

UBE2N antibody (pAb)

Catalog No: 39224

RRID: AB_2793194

Application(s): WB

Reactivity: Human

Quantity: 100 µg

Purification: Affinity Purified

Host: Rabbit

Isotype: IgG

Concentration: 0.5 µg/µl

Molecular Weight: 15 kDa

Background: UBE2N (UBC13, BLU, ubiquitin-conjugating enzyme E2N) heterodimer with UBE2V2 catalyzes the synthesis of non-canonical poly-ubiquitin chains that are linked through 'Lys-63'. This type of poly-ubiquitination does not lead to protein degradation by the proteasome. UBE2N mediates transcriptional activation of target genes and plays a role in the control of progress through the cell cycle and differentiation. UBE2N also plays a role in the error-free DNA repair pathway and contributes to the survival of cells after DNA damage and acts together with the E3 ligase SHPRH in the poly-ubiquitination of PCNA 'Lys-164' upon genotoxic stress, which is required for DNA repair.

Immunogen: This UBE2N antibody was raised against a mixture of synthetic peptides corresponding to amino acid residues 2-19 and 131-148 of human UBE2N.

Buffer: PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

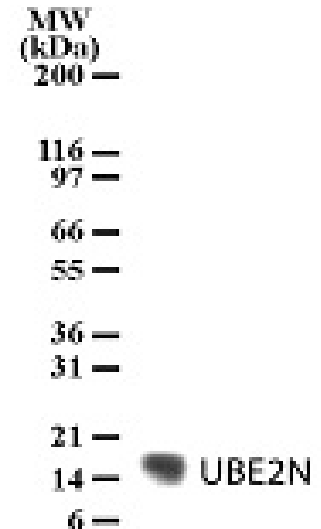
Applications Validated by Active Motif:

WB: 2 µg/ml dilution

For optimal results in Western blotting, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



UBE2N pAb tested by Western blot.
UBE2N detection by Western blot. The analysis was performed using human heart whole-cell extract and UBE2N pAb at a 2 µg/ml dilution.