

UBE2V2 (Mms2) [6His-tagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: UEV2, Methyl Methanesulphonate Sensitive 2, Mms2, Enterocyte Differentiation-Promoting Factor 1, EDPF1

Cat. No. 62-0099-020
Lot. No. 30039

Quantity: 20 µg
Storage: -70°C

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). UBE2V2 is a member of the E2 conjugating enzyme family and cloning of the human gene was first described by Fritsche *et al.* (1997). UBE2V2 shares 90% sequence identity with UBE2V1 in its C-terminal domain (Sancho *et al.*, 1998). The UEV protein Mms2 (yeast homologue of human UBE2V2) forms a heterodimer with yeast Ubc13 (UBE2N) which is recruited to chromatin by the RING finger proteins RAD5 and RAD18 in the RAD6 dependent post-replicative DNA repair pathway (Hofmann and Pickart 1999). These proteins also play a central role in the assembly of K63-linked polyubiquitin chains (Ulrich and Jentsch 2000; Xiao *et al.*, 1998). UEV/Ubc complexes have been implicated in the assembly of Lys63-linked polyubiquitin chains that act as a novel signal in post-replicative DNA repair and IκBα kinase activation. Recent crystal structure analysis provides direct evidence that the Mms2/Ubc13 heterodimer is necessary for DNA repair (Morales *et al.*, 2001).

Continued on page 2

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: ~18 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

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

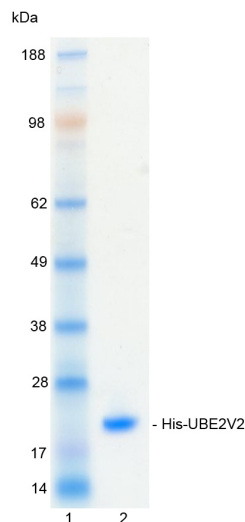
Protein Sequence:

MGSSHHHHHSSGL~~LEVL~~FQGP~~GS~~MAVST
GVKVPNRNFRLLLEEELEGQKGVGDGTVSWGLED
DEDMLTRWTGMIIGPPRTNYENRIYSLKVECG
PKYPEAPPSVRFVTKINMNGINSSGMVDAR
SIPVLAKWQNSYSIKVVLQELRRLMMSKENM
KLPQPPEGQTYNN

Tag (**bold text**): N-terminal His
Protease cleavage site: PreScission™ (~~LEVL~~FQ▼GP)
UBE2V2 (regular text): Start **bold italics** (amino acid residues 1-145)
Accession number: NP_003341.1

Quality Assurance

Purity: 4-12% gradient SDS-PAGE InstantBlue™ staining
Lane 1: MW markers
Lane 2: 1 µg His-UBE2V2



Protein Identification:

Confirmed by mass spectrometry.

Polyubiquitin Chain Formation Assay:

The activity of His-UBE2V2 was validated in a polyubiquitin chain formation assay. Incubation of UBE1, UBE2N and CHIP - with and without His-UBE2V2 - in the presence of ubiquitin and ATP at 30°C was compared at two time points, T₀ and T₆₀ minutes. Polyubiquitin chains were detected by Western blot using a monoclonal anti-ubiquitin conjugate antibody. The ability of His-UBE2V2 to promote the formation of polyubiquitin chains was observed.



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

UBE2V2 (Mms2) [6His-tagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: UEV2, Methyl Methanesulphonate Sensitive 2, Mms2, Enterocyte Differentiation-Promoting Factor 1, EDPF1

Cat. No. 62-0099-020

Lot. No. 30039

Quantity: 20 µg

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:

Fritsche, J., M. Rehli, *et al.* (1997). Molecular cloning of a 1alpha,25-dihydroxyvitamin D3-inducible transcript (DDVit 1) in human blood monocytes. *Biochem Biophys Res Commun* **235**(2): 407-12.

Hofmann RM, Pickart CM (1999). Noncanonical MMS2-encoded ubiquitin-conjugating enzyme functions in assembly of novel polyubiquitin chains for DNA repair. *Cell* **96**, 645-53.

Moraes, T. F., R. A. Edwards, *et al.* (2001). Crystal structure of the human ubiquitin conjugating enzyme complex, hMms2-hUbc13. *Nat Struct Biol* **8**(8): 669-73.

Sancho, E., M. R. Vila, *et al.* (1998). Role of UEV-1, an inactive variant of the E2 ubiquitin-conjugating enzymes, in *in vitro* differentiation and cell cycle behavior of HT-29-M6 intestinal mucosecretory cells. *Mol Cell Biol* **18**(1): 576-89.

Ulrich, H. D. and S. Jentsch (2000). Two RING finger proteins mediate cooperation between ubiquitin-conjugating enzymes in DNA repair. *EMBO J* **19**(13): 3388-97.

Xiao, W., S. L. Lin, *et al.* (1998). The products of the yeast MMS2 and two human homologs (hMMS2 and CROC-1) define a structurally and functionally conserved Ubc-like protein family. *Nucleic Acids Res* **26**(17): 3908-14.



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