

UBA1 recombinant protein

Other names: UBA1A, UBE1, UBE1X, A1S9, A1S9T, AMCX1, SMAX2, A1ST, GXP1, CFAP124, POC20

Cat. No. C23010111

Source: Baculovirus

Lot #: 001

Size: 5 µg/ 10 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, 0.15 M NaCl, 10% glycerol, 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 UBA1 (Ubiquitin-Like Modifier Activating Enzyme 1) protein, produced in Baculovirus.

Protein description

UBA1 (UniProt/Swiss-Prot entry P22314) catalyzes the first step in ubiquitin conjugation to mark cellular proteins for degradation. It activates ubiquitin by adenylating the C-terminal glycine residue with ATP and subsequently linking the adenylated glycine to the side chain of a cysteine residue in E1 thereby forming a ubiquitin-E1 thioester. UBA1 is also required for DNA repair response to replication stress.

Quality control

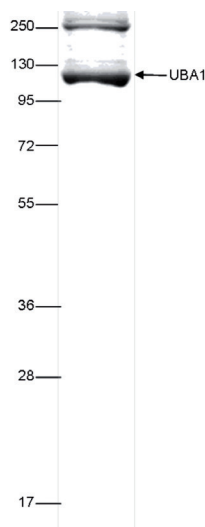


Figure 1.

SDS page of the UBA1 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 28, 2015