

HP1 α recombinant protein

Other names: CBX1, MOD1, M31, P25 β

Cat. No. C23010101

Source: E. coli

Lot #: 001

Size: 25 μ g/ 50 μ l

Concentration: 0.5 μ g/ μ l

Specificity: Human

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

Storage buffer: 20 mM Tris-Cl pH 7.6, 1 mM EDTA, 0.15 M NaCl, 10% glycerol, 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 human N-terminal His tagged HP1 α (Heterochromatin Protein 1 Homolog Alpha) protein, produced in E. coli.

Protein description

HP1 alpha (UniProt/Swiss-Prot entry P45973) is a component of heterochromatin. It recognizes and binds to histone H3 tails methylated at 'Lys-9', leading to epigenetic repression of transcription. HP1 alpha may also interact with lamin B receptor (LBR), thereby contributing to the association of the heterochromatin with the inner nuclear membrane. Further, HP1 alpha is involved in the formation of a functional kinetochore through interaction with MIS12 complex proteins.

Quality control

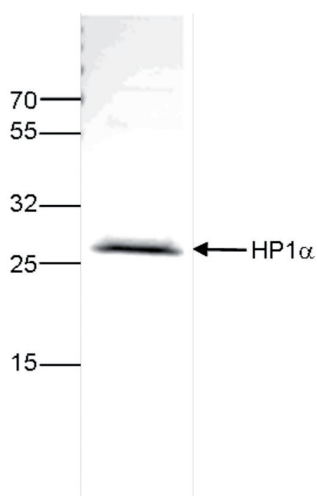


Figure 1.

SDS page of the HP1 α recombinant protein. 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: August 19, 2015