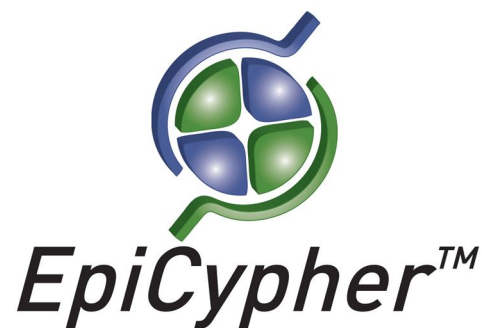


# Histone H3/H4 Tetramer, Recombinant Human



**Catalog No.** 16-0008  
**Lot No.** 17020002  
**Pack Size** 50 µg

## Product Description:

Histone H3/H4 Tetramers made from recombinant histones expressed in *E. coli* (accession numbers: H3-P68431; H4-P62805). Histones H3 and H4 were expressed and purified individually by FPLC, then assembled into tetramers that were further purified using gel filtration chromatography.

## Formulation:

50 µg of H3/H4 Tetramer (0.61 mg/ml) in 10mM Tris-HCl pH 7.5, 2M NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol. Molec. weight: 53,210 Da. Concentration: 11.5 µM.

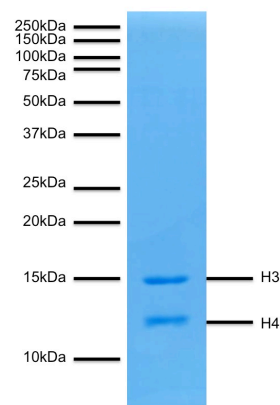
## Storage and Stability:

Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

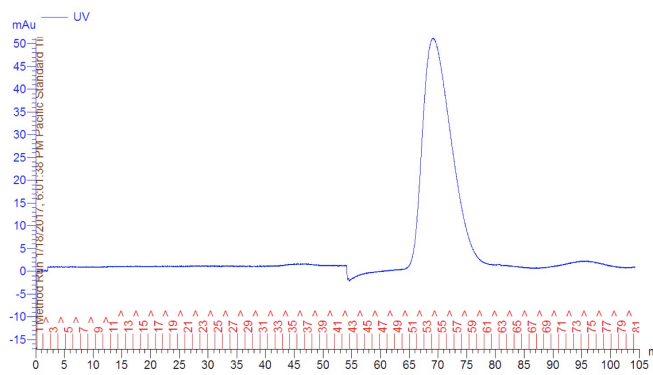
## Application Notes:

Recombinant histone H3/H4 Tetramer is suitable for enzyme assays and nucleosome reconstitution.

## References:



**Protein Gel Data:** Histone H3/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 H3 and H4 are indicated.



**Purification Data:** Chromatogram from gel filtration purification of Histone H3/H4 Tetramer, Recombinant Human. Only fractions corresponding to the intact tetramer (as indicated, black bar) were collected and pooled.