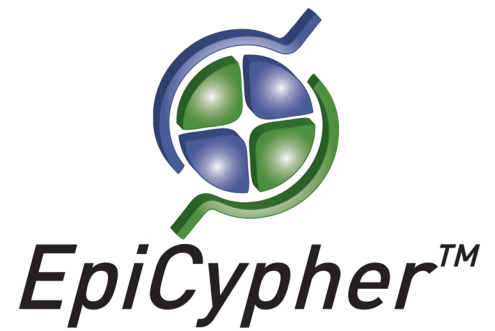


# Histone CENP-A/H4 Tetramer, Recombinant Human

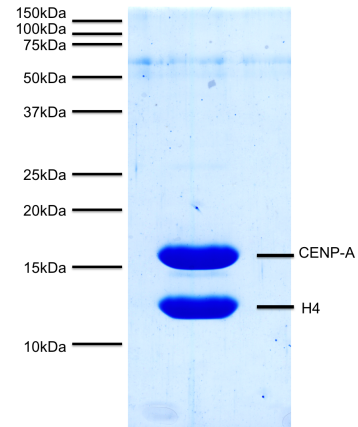
**Catalog No.** 16-0010  
**Lot No.** 16028001  
**Pack Size** 50 µg



**Type** Histone                      **Expressed In** *E. coli*  
**Mol. Wgt.** 54 kDa                      **Epitope Tag** None

## Product Description:

CENP-A/ Histone H4 Tetramers made from recombinant histones expressed in *E. coli* (Centromere protein A, CENPA, CenH3; accession numbers: CENP-A: P49450; H4: P62805). CENP-A and H4 were expressed and purified individually by FPLC, then assembled into tetramers that were further purified using gel filtration chromatography. CENP-A is a specialized variant of H3 and component of modified nucleosomes found at the core centromere.



**Protein Gel Data:** Histone CENP-A/H4 Tetramer, Recombinant Human (1 µg) was run on a PAGE gel and stained with Coomassie blue. The migration and molecular weight of the protein standards and migration of CENP-A and H4 are indicated.

## Formulation:

50 µg of CENP-A/H4 Tetramer (2 mg/ml) in 40mM sodium phosphate buffer, pH 7.2, 1M NaCl, 7mM β-ME, 50%

## 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:

Recombinant histone CENP-A/H4 Tetramer is suitable for enzyme assays and nucleosome reconstitution.

## References:

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