

Human Recombinant Nucleosome

Cat. No. C01030101

Source: E. coli

Lot #: 001

Size: 5 µg/ 20 µl

Concentration: 0.25 µg/µl

Specificity: Human

Purity: Purified using FPLC, >98% purity as determined by SDS-PAGE

Storage buffer: 20 mM Tris-Cl pH 7.9, 1mM EDTA, 0.5 mM PMSF, 1 mM DTT and 40% glycerol.

Storage: Store at -20°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: Recombinant nucleosome, assembled in vitro by salt dialysis using the 147 bp 601 DNA sequence and the four core histones, H2A, H2B, H3 and H4, produced in E. Coli. Suitable for in vitro chromatin assembly and enzyme activity assays

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

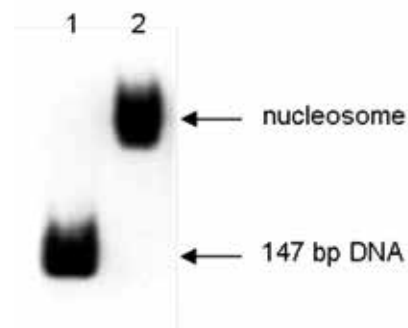


Figure 1.

Native page of the human recombinant nucleosome (lane 2). Lane 1 shows the 147 bp 601 DNA sequence.

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