

## UBE2B Protein

Full length recombinant protein expressed in E. coli cells

**Catalog # U211-30H**

Lot # F736-3

### Product Description

Recombinant full length human UBE2B was expressed in E. coli cells using an N-terminal His tag. The gene accession number is [NM\\_003337](#).

### Gene Aliases

E2-17kDa; HHR6B; HR6B; RAD6B; UBC2

### 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.

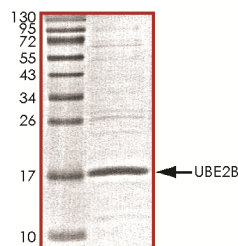
### Scientific Background

UBE2B or ubiquitin-conjugating enzyme E2B is a member of the E2 ubiquitin-conjugating enzyme family which is required for post-replicative DNA damage repair. UBE2B is specifically important during chronic low-dose ultraviolet exposure to prevent counterproductive DNA checkpoint activation and allow cells to proliferate normally (1). The activity of UBE2B is required within the synaptonemal complex and for meiotic recombination in spermatocytes (2). Experimental inactivation of the UBE2B gene in mice causes male infertility.

### References

- Hishida, T. et.al: RAD6-RAD18-RAD5-pathway-dependent tolerance to chronic low-dose ultraviolet light. *Nature* 457: 612-615, 2009.
- Baarends, W. M. et.al: Loss of HR6B ubiquitin-conjugating activity results in damaged synaptonemal complex structure and increased crossing-over frequency during the male meiotic prophase. *Molec. Cell. Biol.* 23: 1151-1162, 2003.

### Purity



The purity of UBE2B was determined to be **>90%** by densitometry. Approx. MW **17 kDa**.

## UBE2B Protein

Full length recombinant protein expressed in E. coli cells

Catalog Number	U211-30H
Specific Lot Number	F736-3
Purity	>90%
Concentration	0.1µg/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.**