

ASF1b recombinant protein

Other names: CIA-II, HAsf1b, HCIA-II

Cat. No. C23010105

Source: E. coli

Lot #: 001

Size: 25 µg/ 50 µl

Concentration: 0.5 µg/µl

Specificity: Human

Purity: Purified using FPLC, >95% purity as determined by SDS-PAGE.

Storage buffer: 20 mM Tris-Cl pH 7.6, 1 mM DTT, 0.15 M NaCl, 10% glycerol.

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 GST tagged ASF1b (Anti-Silencing Function 1B Histone Chaperone) protein, produced in E. coli.

Protein description

ASF1b (UniProt/Swiss-Prot entry Q9NVP2) belongs to the family of histone chaperones and interacts with both histones H3 and H4. It's a key component for nucleosome assembly and disassembly. ASF1b cooperates with chromatin assembly factor 1 (CAF-1) to promote replication-dependent chromatin assembly, but is not involved in replication-independent chromatin assembly. ASF1b is essential for spermatogenesis.

Quality control

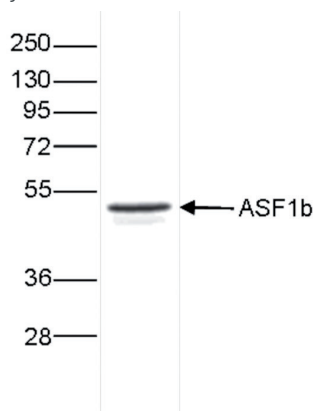


Figure 1.

SDS page of the ASF1b 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 25, 2015