

ASF1a recombinant protein

Other names: CIA, HAsf1, HCIA, HSPC146, CGI-98

Cat. No. C23010104

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 ASF1a (Anti-Silencing Function 1A Histone Chaperone) protein, produced in E. coli.

Protein description

ASF1a (UniProt/Swiss-Prot entry Q9Y294) 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 and promotes both replication-dependent chromatin assembly and with HIRA to promote replication-independent chromatin assembly.

Quality control

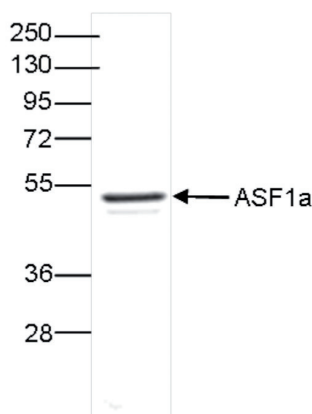


Figure 1.

SDS page of the ASF1a 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 21, 2015