

## Recombinant Mononucleosomes H3.3 (K27M) - biotin

---

**Catalog No:** 81006, 81706

**Lot No:** 17717001

**Expressed In:** *E. coli*

**Quantity:** 20, 1000 µg

**Concentration:** 0.75 µg/µl

**Source:** Human

**Buffer Contents:** Recombinant Mononucleosomes H3.3 (K27M) - biotinylated (20 µg protein + 20 µg DNA) is supplied at a protein concentration of 0.75 µg/µl in 10 mM Tris-HCl, pH 8.0, 1 mM EDTA, 2 mM DTT and 20% glycerol.

**Background:** *In vivo*, the nucleosome is the basic structural unit of chromatin. It consists of about 146 bp of DNA wrapped around a core of eight histones of four different types: H2A, H2B, H3 and H4. Histones are subject to posttranslational modifications, such as methylation, acetylation, phosphorylation, mono-ubiquitination, etc. Histone modifications influence multiple chromatin templated processes such as gene transcription, DNA repair and recombination. Besides the “major” histones, there are some histone variants in specific regions of chromatin or in specific cell types. Histone variants were involved in multiple biology processes including chromosome segregation, DNA repair, transcriptional regulation and mRNA processing. Histone H3.3 point mutations (K27 and G34) are found in 1/3 of pediatric glioblastomas. Up to 78% of diffuse intrinsic pontine gliomas (DIPGs) carry K27M and 36% of non-brainstem gliomas carry either K27M or G34R/V mutations. High-frequency mutation of histone H3 to K36M is seen in chondroblastomas and mutation to G34W/L in giant cell tumors of bone, which are diseases of adolescents and young adults. Histone H3.3 mutations may drive pediatric glioblastoma through upregulation of MYCN.

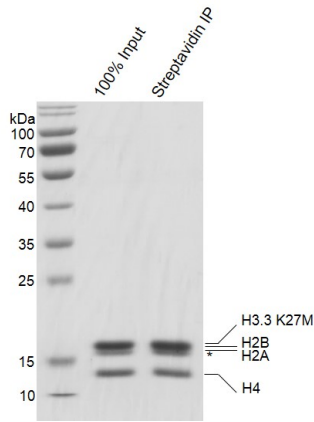
Nucleosomes are more physiologically relevant substrates than histones and histone-derived peptides for *in vitro* studies. More importantly, some histone methyltransferases are significantly more active, as well as specific, when using nucleosomal substrates in HMT assays, such as DOT1L and NSD family enzymes. Nucleosomes are also widely used in histone methyltransferase screening assays to identify small molecular inhibitors for drug discovery.

**Protein Details:** Recombinant Mononucleosomes H3.3 (K27M) – biotinylated, consist of a 167 bp of 601 DNA with 5' biotin tag and two molecules each of histones H2A that includes amino acids 1-130 (end) (accession number NP\_003503.1), H2B that includes amino acids 1-126 (end) (accession number NP\_003509.1), H3.3 that includes amino acids 1-136 (end) (accession number NP\_002098.1) with a point mutation Lys27Met, and H4 that includes amino acids 1-103 (end) (accession number NP\_003539.1). All of these histones were expressed in *E. coli* cells. The molecular weight of histone octamer is ~108 kDa. The recombinant protein is ≥95% by SDS-PAGE gel.

**Application Notes:** Recombinant Mononucleosomes H3.3 (K27M) - biotinylated are suitable for use in the study of enzyme kinetics, inhibitor screening, and selectivity profiling.

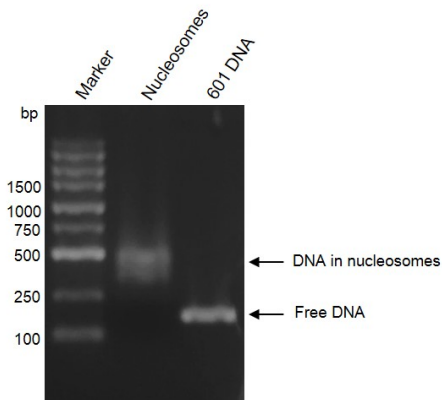
**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

### Streptavidin pull-down for Recombinant Mononucleosomes H3.3 (K27M) - biotin.



Recombinant Mononucleosomes H3.3 (K27M) - biotinylated were pulled down by streptavidin beads. Input nucleosomes (Lane 2) and the nucleosomes (Lane 3) pulled down by streptavidin were run on a 13% SDS-PAGE gel and stained with Commassie blue. The SDS-PAGE gel result shows that almost all of biotinylated mononucleosomes H3.3 (K27M) are pulled down by streptavidin beads. \* indicates streptavidin.

### Recombinant Mononucleosomes H3.3 (K27M) - biotin, DNA gel.



Recombinant Mononucleosomes H3.3 (K27M) - biotinylated were run on a 2% agarose gel and stained with ethidium bromide. Lane 1: DNA marker. Lane 2: Intact mononucleosomes H3.3 (K27M). Lane 3: 601 DNA. Intact mononucleosomes H3.3 (K27M) migrated much higher than free 601 DNA. The agarose gel result shows almost all of 601 DNA wrap histone octamers to form nucleosomes.

### Western Blot analysis for Recombinant Mononucleosomes H3.3 (K27M).

PRC2 Complex:	0	0	1	1	:ug
Nucleosomes H3.3 (WT):	1	0	1	0	
Nucleosomes H3.3 (K27M):	0	1	0	1	



1  $\mu$ g PRC2 Complex (Cat. No. 31387) was incubated with 2  $\mu$ g wild type Mononucleosomes H3.3 or K27M mutant Mononucleosomes H3.3, respectively, in 30  $\mu$ l reaction system containing 50 mM Tris-HCl, pH 8.6, 0.02% Triton X-100, 2 mM MgCl<sub>2</sub>, 1 mM TCEP, 50  $\mu$ M SAM for 3 h at room temperature. 6  $\mu$ l reaction products were loaded and run on a 12.5% SDS-PAGE gel. Western blot was used to detect the generation of reaction products (Anti-H3K27me3, Cat. No. 39155). The Western blot result shows that K27M mutation inhibits the activity of PRC2 Complex.