

Recombinant Histone H3

Cat. No. C23010003

Source: E. coli

Lot #: 001

Size: 100 µg/ 200 µl

Concentration: 0.5 µg/µl

Specificity: Human

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

Storage buffer: 20 mM sodium phosphate pH 7.0, 0.3 M NaCl, 1mM EDTA, 0.5 mM PMSF and 1 mM DTT.

Storage: Store at -80°C; guaranteed stable for 2 years from date of receipt when stored properly.

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

Description: Full length recombinant histone H3, produced in E. coli.

Protein description

Histones are the main constituents of the protein part of chromosomes of eukaryotic cells. They are rich in the amino acids arginine and lysine and have been greatly conserved during evolution. Histones pack the DNA into tight masses of chromatin. Two core histones of each class H2A, H2B, H3 and H4 assemble and are wrapped by 146 base pairs of DNA to form one octameric nucleosome. 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

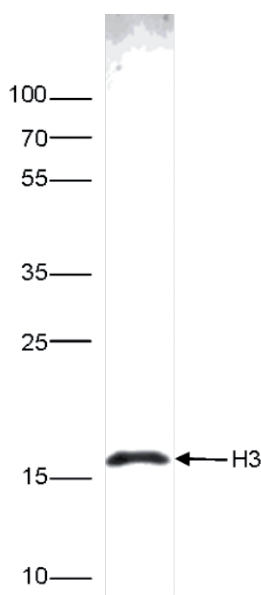


Figure 1.

SDS page of the Recombinant Histone H3. The position of the protein of interest 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 1, 2015