# UBE2B (HR6B) [6His-tagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: HHR6B, HR6B, RAD6B, Ubiquitin carrier protein B,

Ubiquitin protein ligase B

Cat. No. 62-0072-100 Quantity: 100 µg -70°C Lot. No. 1823 Storage:

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



**CERTIFICATE OF ANALYSIS Page 1 of 2** 

# **Background**

The enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including regulated and targeted proteasomal degradation of substrate proteins. Three classes of enzymes are involved in the process of ubiquitylation; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). UBE2B is a member of the E2 ubiquitin-conjugating enzyme family and cloning of the human gene was first described by Koken et al. (1991). UBE2B shares 70% identity with its veast homologue but lacks the acidic C-terminal domain. The ring finger proteins RAD5 and RAD18 interact with UBE2B and other members of the RAD6 pathway (Notenboom et al., 2007; Ulrich and Jentsch, 2000). In complex UBE2B and RAD18 trigger replication fork stalling at DNA damage sites during the post replicative repair process (Tsuji et al., 2008). Null mutations of the UBE2B gene in mice are associated with structural abnormalities in sperm and SNP analysis of human UBE2B variants has provided evidence for association of this gene with male infertility (Escalier et al., 2003; Suryavathi et al., 2008).

Continued on page 2

# Physical Characteristics

Species: human

Source: E. coli expression

Quantity: 100 µg

Concentration: 1 mg/ml

Formulation: 50 mM HEPES pH 7.5. 150 mM sodium chloride, 2 mM dithiothreitol, 10% glycerol

Molecular Weight: ~20 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

Stability/Storage: 12 months at -70°C;

aliquot as required

## **Protein Sequence:**

MYHHHHHSSGLEVLFQGPLGSSTPAR RRLMRDFKRLQEDPPVGVSGAPSEN | NIMQWNAVIFGPEGTPFEDGT FKLVIEFSEEYPNKPPTVRFLSKM FHPNVYADGSICLDILQNRWSPTYD VSSILTSIQSLLDEPNPNSPANSQAAQ LYQENKREYEKRVSAIVEQSWNDS

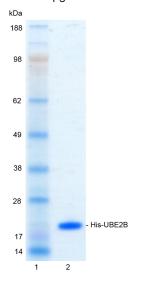
Tag (bold text): N-terminal His Protease cleavage site: PreScission™ (LEVLFQ ▼ GP) UBE2B (regular text): Start bold italics (amino acid residues

Accession number: NP\_003328

# **Quality Assurance**

### **Purity:**

4-12% gradient SDS-PAGE InstantBlue™ staining Lane 1: MW markers Lane 2: 1 µg His-UBE2B



### **Protein Identification:**

Confirmed by mass spectrometry.

## **E2-Ubiquitin Thioester Loading Assay:**

The activity of His-UBE2B was validated by loading E1 UBE1 activated ubiquitin onto the active cysteine of the His-UBE2B E2 enzyme via a transthiolation reaction. Incubation of the UBE1 and His-UBE2B enzymes in the presence of ubiquitin and ATP at 30°C was compared at two time points,  $T_0$  and  $T_{10}$  minutes. Sensitivity of the ubiquitin/His-UBE2B thioester bond to the reducing agent DTT was confirmed.



## **ORDERS / SALES SUPPORT**

International: +1-617-245-0003 US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233)

Email: sales.support@ubiquigent.com

## **UK HQ and TECHNICAL SUPPORT**

International: +44 (0) 1382 381147 (9AM-5PM UTC) US/Canada: +1-617-245-0020 (9AM-5PM UTC) Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© **Ubiquigent 2011**. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0

# UBE2B (HR6B) [6His-tagged]

# E2 - Ubiquitin Conjugating Enzyme

Alternate Names: HHR6B, HR6B, RAD6B, Ubiquitin carrier protein B,

Ubiquitin protein ligase B

**Cat. No. 62-0072-100** Quantity: 100 μg **Lot. No. 1823** Storage: -70°C

FOR RESEARCH USE ONLY NOT FOR USE IN HUMANS



**CERTIFICATE OF ANALYSIS Page 2 of 2** 

# **Background**

Continued from page 1

#### References:

Escalier D, Bai XY, Silvius D, Xu PX, Xu X (2003) Spermatid nuclear and sperm periaxonemal anomalies in the mouse Ube2b null mutant. *Mol Reprod Dev* **65**, 298-308.

Koken MH, Reynolds P, Jaspers-Dekker I, Prakash L, Prakash S, Bootsma D, Hoeijmakers JH (1991) Structural and functional conservation of two human homologs of the yeast DNA repair gene RAD6. *Proc Natl Acad Sci U S A* **88**, 8865-9.

Notenboom V, Hibbert RG, van Rossum-Fikkert SE, Olsen JV, Mann M, Sixma TK (2007) Functional characterization of Rad18 domains for Rad6, ubiquitin, DNA binding and PCNA modification. *Nucleic Acids Res* **35**, 5819-30.

Suryavathi V, Khattri A, et al. (2008) Novel variants in UBE2B gene and idiopathic male infertility. J Androl 29, 564-71.

Tsuji Y, Watanabe K, et al. (2008) Recognition of forked and single-stranded DNA structures by human RAD18 complexed with RAD6B protein triggers its recruitment to stalled replication forks. *Genes Cells* 13, 343-54.

Ulrich HD, Jentsch S (2000) Two RING finger proteins mediate cooperation between ubiquitin-conjugating enzymes in DNA repair. *EMBO J* **19**, 3388-97.



Dundee, Scotland, UK

## **ORDERS / SALES SUPPORT**

International: +1-617-245-0003

US Toll-Free: 1-888-4E1E2E3 (1-888-431-3233) Email: sales.support@ubiquigent.com

## **UK HQ and TECHNICAL SUPPORT**

| International: +44 (0) 1382 381147 (9AM-5PM UTC) | US/Canada: +1-617-245-0020 (9AM-5PM UTC) | Email: tech.support@ubiquigent.com

Email services@ubiquigent.com for enquiries regarding compound profiling and/or custom assay development services.

© Ubiquigent 2011. Unless otherwise noted, Ubiquigent, Ubiquigent logo and all other trademarks are the property of Ubiquigent, Ltd.

Limited Terms of Use: For research use only. Not for use in humans or for diagnostics. Not for distribution or resale in any form, modification or derivative OR for use in providing services to a third party (e.g. screening or profiling) without the written permission of Ubiquigent, Ltd.

Lot-specific COA version tracker: v1.0.0