

## Histone CENP-A/H4 Tetramer, Recombinant Human

|                      |             |                |                |
|----------------------|-------------|----------------|----------------|
| <b>Catalog No</b>    | 16-0010     | <b>Species</b> | Human          |
| <b>Lot No</b>        | 23060007-02 | <b>Source</b>  | <i>E. coli</i> |
| <b>Pack Size</b>     | 50 µg       | <b>Tag</b>     | None           |
| <b>Concentration</b> | 19 µM       | <b>MW</b>      | 54 kDa         |

### DESCRIPTION

Human CENP-A/Histone H4 Tetramers made from recombinant histones expressed in *E. coli*. 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 a component of modified nucleosomes found at the core centromere.

### TECHNICAL INFORMATION

|                    |   |
|--------------------|---|
| <b>Storage</b>     | DO NOT DEEP FREEZE (-80°C)!! Stable for six months at -20°C from date of receipt. For best results, aliquot and avoid freeze/thaws. |
| <b>Formulation</b> | 1.0 mg/mL CENP-A/H4 Tetramer in 48.7 µL 10 mM sodium phosphate buffer pH 6.9, 1 M NaCl, 7 mM β-ME, 0.5 mM PMSF, and 50% glycerol    |

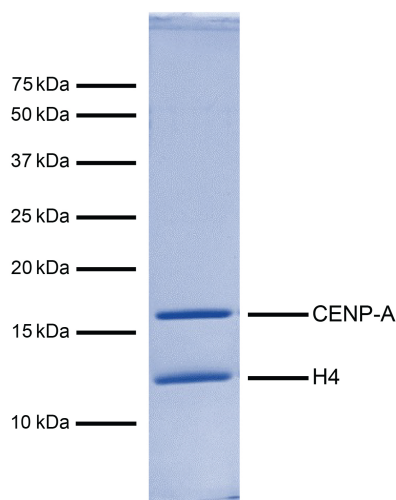
### APPLICATION NOTES

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

### GENE & PROTEIN INFORMATION

|                   |   |
|-------------------|---|
| <b>UniProt ID</b> | CENP-A - P49450 (alt. names: Centromere protein A, CENPA, CenH3)<br>H4 - P62805 |
|-------------------|---|

### VALIDATION DATA



**FIGURE 1 Protein gel data.** Coomassie stained SDS-PAGE gel of proteins in CENP-A/H4 Tetramer (1 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the histones (CENP-A and H4) are indicated.