

HeLa Polynucleosomes, Purified

Catalog No. 16-0003
Lot No. 13059001
Pack Size 50 µg

Product Description:

Human polynucleosomes purified from HeLa cells. 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. HeLa Polynucleosomes are predominantly trimers, with some dimers and tetramers.

Formulation:

Purified HeLa polynucleosomes at a concentration of 6 mg/ml (DNA + protein) in 20 mM HEPES, pH 7.5, 1 mM EDTA.

Storage and Stability:

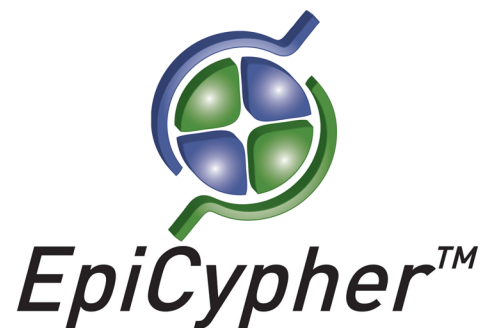
Stable for six months at -80°C from date of receipt. For best results, aliquot and avoid multiple freeze/thaws.

Application Notes:

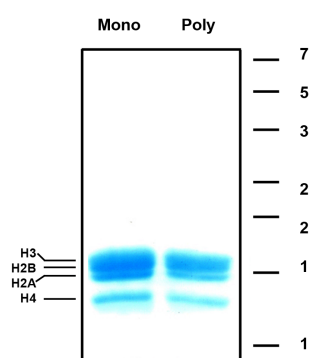
HeLa polynucleosomes are supplied concentrated to maintain their stability. Use the included **Dilution Buffer 1** to dilute them to your desired final concentration. HeLa poly-nucleosomes are suitable for use in enzyme assays such as acetylation or methylation, chromatin binding assays, or for use as a positive control in Western blotting. Use 1-2 µg per reaction.

References:

Kuo AJ et al (2011). *Mol Cell* 44: 609-620.
Matthews AG et al (2007). *Nature* 450: 1106-1110.
Shi X et al (2006). *Nature* 442: 96-99.



DNA Gel Data: DNA was purified from HeLa Mono-nucleosomes (Mono, Catalog No. 16-0002) and HeLa Polynucleosomes (Poly) and run on an agarose gel to show the size of nucleosomal DNA compared to molecular weight markers (base pairs). Note that mononucleosomal DNA runs at ~147 base pairs.



Protein Gel Data: Coomassie stained PAGE gel of proteins in HeLa Mononucleosomes (Mono, Catalog No. 16-0002) and HeLa Polynucleosomes (Poly) to demonstrate the purity of the histones in the preparations. Sizes of molecular weight markers (kDa) and position of the core histones are indicated.