

Histone Octamer (H3.1ΔN32), Human Recombinant

Catalog No. 16-8016
Lot No. 17255001
Pack Size 50 µg



EpiCypher®

Product Description:

Histone Octamer assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3.1ΔN32 and H4. Accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805), with the amino acid sequence of H3 beginning with glycine 33 (amino acids 1-32 are deleted).

Formulation:

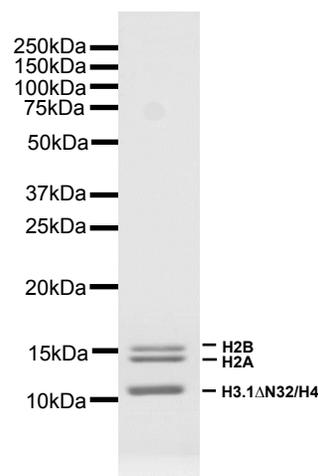
Purified recombinant histone octamer (50µg) in 20.8µl 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 2M NaCl, 2 mM DTT, & 20% glycerol. Concentration of histone octamer is 23.54 µM. Histone octamer molecular weight = 101,934.6 Da.

Storage and Stability:

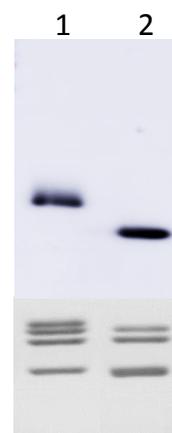
Stable for six (6) months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

Histone Octamer (H3.1ΔN32), 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 (H3.1ΔN32), Human Recombinant (1 µg) demonstrates the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3.1ΔN32 and H4) are indicated. H3.1ΔN32 and H4 co-migrate.



Western Blot Data: Western Analysis of Histone octamer (H3.1ΔN32), Human Recombinant. Top Panel: Wild type (Lane 1) and H3.1ΔN32-containing octamers (Lane 2) were probed with an anti-H3 COOH-terminal antibody and analyzed via ECL readout. Bottom Panel: Detail from Coomassie stained gel showing histones from H3.1 WT (Lane 1) and H3.1ΔN32 histone octamer (Lane 2).

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