

# Histone Octamer, Recombinant Xenopus

**Catalog No.** 16-0007  
**Lot No.** 15036001  
**Pack Size** 1 mg

## Product Description:

Recombinant histone octamers (two each of histones H2A, H2B, H3 and H4) made from recombinant *Xenopus* histones expressed in *E. coli* (accession numbers: H2A-P06897; H2B-P02281; H3-Q92133; H4-P62799). 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.

## Formulation:

Gel filtration-purified recombinant histone octamers (1.0 mg/ml) in 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2 M NaCl, 2 mM DTT, 20% glycerol.

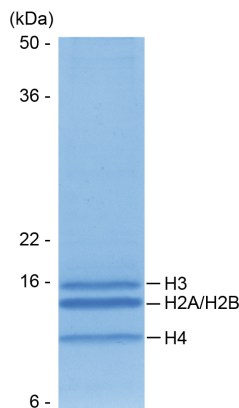
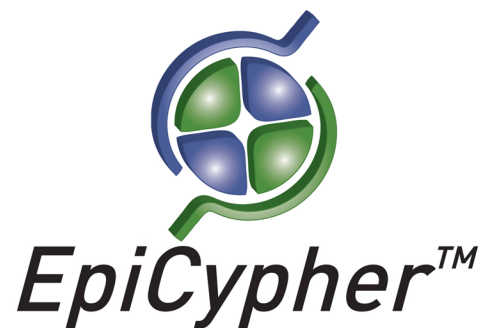
## Storage and Stability:

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

## Application Notes:

Histone Octamer, Recombinant Xenopus 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:



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