

Chicken Polynucleosomes, Purified

Catalog No 16-0004
Lot No 19161002
Pack Size 50 µg



EpiCypher®

Product Description:

Polynucleosomes purified from chicken erythrocytes. The nucleosome is the basic subunit of chromatin consisting of the histone octamer (two each of the four core histones, H2A, H2B, H3 and H4) wrapped by 147 base pairs of DNA. As shown by the size of the DNA species, Chicken Polynucleosomes are predominantly dimers, trimers and tetramers.

Formulation:

Purified Chicken Polynucleosomes at a concentration of 1.8 mg/mL (DNA + protein) in 10 mM Tris-HCl pH 8.0, 1 mM EDTA and 10% glycerol.

Application Notes:

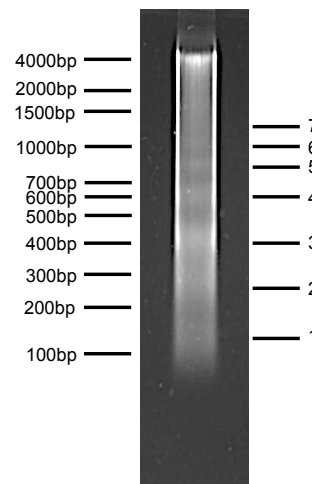
Chicken Polynucleosomes are suitable for use in enzyme assays such as acetylation, methylation or phosphorylation, chromatin binding assays, or for use as a positive control in Western blotting.

Storage and Stability:

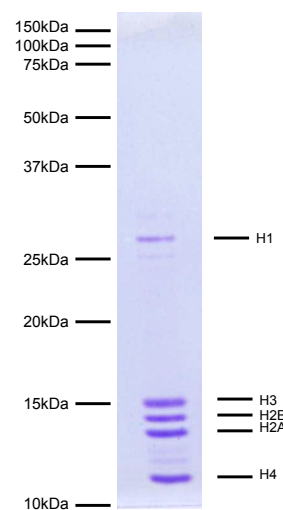
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

References:

Adhvaryu KK et al (2005) Eukaryot Cell 4:1455-1564.
Kizer KO et al (2005) Mol Cell Biol 25:3305-3316.
Morris SA et al (2005) Eukaryot Cell 4:1446-1454.

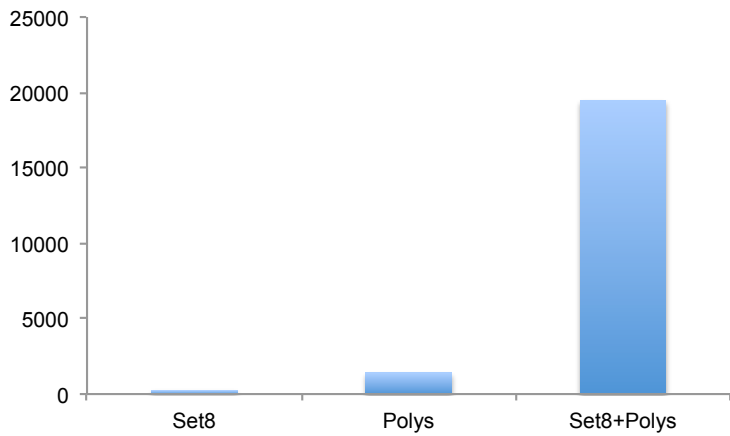


DNA Gel Data: DNA was purified from Chicken Polynucleosomes (4 µg) and run on an agarose gel to show the size of nucleosomal DNA compared to molecular weight markers (base pairs). Note that while mononucleosomal DNA runs at 147 base pairs, oligonucleosomal DNA is roughly 200 base pairs per nucleosome.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Chicken Polynucleosomes (1.5 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers (kDa) and positions of the core histones (H2A, H2B, H3 and H4) and histone H1 are indicated.

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



Histone Methyltransferase Data: Chicken Polynucleosomes were used in a radioactive histone methyltransferase assay with recombinant SET8. Radioactive SAM was incubated with Chicken Polynucleosomes (Polys, 1.5 μ g), recombinant SET8 (SET8, 1 μ g), or both (SET8+Polys).

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