

## Histone Octamer, Recombinant Human

Catalog No	16-0001	Species	Human
Lot No	24342016-03	Source	<i>E. coli</i>
Pack Size	50 µg	Tag	None
Concentration	9.9 µM	MW	108,545 Da

### DESCRIPTION

Human histone octamers (two each of histones H2A, H2B, H3 and H4) are made from recombinant histones expressed in *E. coli*. Octamers are the protein component of nucleosomes, the basic subunit of chromatin. Assembling this product with DNA bearing a 601 Widom sequence (e.g. EpiCypher 18-0005, 18-0006) produces a nucleosome. The 601 Widom sequence, originally discovered by Lowary and Widom [1], has high affinity for octamers.

### TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2 M NaCl, 2 mM DTT, 20% glycerol

### APPLICATION NOTES

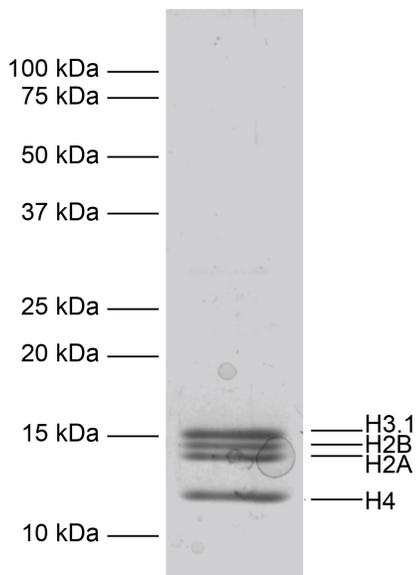
Human recombinant histone octamers can be used for chromatin reconstitution experiments or as substrates for histone modifying enzymes. See Luger et al. [2] for recommended nucleosome reconstitution protocol. Histone octamers should not be stored under conditions where the salt concentration is lower than 800 mM. **Please see lot-specific Technical Datasheets for product concentration.**

### 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] Luger et al. *Methods Mol. Biol.* (1999). PMID: 10804500



**FIGURE 1 Protein gel data.** Histone Octamer, Recombinant Human (1  $\mu$ g) run on an SDS-PAGE gel and stained with Coomassie blue to demonstrate the purity of the preparation. The individual histones are indicated.