
UBE2V1 Polyclonal Antibody

(Catalog # A63744)

Background

Has no ubiquitin ligase activity on its own. The UBE2V1-UBE2N heterodimer catalyzes the synthesis of non-canonical poly-ubiquitin chains that are linked through Lys-63. This type of poly-ubiquitination activates IKK and does not seem to involve protein degradation by the proteasome. Plays a role in the activation of NF-kappa-B mediated by IL1B, TNF, TRAF6 and TRAF2. Mediates transcriptional activation of target genes. Plays a role in the control of progress through the cell cycle and differentiation. Plays a role in the error-free DNA repair pathway and contributes to the survival of cells after DNA damage. Promotes TRIM5 capsid-specific restriction activity and the UBE2V1-UBE2N heterodimer acts in concert with TRIM5 to generate 'Lys-63'-linked polyubiquitin chains which activate the MAP3K7/TAK1 complex which in turn results in the induction and expression of NF-kappa-B and MAPK-responsive inflammatory genes.

Description

UBE2V1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q13404

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Ubiquitin-conjugating enzyme E2 variant 1 protein (2-147AA)

Storage

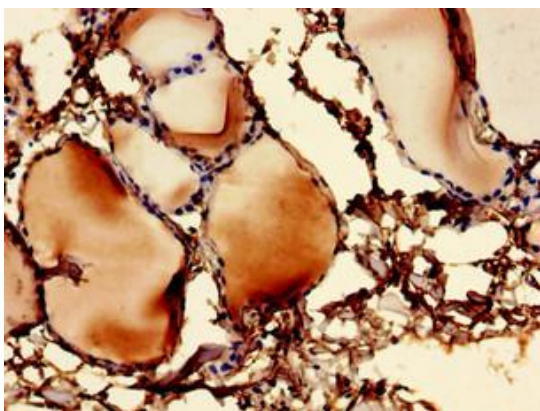
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Ubiquitin-conjugating enzyme E2 variant 1, UEV-1, CROC-1, TRAF6-regulated IKK activator 1 beta Uev1A, UBE2V1, CROC1, UBE2V, UEV1, P/OKcl.19

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human thyroid tissue at dilution of 1:100