

Nucleosome, Recombinant Human, H4 Tetraacetyl

Catalog No. 16-0313
Lot No. 17031002
Pack Size 50 µg



Product Description:

Mononucleosomes assembled from recombinant human histones expressed in *E. coli* (two each of histones H2A, H2B, H3 and H4; accession numbers: H2A-P04908; H2B-O60814; H3.1-P68431; H4-P62805) wrapped by 147 base pairs of 601 positioning sequence DNA. Histone H4 (created by a proprietary semi-synthetic method) contains N-term. α-acetylation and acetyl-lysine at positions 5, 8, 12 and 16. The nucleosome is the basic subunit of chromatin. The 601 sequence, identified by Lowary and Widom, is a 147-base pair sequence that is useful for nucleosome assembly and contains a 5' biotin TEG group.

Formulation:

Recombinant H4 Tetraacetyl mononucleosomes (27.3 µg protein by mass, 50.0 µg protein+DNA in 42.0 µl) in 10 mM Tris-HCl pH 7.5, 1 mM EDTA, 25 mM NaCl, 2 mM DTT, & 20% glycerol. Concentration of nucleosomes is 5.92 µM. Nucleosome molecular weight = 201,031 Da.

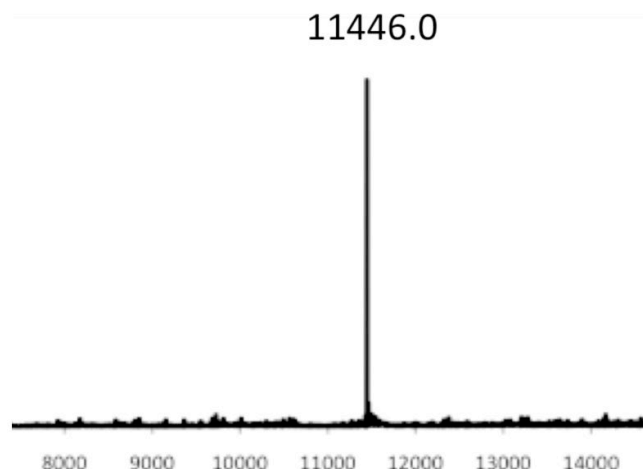
Storage and Stability:

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

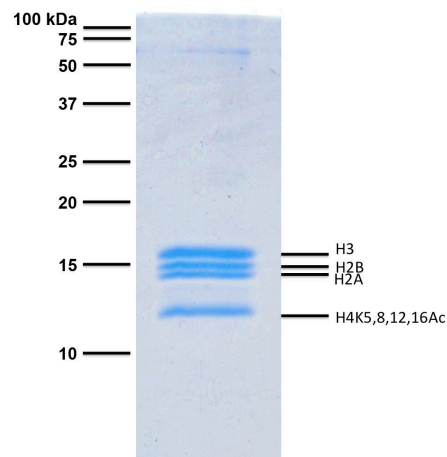
Application Notes:

Nucleosome, Recombinant Human, H4 Tetraacetyl are highly purified and are suitable for use as substrates in enzyme screening assays or for effector protein (especially bromodomain) binding experiments. **Nucleosome, Recombinant Human, H4 Tetraacetyl from EpiCypher does not contain free DNA which could alter assayed activities.**

References Using this Product:

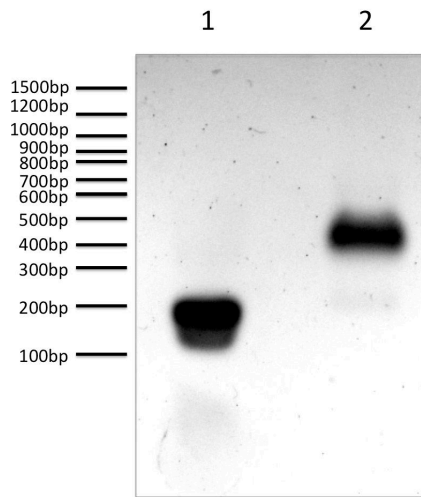


Mass Spec Data: Synthetic H4 Tetraacetyl protein analyzed by ESI-TOF mass spectrometry. Expected mass = 11446.0 Da. Determined mass = 11446.0 Da.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Nucleosome, Recombinant Human, H4 Tetraacetyl (1 µg) to demonstrate the purity of the histones in the preparation. Sizes of molecular weight markers and positions of the core histones (H2A, H2B, H3 and H4 Tetraacetyl) are indicated.

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



DNA Gel Data: Nucleosome, Recombinant Human, H4 Tetraacetyl
run on an agarose gel and stained with ethidium bromide to
visualize DNA.

Lane 1: Free DNA extracted from nucleosomes (200 ng).

Lane 2: Intact nucleosomes (400 ng).