

## H3S28ar1 Recombinant Nucleosome, Biotinylated

<b>Catalog No</b>	16-0420	<b>Species</b>	Human
<b>Lot No</b>	25038001-02	<b>Source</b>	<i>E. coli</i> & synthetic DNA
<b>Pack Size</b>	25 µg	<b>Tag</b>	Biotinylated
<b>Concentration</b>	4.58 µM	<b>MW</b>	200,728.3 Da

### DESCRIPTION

ADP-ribosylation (ADPr) is a reversible histone modification that regulates chromatin structure, transcription, DNA replication, and DNA damage repair [1]. Occurring as either mono- or poly-ADP-ribose, ADPr supports recruitment of DNA repair proteins and chromatin remodelers [1]. Recombinant ADPr nucleosomes provide a defined platform for studying biological, chromatin-based functions of ADP-ribosylation.

#### Product Highlights:

- **Physiologically relevant binding** – Nucleosomes provide a physiologically relevant substrate for studying chromatin biology, capturing the multivalent binding environment that many chromatin-associated proteins require.
- **Built for versatility** – Designed for broad biochemical assays, these nucleosome substrates enable robust, quantitative analysis of protein binding, enzymatic activity, and chromatin interactions in a defined relevant format.
- **Research ready performance** – Built on deep nucleosome design expertise: EpiCypher's designer nucleosomes are supported by > 1,000 publications across diverse research applications including immunology, cancer, and molecular biology.

H3S28ar1 (histone H3 serine 28 mono-ADPr) Recombinant Nucleosome, Biotinylated consists of 147 base pairs of 601 sequence DNA [2] wrapped around an octamer of core histone proteins (two each of H2A, H2B, H3.2, and H4) to form a nucleosome, the basic repeating unit of chromatin. The DNA contains a 5' biotin-TEG group. H3S28ar1 nucleosome contains H3.2 with mono-ADP-ribosylated serine at position 28. Histone H3.2 also contains a Cys to Ala substitution at position 110. H3S28ar1 is ADP-ribosylated by poly-ADPr polymerase 1 (PARP1) and PARP2, and ADPr removal is achieved by the glycohydrolase ARH3 [1].

### TECHNICAL INFORMATION

<b>Storage</b>	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
<b>Formulation</b>	0.92 mg/mL mononucleosome in 27.2 µL 10 mM Tris HCl pH 7.5, 25 mM NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol (13.6 µg protein, 25 µg DNA + protein).

### APPLICATION NOTES

H3S28ar1 nucleosome is highly purified and suitable for a variety of applications, including use as a substrate in enzyme assays, high-throughput screening and inhibitor testing, chromatin binding studies, protein-protein interaction assays, structural studies, and in effector protein binding experiments. For a corresponding unmodified control, we recommend EpiCypher 16-0006.

### GENE & PROTEIN INFORMATION

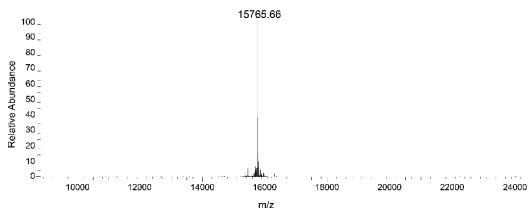
<b>UniProt ID</b>	H2A - P04908 (alt. names: H2A type 1-B/E, H2A.2, H2A/a, H2A/m) H2B - O60814 (alt. names: H2B K, HIRA-interacting protein 1) H3.2 - Q71DI3 H4 - P62805
-------------------	--

### REFERENCES

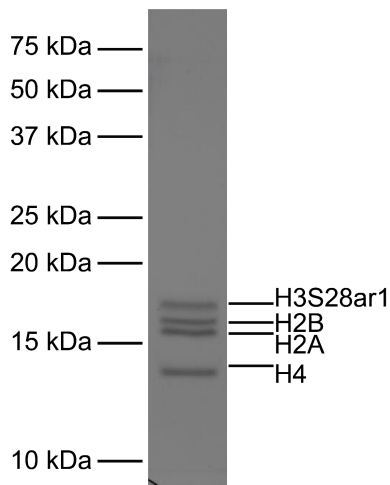
- [1] Longarini & Matic *DNA Repair*. (2022). PMID: 35963141  
[2] Lowary & Widom *J. Mol. Biol.* (1998). PMID: 9514715



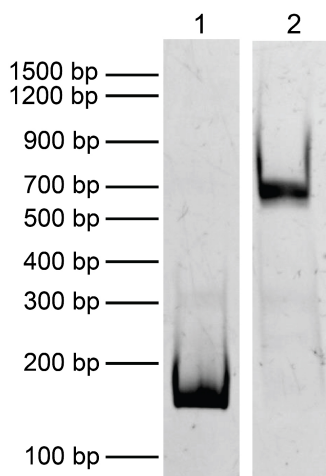
**FIGURE 1 Western blot data.** Western analysis of H3S28ar1 nucleosome. All lanes were resolved on a single gel. **Top Panel:** Unmodified (EpiCypher 16-0006; Lane 1) and H3S28ar1 (Lane 2) nucleosomes were probed with an anti-Poly/Mono-ADP Ribose antibody and analyzed via enhanced chemiluminescence (ECL). Only the H3S28ar1 sample produced a detectable signal. **Bottom Panel:** Detail from Coomassie stained gel showing unmodified (Lane 1) and H3S28ar1 (Lane 2) nucleosomes.



**FIGURE 2 Mass spec data.** H3S28ar1 histone analyzed by high resolution mass spectrometry. Expected mass = 15765.88 Da. Determined mass = 15765.66 Da.



**FIGURE 3 Protein gel data.** Coomassie stained SDS-PAGE gel of proteins in H3S28ar1 nucleosome (2  $\mu$ g) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3S28ar1, and H4) are indicated.



**FIGURE 4 DNA gel data.** H3S28ar1 nucleosome resolved via native PAGE and stained with ethidium bromide to visualize DNA. All lanes were resolved on a single gel. **Lane 1:** Free DNA (EpiCypher 18-0005; 100 ng). **Lane 2:** Intact H3S28ar1 nucleosomes (400 ng).