

## 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.

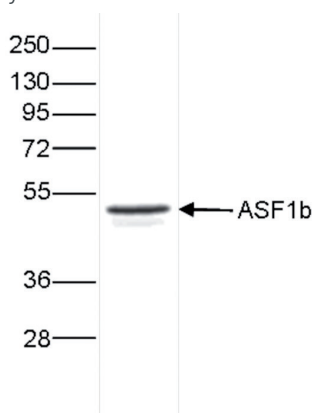
**Precautions:** This product is for research use only. Not for use in diagnostic or therapeutic procedures.

**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