

Chicken Mononucleosomes, Purified

Catalog No. 16-0019
Lot No. 17248001
Pack Size 50 µg

Product Description:

Mononucleosomes 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.

Formulation:

Purified Chicken Mononucleosomes at a concentration of 0.52 mg/ml (DNA + protein) in 10 mM Tris-HCl pH 8.0, 1 mM EDTA and 10% glycerol. Approx. MW = 225,000 Da
Approx. concentration = 2.3 µM.

Concentration may differ from previous lots.

Storage and Stability:

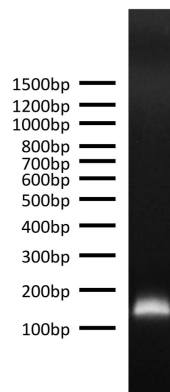
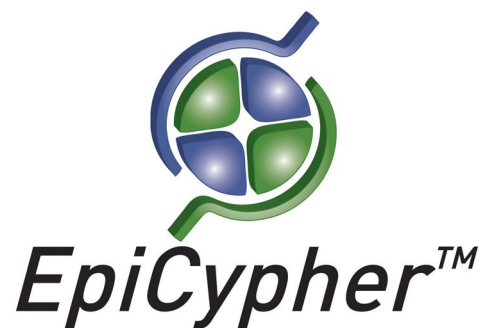
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

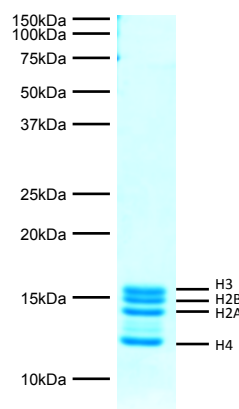
Chicken Mononucleosomes are suitable for use in enzyme assays such as acetylation, methylation or phosphorylation, chromatin binding assays, or for use as a positive control in Western blotting.

References:

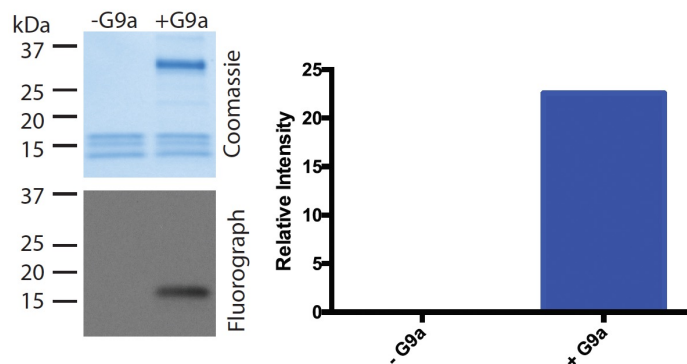
Adhvaryu KK *et al* (2005). *Eukaryot Cell* 4: 1455-1564.
Kizer KO *et al* (2005). *Mol Cell Biol* 25: 3305-3316.
Morris SA *et al* (2005). *Eukaryot Cell* 4: 1446-1454.



DNA Gel Data: DNA was purified from Chicken Mononucleosome (1 µg) and run on an agarose gel to show the size of nucleosomal DNA compared to molecular weight markers (base pairs). Note that the mononucleosomal DNA runs at ~150 base pairs.



Protein Gel Data: Coomassie stained PAGE gel of proteins in Chicken Mononucleosomes (2 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) are indicated.



Histone Methyltransferase Data: Chicken Mononucleosomes were used in a radioactive histone methyltransferase assay with recombinant G9a. Radioactive SAM was incubated with Chicken Mononucleosomes in the absence (-) or presence (+) of recombinant G9a and run on an acrylamide gel. HMTase activity co-purifying with the nucleosomes is undetectable. **Left Top:** Coomassie stained gel. **Left Bottom:** autoradiogram of gel. **Right:** Densitometry quantification of autoradiogram.

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