

Histone Octamer (H4 Δ N15), Human Recombinant



EpiCypher™

Catalog No. 16-8018
Lot No. 17255001
Pack Size 50 μ g

Product Description:

Histone octamer assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4 Δ N15. Accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805) with the amino acid sequence of H4 beginning with lysine 16 (amino acids 1-15 are deleted).

Formulation:

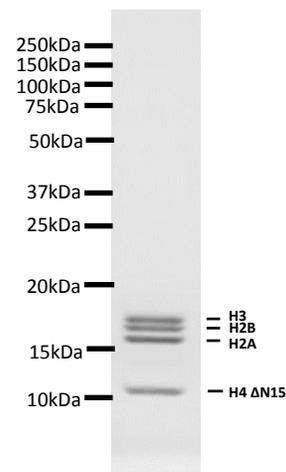
Purified recombinant histone octamers (50 μ g) in 104 μ l 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2M NaCl, 2 mM DTT, & 20% glycerol. Concentration of histone octamer is 4.53 μ M. Histone octamer molecular weight = 106,037.4Da.

Storage and Stability:

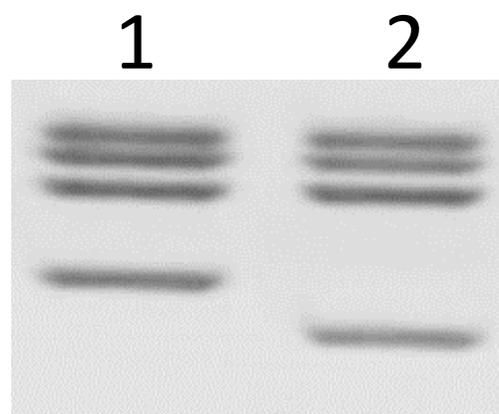
Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Histone Octamer (H4 Δ N15), Human Recombinant are highly purified and suitable for use as substrates in enzyme screening assays, structural studies, or effector protein binding experiments. The N-terminal deletion allows for the study of the role of the N-terminus in many aspects of chromatin biology.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Histone Octamer (H4 Δ N15), Human Recombinant (1 μ g) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4 Δ N15) are indicated.



Detailed Protein Gel Data: Coomassie stained PAGE gel of proteins in wildtype histone octamers (Lane 1) resolved alongside proteins in Histone Octamer (H4 Δ N15), Human Recombinant (Lane 2). The faster migration of the H4 species observed in Lane 2 as compared to Lane 1 denotes the deletion of the amino acids from the N-terminus of the histone protein.

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