

Anti-UBA1 (UBE1)

Rabbit Polyclonal Antibody

Catalog # U201-62R

Lot # B3216-90

Cited Applications

ELISA, IHC, WB

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

Specificity

Recognizes the human Ubiquitin Activating Enzyme E1 protein

Cross Reactivity

Human

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

The antibody was produced against recombinant protein corresponding to full-length human Ubiquitin Activating Enzyme E1

Formulation

0.02 M Potassium Phosphate, 0.15 M Sodium Chloride, pH 7.2 + 0.01% (w/v) Sodium Azide

Stability

1yr at -20°C from date of shipment

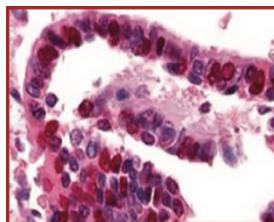
Scientific Background

UBA1 or ubiquitin-like modifier activating enzyme 1 is an E1 enzyme that initiates the activation and conjugation of ubiquitin-like proteins and catalyzes the first step in ubiquitin conjugation to mark cellular proteins for degradation (1). UBE1 is able to transfer ubiquitin to a wide range of E2 substrates. UBA1 protein may participate in DNA topoisomerase-2 activity (2). UBA1 complements an X-linked mouse temperature-sensitive defect in DNA synthesis, and thus may function in DNA repair.

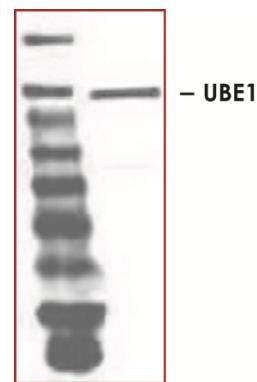
References:

1. Jin, J.et.al: Dual E1 activation systems for ubiquitin differentially regulate E2 enzyme charging. *Nature* 447: 1135-1138, 2007.
2. Zackenhaus, E.et.al: Identification of human gene complementing ts A1S9 mouse L-cell defect in DNA replication following DNA-mediated gene transfer. *Somat. Cell Molec. Genet.* 14: 371-379, 1988.

Sample Data



Immunohistochemical analysis of formalin-fixed, paraffin embedded human lung tissue using Anti-UBA1 (UBE1) (10 µg/ml)human lung tissue.



Western blot using Anti-UBA1 (UBE1) (1:1,000 dilution) shows detection of a band at ~118 kDa corresponding to UBE1 (lane 1 800 nm channel). Load: ~35 µg of an A431 whole cell lysate.

Anti-UBA1 (UBE1)

Rabbit Polyclonal Antibody

Catalog #	U201-62R
Lot #	B3216-90
Purification	Multistep procedure
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
www.signalchem.com

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