

## Histone H3/H4 Tetramer, Recombinant Human

<b>Catalog No</b>	16-0008	<b>Species</b>	Human
<b>Lot No</b>	19228003	<b>Source</b>	<i>E. coli</i>
<b>Pack Size</b>	50 µg	<b>Tag</b>	None
<b>Concentration</b>	25.6 µM	<b>MW</b>	53,210 Da

### DESCRIPTION

Histone H3/H4 Tetramers made from recombinant histones expressed in *E. coli*. Histones H3 and H4 were expressed and purified individually, then assembled into tetramers that were further purified using gel filtration chromatography.

### TECHNICAL INFORMATION

<b>Storage</b>	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
<b>Formulation</b>	1.36 mg/mL H3/H4 Tetramer in 36.8 µL 10 mM Tris-HCl pH 7.5, 2M NaCl, 1 mM EDTA, 2 mM DTT, 20% glycerol.

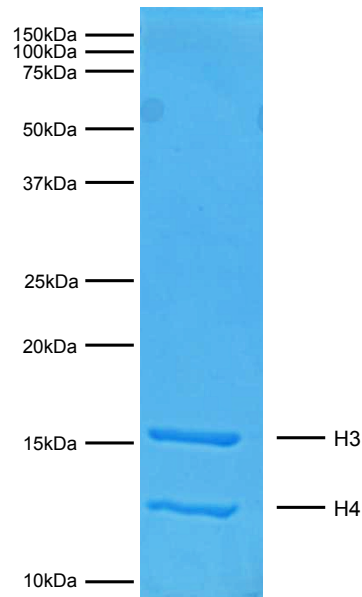
### APPLICATION NOTES

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

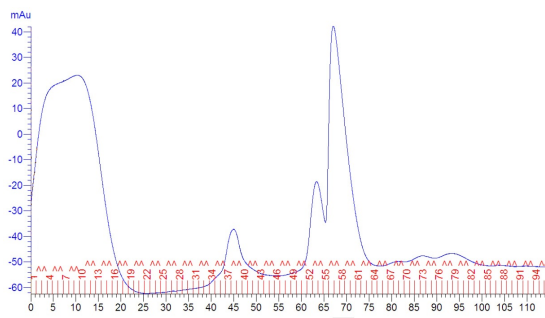
### GENE & PROTEIN INFORMATION

<b>UniProt ID</b>	H3 - P68431 H4 - P62805
-------------------	----------------------------

VALIDATION DATA



**FIGURE 1 Protein gel data.** Histone H3/H4 Tetramer, Recombinant Human (1  $\mu$ g) 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.



**FIGURE 2 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.