

UBE2D1 (UbcH5a) [untagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: E2(17)KB 1, EC 6.3.2.19, SFT, Stimulator of Fe transport, homolog of UBC4/5, UbcH5, UbcH5A, Ubiquitin protein ligase, Ubiquitin-conjugating enzyme E2-17 kDa 1, Ubiquitin-conjugating enzyme UbcH5A

Cat. No. 62-0010-100

Lot. No. 1458

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 ubiquitylation; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). UBE2D1 is a member of the E2 ubiquitin-conjugating enzyme family and cloning of the human gene was first described by Scheffner *et al.* (1994). UBE2D1 shares 89% sequence identity with its Drosophila homologue and mediates E6/UBE3A (E6AP)-induced ubiquitylation of p53 (Jensen *et al.*, 1995; Scheffner *et al.*, 1994). Ubiquitylation of the yeast PTS1 import receptor (pex5p) has been demonstrated in an *in vitro* assay in the presence of the human UBE2D1 in combination with the ring domain of the yeast E3 ligase pex10p (Williams *et al.*, 2008). Sequence encoding the stimulated Iron transport gene SFT overlaps with intron 7 and exon 6 of UBE2D1, and RT/PCR has shown significantly upregulated levels of UBE2D1 in livers of iron-overloaded patients with hereditary hemochromatosis (Gehrke *et al.*, 2003).

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:

GSHMASMTGGQQMGRGSALKRIQKELSDLQRD
PPAHCSAGPVGDDLFHWQATIMGPPDSAYQG
GVFFLTVHFPTDYPFKPKIAFTTKIYHPNIN
SNGSICLDILRSQWSPALTVSKVLLSICSL
CDPNPDDPLVPDIAQIYKSDKEKYNRHARE
WTQKYAM

The residues underlined remain after cleavage and removal of the purification tag.

UBE2D1 (regular text): Start **bold italics** (amino acid residues 2-147)

Accession number: NP_003329

Quality Assurance

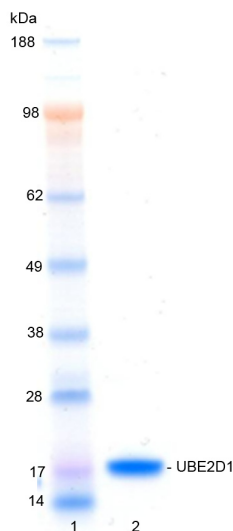
Purity:

4-12% gradient SDS-PAGE

InstantBlue™ staining

Lane 1: MW markers

Lane 2: 1 µg UBE2D1



Protein Identification:

Confirmed by mass spectrometry.

E2-Ubiquitin Thioester Loading Assay:

The activity of UBE2D1 was validated by loading E1 UBE1 activated ubiquitin onto the active cysteine of the UBE2D1 E2 enzyme via a transthioylation reaction. Incubation of the UBE1 and UBE2D1 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/UBE2D1 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

UBE2D1 (UbcH5a) [untagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: E2(17)KB 1, EC 6.3.2.19, SFT, Stimulator of Fe transport, homolog of UBC4/5, UbcH5, UbcH5A, Ubiquitin protein ligase, Ubiquitin-conjugating enzyme E2-17 kDa 1, Ubiquitin-conjugating enzyme UbcH5A

Cat. No. 62-0010-100

Lot. No. 1458

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

References:

Gehrke SG, Riedel HD, Herrmann T, Hadaschik B, Bents K, Veltkamp C, Stremmel W (2003) UbcH5A, a member of human E2 ubiquitin-conjugating enzymes, is closely related to SFT, a stimulator of iron transport, and is up-regulated in hereditary hemochromatosis. *Blood* **101**, 3288-93.

Jensen JP, Bates PW, Yang M, Vierstra RD, Weissman AM (1995) Identification of a family of closely related human ubiquitin conjugating enzymes. *J Biol Chem* **270**, 30408-14.

Scheffner M, Huibregtse JM, Howley PM (1994) Identification of a human ubiquitin-conjugating enzyme that mediates the E6-AP-dependent ubiquitination of p53. *Proc Natl Acad Sci USA* **91**, 8797-801.

Williams C, van den Berg M, Geers E, Distel B (2008) Pex10p functions as an E3 ligase for the Ubc4p-dependent ubiquitination of Pex5p. *Biochem Biophys Res Commun* **374**, 620-4.



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