

# Histone Octamer, Recombinant Human

**Catalog No.** 16-0001-0.5 mg

**Lot No.** 14283003

**Pack Size** 500 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; H3.2-Q71DI3; 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 (octamer only in this product).

## Formulation:

Gel filtration-purified recombinant human histone octamers (0.56 mg/ml) in 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2 M NaCl, 2 mM DTT, 50% glycerol.

## Storage and Stability:

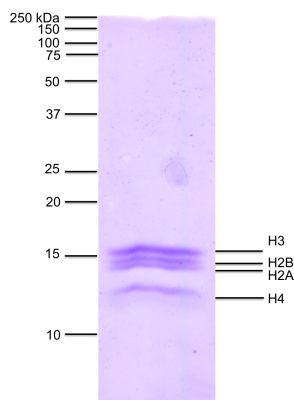
Stable for six months at -20°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

## Application Notes:

Human recombinant histone octamers can be used for chromatin reconstitution experiments, or as substrates for histone modifying enzymes. See product page on web for links to protocols for chromatin reconstitution. Histone octamers should not be stored under conditions where the salt concentration is lower than 800 mM.

## References:

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



**Protein Gel Data:** Histone Octamer, Recombinant Human (2 µg) run on a PAGE gel and stained with Coomassie blue to demonstrate the purity of the preparation. The individual histones are indicated.