

## Human Native Nucleosome

**Cat. No.** C01030102

**Source:** E. coli

**Lot #:** 001

**Size:** 20 µg/ 100 µl

**Concentration:** 0.2 µg/µl

**Specificity:** Human

**Purity:** Purified using glycerol gradient and ion exchange chromatography, >98% purity as determined by SDS-PAGE.

**Storage buffer:** PBS containing 0.5 mM PMSF and 50% glycerol.

**Storage:** Store at -80°C; guaranteed stable for 1 year from date of receipt when stored properly.

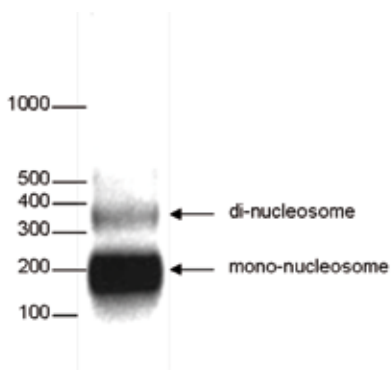
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**Description:** Human native nucleosomes purified from HeLa cells, digested with micrococcal nuclease.

### Protein description

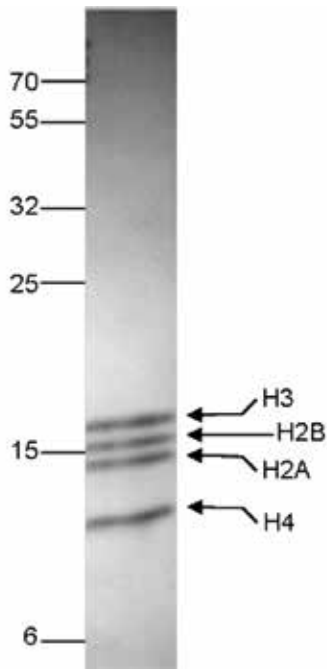
The nucleosome is the smallest part of the chromosome. It consists of two core histones of each class H2A, H2B, H3 and H4 which assemble and are wrapped around 146 base pairs of DNA. Histones are the main constituents of the protein part of chromosomes of eukaryotic cells. They pack the DNA into tight masses of chromatin. Histones are rich in the amino acids arginine and lysine and have been greatly conserved during evolution. Histone tails undergo numerous post-translational modifications, which either directly or indirectly alter chromatin structure to facilitate transcriptional activation or repression or other nuclear processes.

### Quality control



**Figure 1.**

Agarose gel analysis of the human native nucleosome. The position of the mono- and di-nucleosomes is indicated on the right; the marker (in bp) is shown on the left.



**Figure 2.**

SDS-PAGE gel analysis of the human native nucleosome. The position of the four core histones is indicated on the right; the marker (in kDa) is shown on the left.

**Diagenode sa. BELGIUM | EUROPE**

LIEGE SCIENCE PARK  
Rue Bois Saint-Jean, 3  
4102 Seraing (Ougrée) - Belgium  
Tel: +32 4 364 20 50  
Fax: +32 4 364 20 51  
orders@diagenode.com  
info@diagenode.com

**Diagenode Inc. USA | NORTH AMERICA**

400 Morris Avenue, Suite 101  
Denville, NJ 07834 - USA  
Tel: +1 862 209-4680  
Fax: +1 862 209-4681  
orders.na@diagenode.com  
info.na@diagenode.com

**Last update:** September 10, 2015