

## UBE2V2 Protein

Recombinant protein expressed in E.coli cells

**Catalog # U239-30H**

Lot # J689-3

### Product Description

Recombinant human UBE2V2 (2-end) was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM\\_003350](#).

### Gene Aliases

DDVit-1; DDVIT1; EDAF-1; EDPF1; MMS2; UEV-2; UEV2

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

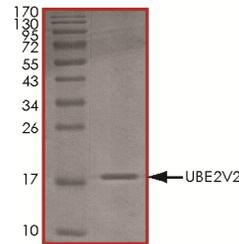
### Scientific Background

UBE2V2 or ubiquitin-conjugating enzyme E2 variant 2 constitutes a distinct subfamily within the E2 protein family which is critical for the catalytic activity of E2s and also shares homology with ubiquitin-conjugating enzyme E2 variant 1 and yeast MMS2 gene product. UBE2V2 may be involved in the differentiation of monocytes and enterocytes (1). UBE2V2 functions in assembly of novel polyubiquitin chains for DNA repair (2).

### References

1. Fritsche, J.et.al: Molecular cloning of a 1-alpha,25-dihydroxyvitamin D3-inducible transcript (DDVit 1) in human blood monocytes. *Biochem. Biophys. Res. Commun.* 235: 407-412, 1997.
2. Hofmann, R. M.et.al: Noncanonical MMS2-encoded ubiquitin-conjugating enzyme functions in assembly of novel polyubiquitin chains for DNA repair. *Cell* 96: 645-653, 1999.

### Purity



The purity of UBE2V2 was determined to be **>95%** by densitometry. Approx. MW **18 kDa**.

## UBE2V2 Protein

Recombinant protein expressed in E. coli cells

Catalog Number	U239-30H
Specific Lot Number	J689-3
Purity	>95%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

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.**