

UBE2L6 (UbcH8) [untagged]

E2 - Ubiquitin or ISG15 Conjugating Enzyme

Alternate Names: MGC40331, Retinoic acid induced gene B protein, RIG-B, UbcH8, Ubiquitin conjugating enzyme UbcH8

Cat. No. 62-0044-100
Lot. No. 1471

Quantity: 100 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 ubiquitination; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). UBE2L6 is a member of the E2 conjugating enzyme family, identification and cloning of the gene from a yeast 2 hybrid screen was first described by Kumar *et al.* (1997). Human UBE2L6 has been mapped to chromosome 11q12 and shares 46% homology with UBE2L3 (Recombinant UBE2L6 forms thiol ester complexes with ubiquitin *in vitro* and is able to transfer ubiquitin to E6AP (Ardley *et al.*, 2000). UBE2L6 has also been shown to conjugate the ubiquitin like protein (Ubl) ISG15. *In vitro* assays and RNA interference have shown that of the ubiquitin E2 enzymes tested UBE2L6 is a major E2 enzyme for ISG15 conjugation (Li *et al.*, 2006; Zhao *et al.*, 2004). The nuclear matrix protein Msx2-interacting nuclear target protein (MINT) regulates the expression of key transcriptional effectors in diverse signalling pathways and has been identified as a binding partner of UBE2L6 by co-immunoprecipitation and GST pull down (Li *et al.*, 2006). Degradation of the chromosomal translocation products AML1-ETO and PML-RAR alpha which contribute to the pathogenesis of leukaemias is mediated by the E2 en-

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: ~18 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

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

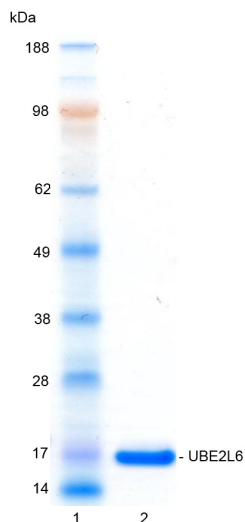
Protein Sequence:

GPGSMMASMRVVKELEDLQKKPPPYLRNLSSD
DANVLVWHALLLPDQPPYHLKAFNLRI SFPEYP
FKPPMIKF~~TT~~KIYHPNVDENGQICLP IISSEN
WKPC~~TK~~T~~CQ~~VLEALNVLVNRPNIREPLRMD
LADLLTQNP~~EL~~FRKNAEEFTLRF~~G~~VDRPS

The residues underlined remain after cleavage and removal of the purification tag.
UBE2L6 (regular text): Start **bold italics** (amino acid residues 1-153)
Accession number: NP_004214

Quality Assurance

Purity: 4-12% gradient SDS-PAGE InstantBlue™ staining
Lane 1: MW markers
Lane 2: 1 µg UBE2L6



Protein Identification:

Confirmed by mass spectrometry.

E2-Ubiquitin Thioester Loading Assay:

The activity of UBE2L6 was tested by loading E1 UBE1 activated ubiquitin onto the active cysteine of the UBE2L6 E2 enzyme via a transthioester reaction. Incubation of the UBE1 and UBE2L6 enzymes in the presence of ubiquitin and ATP at 30°C was compared at two time points, T₀ and T₁₀ minutes. Under these conditions tested no UBE2L6 ubiquitin thioester loading 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

UBE2L6 (UbcH8) [untagged]

E2 - Ubiquitin or ISG15 Conjugating Enzyme

Alternate Names: MGC40331, Retinoic acid induced gene B protein, RIG-B, UbcH8, Ubiquitin conjugating enzyme UbcH8

Cat. No. 62-0044-100
Lot. No. 1471

Quantity: 100 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

zyme UBE2L6 and the E3 ligase SIAH-1. Thus, UBE2L6 could be a potential therapeutic target in the treatment of leukaemia (Kramer *et al.*, 2008).

References:

Ardley HC, Rose SA, Tan N, Leek JP, Markham AF, Robinson PA (2000) Genomic organization of the human ubiquitin-conjugating enzyme gene, UBE2L6 on chromosome 11q12. *Cytogenet Cell Genet* **89**, 137-40.

Kramer OH, Muller S, Buchwald M, Reichardt S, Heinzel T (2008) Mechanism for ubiquitylation of the leukemia fusion proteins AML1-ETO and PML-RARalpha. *FASEB J* **22**, 1369-79.

Kumar S, Kao WH, Howley PM (1997) Physical interaction between specific E2 and Hect E3 enzymes determines functional cooperativity. *J Biol Chem* **272**, 13548-54.

Li J, Wang J, Yang X, Qin H, Dong X, Zhu Y, Liang L, Liang Y, Han H (2006) The Spen homolog Msx2-interacting nuclear target protein interacts with the E2 ubiquitin-conjugating enzyme UbcH8. *Mol Cell Biochem* **288**, 151-7.

Zhao C, Beaudenon SL, Kelley ML, Waddell MB, Yuan W, Schulman BA, Huibregtse JM, Krug RM (2004) The UbcH8 ubiquitin E2 enzyme is also the E2 enzyme for ISG15, an IFN-alpha/beta-induced ubiquitin-like protein. *Proc Natl Acad Sci U S A* **101**, 7578-82.



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