

Anti-UBC9

Rabbit Polyclonal Antibody

Catalog # U19-63R

Lot # R027-6

Cited Applications

- Western blot (1:1000 dilution)

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

Specificity

Recognizes the UBC9 protein

Cross Reactivity

- Western blot UBC9 from human and mouse cells

UBC9 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide of c-terminal amino acid sequence of human UBC9 (P63279)

Formulation

TBS, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

Scientific Background

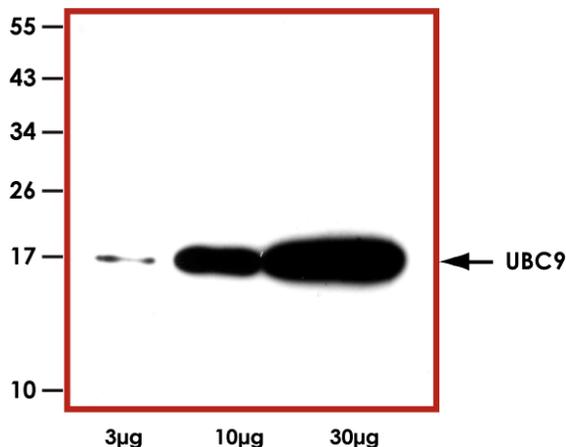
UBC9 is a member of the E2 ubiquitin-conjugating enzyme family that is involved in ubiquitin-dependent protein degradation system (1). Ubiquitination involves at least three classes of enzymes: ubiquitin-activating enzymes, ubiquitin-conjugating enzymes and ubiquitin-protein ligases. UBC9 can function as a co-factor of PLAGL2, uncoupling its enzymatic activity to mediate PLAGL2 interactive SP-C promoter activity (2). UBC9 also plays an important role in the up-regulation of the bone morphogenetic protein signaling pathway.

References

1. Yasugi, T. et.al: Identification of the structural and functional human homolog of the yeast ubiquitin conjugating enzyme UBC9. *Nucleic Acids Res.* 24: 2005-2010, 1996.
2. Wang, Z.-Y. et.al: Molecular cloning of the cDNA and chromosome localization of the gene for human ubiquitin-conjugating enzyme 9. *J. Biol. Chem.* 271: 24811-24816, 1996.

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

Sample Data



Representative western blot with Anti-UBC9 (1:1000) using 3 µg, 10 µg and 30 µg of Jurkat cell lysate.

Anti-UBC9

Rabbit Polyclonal Antibody

Catalog Number

U19-63R

Specific Lot Number

R027-6

Purification

Affinity chromatography

Concentration

1.0 µg/µl

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.