

Histone CENP-A/H4 Tetramer, Recombinant Human

| | | | |
|----------------------|-------------|----------------|----------------|
| Catalog No | 16-0010 | Species | Human |
| Lot No | 25338008-01 | Source | <i>E. coli</i> |
| Pack Size | 50 µg | Tag | None |
| Concentration | 20.4 µM | MW | 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 defining component of centromere-specific nucleosomes at the core centromere.

TECHNICAL INFORMATION

| | |
|--------------------|---|
| Storage | 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. |
| Formulation | 1.1 mg/mL CENP-A/H4 Tetramer in 45.5 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

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

VALIDATION DATA

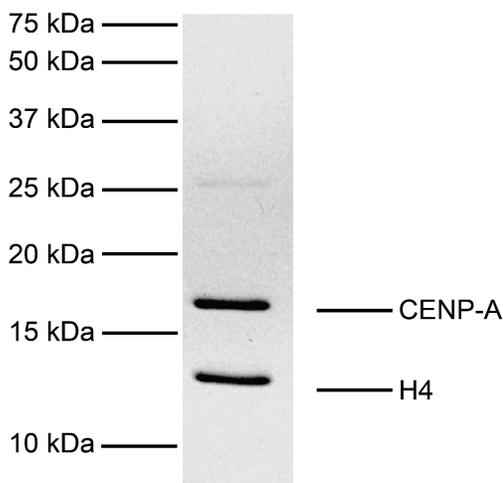


FIGURE 1 Protein gel data. Coomassie stained SDS-PAGE gel of proteins in CENP-A/H4 Tetramer (1 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the histones (CENP-A and H4) are indicated.