

UBLE1A (SAE1) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # U208-30G

Lot # F460-3

Product Description

Recombinant full-length human UBLE1A (SAE1) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM_005500](#).

Gene Aliases

SAE1; SUA1; HSPC140; AOS1; FLJ3091; HSPC140

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

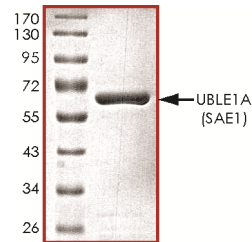
Scientific Background

UBLE1A (also known as SAE1) or SUMO1 activating enzyme subunit 1 is involved in regulating protein structure and intracellular localization. SAE1 and UBA2 form a heterodimer that functions as a SUMO-activating enzyme for the sumoylation of proteins (1). The SAE1/SAE2 dimer functions in SUMO1 activation in a manner analogous to the single E1 ubiquitin-activating enzymes. The SAE2 inactivation may be a therapeutic strategy in MYC-driven cancers (2).

References

- Okuma, T. et.al: In vitro SUMO-1 modification requires two enzymatic steps, E1 and E2. *Biochem. Biophys. Res. Commun.* 254: 693-698, 1999.
- Kessler, J. D. et.al: A SUMOylation-dependent transcriptional subprogram is required for Myc-driven tumorigenesis. *Science* 335: 348-353, 2012.

Purity



The purity of UBLE1A (SAE1) was determined to be **>95%** by densitometry. Approx. MW **65kDa**.

UBLE1A (SAE1) Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	U208-30G
Specific Lot Number	F460-3
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr At -70°C from date of shipment
Storage & Shipping	Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.