

## UBE2W (UBC16) Protein

Full length recombinant protein expressed in E.coli cells

**Catalog # U240-30G**

Lot # J703-4

### Product Description

Recombinant full length human UBE2W was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [BC010900](#).

### Gene Aliases

hUBC-16; UBC-16; UBC16, FLJ11011

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . 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.

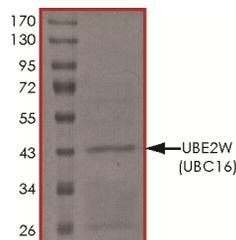
### Scientific Background

UBE2W or ubiquitin-conjugating enzyme E2W (putative) encodes a nuclear-localized ubiquitin-conjugating enzyme (E2) (1). Ubiquitin-conjugating enzyme (E2) along with ubiquitin-activating (E1) and ligating (E3) enzymes coordinates the addition of an ubiquitin moiety to existing proteins. UBE2W interacts with FANCL and regulates the monoubiquitination of Fanconi anemia protein FANCD2 and may be important in the repair of DNA damage (2). UBE2W is highly expressed in heart, brain, liver, and pancreas.

### References

1. Yin, G. et.al: Cloning, characterization and subcellular localization of a gene encoding a human ubiquitin-conjugating enzyme (E2) homologous to the Arabidopsis thaliana UBC-16 gene product. *Front. Biosci.* 11: 1500-1507, 2006.
2. Zhang, Y. et.al: UBE2W interacts with FANCL and regulates the monoubiquitination of Fanconi anemia protein FANCD2. *Molec. Cells* 31: 113-122, 2011.

### Purity



The purity of UBE2W (UBC16) was determined to be **>75%** by densitometry. Approx. MW **45 kDa**.

## UBE2W (UBC16) Protein

Full length recombinant protein expressed in E. coli cells

Catalog Number	U240-30G
Specific Lot Number	J703-4
Purity	>75%
Concentration	0.1 µg/µl
Stability	1yr at $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.**