

Chicken Mononucleosomes, Purified

Catalog No	16-0019	Species	Chicken
Lot No	23212003-08	Source	Chicken erythrocytes
Pack Size	50 µg	Tag	None
Concentration	4.8 µM	MW	225,000 Da

DESCRIPTION

Mononucleosomes purified from chicken erythrocytes. The nucleosome is the basic repeating unit of chromatin, which consists of 147 base pairs of DNA wrapped around an octamer core of histone proteins (two each of H2A, H2B, H3, and H4).

TECHNICAL INFORMATION

Storage	Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid freeze/thaws.
Formulation	1.09 mg/mL purified Chicken Mononucleosome in 45.9 µL 10 mM Tris-HCl pH 8.0, 1 mM EDTA, and 10% glycerol.

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.

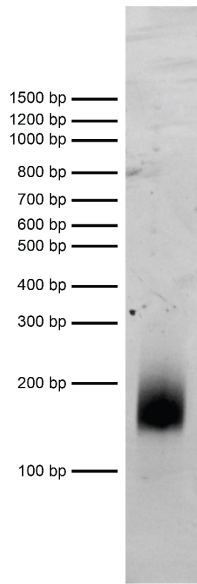


FIGURE 1 DNA gel data. Purified chicken mononucleosome DNA (100 ng) resolved via native PAGE gel and stained with ethidium bromide.

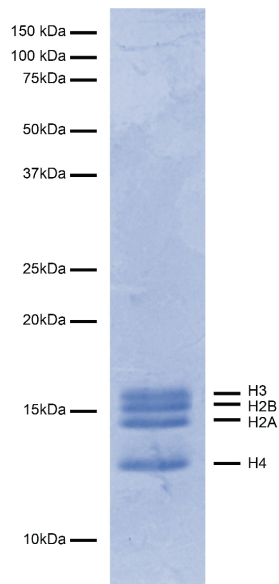


FIGURE 2 Protein gel data. Coomassie stained SDS-PAGE gel of proteins in Chicken Mononucleosomes (2 µg) demonstrates the purity of histones in the preparation. Sizes of molecular weight markers and position of the core histones (H2A, H2B, H3 and H4) are indicated.

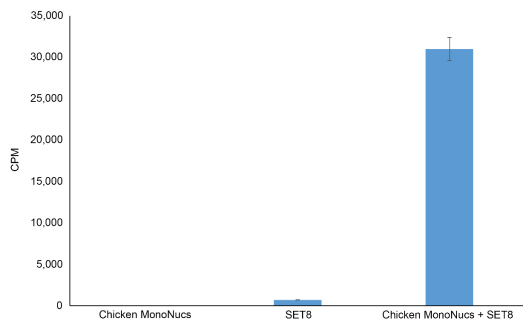


FIGURE 3 Enzyme activity data. HMTase assay with Chicken Mononucleosomes (1 µg) and SET8 (Chicken MonoNucs + SET8; 1 µg) using a standard radiometric filter binding assay protocol. Controls include the mononucleosomes alone (Chicken MonoNucs) or SET8 alone (SET8). SET8 is aa 1-352 of UniProt# Q9NQR1.