

Nucleosome, Recombinant Human, Acidic Patch Mutant H2AE92K Biotinylated

Catalog No	16-0030	Species	Human
Lot No	22224002-01	Source	<i>E. coli</i> & synthetic DNA
Pack Size	50 µg	Tag	Biotinylated
Concentration	4.9 µM	MW	199,741 Da

DESCRIPTION

Recombinant mononucleosomes consist of 147 base pairs of DNA wrapped around an octamer core of histone proteins (two each of H2A, H2B, H3.1, and H4) to form a nucleosome, the basic repeating unit of chromatin. The 147 bp 601 sequence containing a 5' biotin-TEG group has high affinity for histone octamers and is useful for nucleosome assembly [1]. Histone H2A contains a glutamate-to-lysine (E-to-K) substitution at position 92 (H2AE92K). H2AE92 is among key residues forming a negatively charged region on the nucleosome surface named the "acidic patch". The acidic patch is a conserved interaction hub for neighboring nucleosomes and nucleosome binding proteins, often via salt bridges with arginine anchors, and is functionally critical in chromatin condensation and chromatin remodeling [2-4]. H2AE92 resides in the H2A C-terminal extension and is associated with nucleosome binding factors such as histone H4 N-terminal tail, LANA, RCC1, IL-33, SIR3, HMGN2, and remodeling ATPases [2-4]. H2AE92K disrupts binding with SMARCB1 [5].

TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	0.98 mg/mL mononucleosome in 51.1 µL 10 mM Tris-HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. (27.3 µg protein, 50 µg DNA + protein)

APPLICATION NOTES

H2AE92K mononucleosome is highly purified and suitable for a variety of applications to test the effect of acidic patch mutation on enzymatic activity or chromatin binding. The biotinylated DNA enables affinity binding applications.

GENE & PROTEIN INFORMATION

UniProt ID	H2A - P04908 (alt. names: H2A type 1-B/E, H2A.2, H2A/a, H2A/m) H2B - O60814 (alt. names: H2B K, HIRA-interacting protein 1) H3.1 - P68431 (alt. names: H3, H3/a, H3/b, H3/c, H3/d) H4 - P62805
-------------------	---

REFERENCES

- [1] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715
- [2] Kalashnikova et al. *Soc. Interface* (2013). PMID: 23446052
- [3] Levendosky & Bowman *eLife* (2019). PMID: 31094676
- [4] Gamarra et al. *eLife* (2018). PMID: 29664398
- [5] Valencia et al. *Cell* (2019). PMID: 31759698

VALIDATION DATA

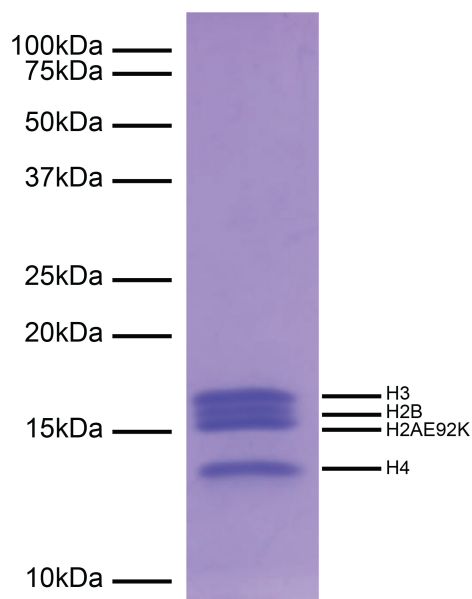


FIGURE 1 Protein gel data. Coomassie stained PAGE gel of proteins in H2AE92K mononucleosome (1 μ g) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2AE92K, H2B, H3 and H4) are indicated.

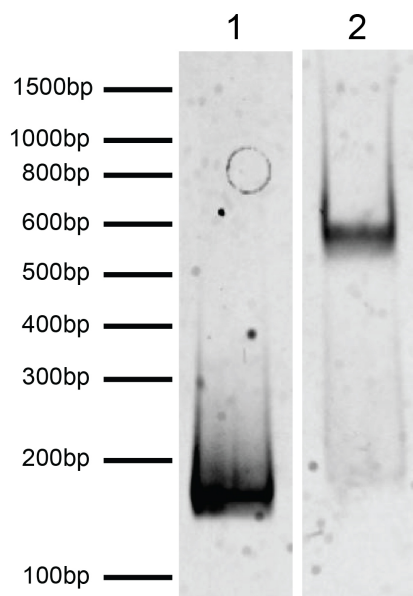


FIGURE 2 DNA gel data. H2AE92K mononucleosome resolved by native PAGE and stained with ethidium bromide to visualize DNA. **Lane 1:** Free DNA (EpiCypher 18-0005; 100 ng). **Lane 2:** Intact H2AE92K mononucleosomes (400 ng).

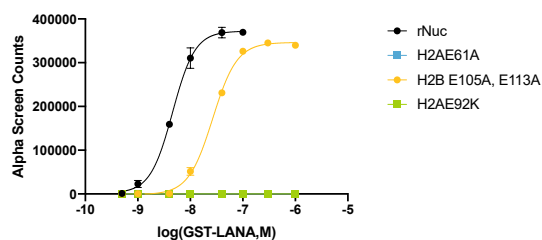


FIGURE 3 Functional binding assay. The presence of acidic patch mutations disrupts LANA peptide binding to recombinant nucleosomes (WT control, EpiCypher 16-0006; H2AE61A, EpiCypher 16-0029; H2AE92K, EpiCypher 16-0030; H2BE105A, E113A, EpiCypher 16-0031). The binding of GST-tagged LANA peptide to biotinylated recombinant nucleosomes was assessed by AlphaLISA assay using Streptavidin Donor Beads and Glutathione Acceptor Beads (PerkinElmer). The presence of H2A acidic patch mutations completely blocks LANA binding, while H2B mutations cause a decrease in LANA binding affinity.