

Histone Octamer, Recombinant Human



EpiCypher®

Catalog No. 16-0001

Lot No. 21084013-04

Pack Size 50 µg

Product Description:

Human histone octamers (two each of histones H2A, H2B, H3 and H4) made from recombinant histones expressed in *E. coli* (accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805). The histone octamer is the protein component of the nucleosome, the basic subunit of chromatin. A nucleosome consists of a histone octamer wrapped with 147 base pairs of DNA (note that there is only histone octamer in this product).

Formulation:

Gel filtration-purified recombinant human histone octamers (0.735 mg/mL) in 68.0 µL of 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2 M NaCl, 2 mM DTT, 20% glycerol. MW: 109,536 Da. Molarity: 6.71 µM.

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:

Human recombinant histone octamers can be used for chromatin reconstitution experiments, or as substrates for histone modifying enzymes. See Luger et al. (1999) for recommended nucleosome reconstitution protocol. Histone octamers should not be stored under conditions where the salt concentration is lower than 800 mM.

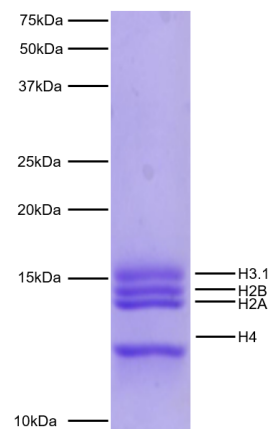
Please note that the concentration may vary compared to previous lots.

References:

Luger et al (1999). *Methods Enzymol.* 304:3-19.

References using this product:

Cao et al (2013). *Epigenetics* 8:477-485.



Protein Gel Data: Histone Octamer, Recombinant Human (1 µg) run on a PAGE gel and stained with Coomassie blue to demonstrate the purity of the preparation. The individual histones are indicated.

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