

UBE2R2 (CDC34B) [6His-tagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: CDC34B, EC 6.3.2.19, FLJ20419, MGC10481, UBC3B

Cat. No. 62-0094-020

Lot. No. 2121

Quantity: 20 µg

Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 1

Background

The enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including the 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). UBE2R2 is a member of the E2 conjugating enzyme family and cloning of the human gene was first described by Semplici *et al.* (2002). Site directed mutagenesis studies have shown that serine 233 in the C-terminal domain of UBE2R2 is the site at which CK2-dependent phosphorylation occurs (Semplici *et al.*, 2002). *In vitro* binding experiments have also demonstrated that phosphorylated UBE2R2 and UBE2R1 bind specifically to the F-box protein beta-TRCP, which results in enhanced degradation of beta-catenin (a substrate of the Beta Transducin Repeat Containing protein (BTRC) (Semplici *et al.*, 2002).

References:

Semplici F, Meggio F, Pinna LA, Oliviero S (2002) CK2-dependent phosphorylation of the E2 ubiquitin conjugating enzyme UBC3B induces its interaction with beta-TrCP and enhances beta-catenin degradation. *Oncogene* 21, 3978-87.

Physical Characteristics

Species: human

Source: *E. coli* expression

Quantity: 20 µg

Concentration: 1 mg/ml

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

Molecular Weight: ~31 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

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

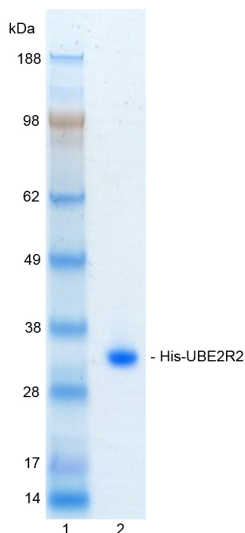
Protein Sequence:

M G S S H H H H H S S G L V P R G S H M A S M T G
G Q Q M G R G S M A Q Q Q M T S S Q K A L M L E L K
S L Q E E P V E G F R I T L V D E S D L Y N W E V A I F G P
P N T L Y E G G Y F K A H I K F P I D Y P S P T F R F L T K
M W H P N I Y E N G D V C I S I L H P P V D D P Q S G E L P S
E R W N P T Q N V R T I L L S V I S L L N E P N T F S P A N V
D A S V M F R K R W R D S K G K D K E Y A E I I R K Q V S A T
K A E A E K D G V K V P T T L A E Y C I K T K V P S N D N S
S D L L Y D D L Y D D D I D D E D E E E E D A D C Y D D D D S
G N E E S

Tag (**bold text**): N-terminal His
Protease cleavage site: PreScission™ (LVPR▼GS)
UBE2R2 (regular text): Start **bold italics** (amino acid residues 1-238)
Accession number: AAH04862

Quality Assurance

Purity: 4-12% gradient SDS-PAGE InstantBlue™ staining
Lane 1: MW markers
Lane 2: 1 µg His-UBE2R2



Protein Identification:

Confirmed by mass spectrometry.

E2-Ubiquitin Thioester Loading Assay:

The activity of His-UBE2R2 was validated by loading E1 UBE1 activated ubiquitin onto the active cysteine of the His-UBE2R2 E2 enzyme via a transthiolation reaction. Incubation of the UBE1 and His-UBE2R2 enzymes in the presence of ubiquitin and ATP at 30°C was compared at two time points, T₀ and T₁₀ minutes. Sensitivity of the ubiquitin/His-UBE2R2 thioester bond to the reducing agent DTT was confirmed.



www.ubiquigent.com
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