effector protein binding experiments.

## References:

- [1] Chan et al. (2013) *Genes Dev* 27:985-990.
- [2] Lowary PT and Widom J (1998) *J Mol Biol* 276:19-42.

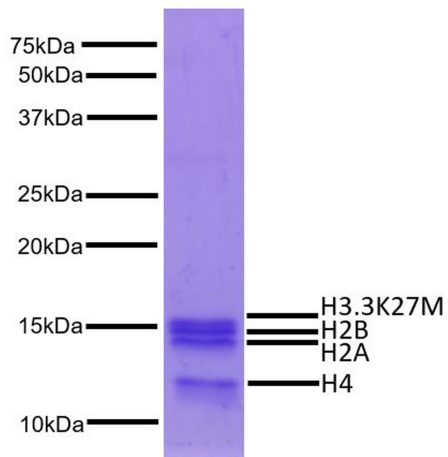


**Western Blot Data:** Western analysis of Mononucleosomes (H3.3K27M), Recombinant Human. **Top Panel:** Wild-type H3.3 (Lane 1) and H3.3K27M nucleosomes (Lane 2) were probed with an H3K27M antibody. Only the H3.3K27M sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel of Western blot.



**DNA Gel Data:** H3.3K27M Mononucleosomes resolved via Native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** Free 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.



**Protein Gel Data:** Coomassie stained PAGE gel of proteins in H3.3K27M Mononucleosomes (1  $\mu$ g) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3.3K27M and H4) are indicated.

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