

Histone CENP-A/H4 Tetramer, Recombinant Human

Catalog No 16-0010

Lot No 22039006-01

Pack Size 50 µg



EpiCypher®

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.

Formulation:

50 µg of CENP-A/H4 Tetramer (1 mg/mL) in 10 mM sodium phosphate buffer pH 6.9, 1 M NaCl, 7 mM β-ME, 0.5 mM PMSF, and 50% glycerol.

Storage and Stability:

DO NOT DEEP FREEZE (-80°C)!! Stable for six months at -20°C from date of receipt. For best results, aliquot for single use.

Application Notes:

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

PAGE gel (Figure 1) serves to confirm that both histone subunits are present in equimolar amounts and does not represent the size of the intact tetramer.

References:

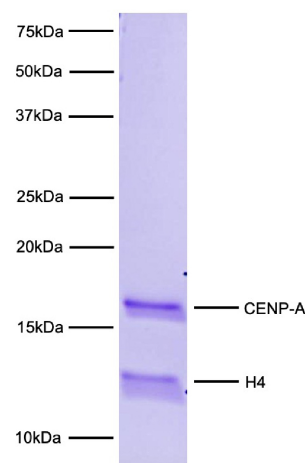


Figure 1: Protein gel data. Coomassie stained PAGE gel of proteins in CENP-A/H4 Tetramer (1 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and positions of the histones (CENP-A and H4) are indicated.

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