

## Anti-UBLE1A (SAE1)

Rabbit Polyclonal Antibody

**Catalog # U208-63R**

Lot # O2121-29

### Cited Applications

WB, ELISA

*Ideal working dilutions for each application should be empirically determined by the investigator.*

### Specificity

Recognizes the UBLE1A (SAE1) protein

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

UBLE1A antibody was raised against a 20 amino acid synthetic peptide near the carboxy terminus of human UBLE1A

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

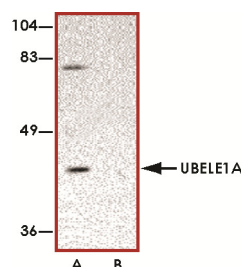
### Scientific Background

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.

### Sample Data



Western blot analysis of UBLE1A in SK-N-SH lysate with UBLE1A antibody at 1 µg/ml in (A) the absence and (B) the presence of blocking peptide.

## Anti-UBLE1A (SAE1)

Rabbit Polyclonal Antibody

Catalog Number

U208-63R

Specific Lot Number

O2121-29

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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

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