

UBE2E3 (UbcH9) [GST-tagged]

E2 – Ubiquitin Conjugating Enzyme

Alternate Names: UbcH9, UbcM2, Ubiquitin conjugating enzyme UbcH9

Cat. No. 62-0021-020
Lot. No. 1392

Quantity: 20 µg
Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS

Background

The enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including regulated and targeted proteosomal 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). UBE2E3 is a member of the E2 ubiquitin-conjugating enzyme family and cloning of the gene was first described by Ito *et al.* (1999). UBE2E3 binds to the RING-finger proteins ARA54 and RNF8, thought to act as E3 ligases in the ubiquitylation of nuclear proteins (Ito *et al.*, 2001). The epithelial Na⁺ channel (ENaC) is regulated by UBE2E3 and the E3 ligase NEDD4.2. UBE2E3 interacts with NEDD4.2 via its UBC domain and ubiquitylation of ENaC occurs by NEDD4.2 binding the PY motifs of its α , β and γ subunits (Debonneville and Staub, 2004). NEDD4.2 is a negative regulator of ENaC and deletions in the PY motifs of the β and γ subunits of ENaC cause Liddle's syndrome, an inherited form of hypertension. The loss of NEDD4.2 binding sites in mutated ENaC causes an increase in channel number at the cell surface and increased Na⁺ reabsorption by the distal nephron, resulting in hypertension (Abriel *et al.*, 1999).

References:

Abriel H, Loffing J, Rebhun JF, Pratt JH, Schild L, Horisberger JD, Rotin D, Staub O (1999) Defective regulation of the epithelial Na⁺ channel by Nedd4 in Liddle's syndrome. *J Clin Invest* **103**, 667-73.

Debonneville C, Staub O (2004) Participation of the ubiquitin-conjugating enzyme UBE2E3 in Nedd4-2-dependent regulation of the epithelial Na⁺ channel. *Mol Cell Biol* **24**, 2397-409.

Ito K, Adachi S, Iwakami R, Yasuda H, Muto Y, Seki N, Okano Y (2001) N-Terminally extended human ubiquitin-conjugating enzymes (E2s) mediate the ubiquitination of RING-finger proteins, ARA54 and RNF8. *Eur J Biochem* **268**, 2725-32.

Ito K, Kato S, Matsuda Y, Kimura M, Okano Y (1999) cDNA cloning, characterization, and chromosome mapping of UBE2E3 (alias UbcH9), encoding an N-terminally extended human ubiquitin-conjugating enzyme. *Cytogenet Cell Genet* **84**, 99-104.

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: ~50 kDa

Purity: >90% by InstantBlue™ SDS-PAGE

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

Protein Sequence:

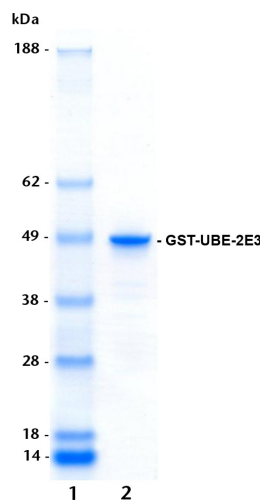
MSPILGYWKIKGLVQPTRLLEYLEEKYEEH
LYERDEGDKWRNKKFELGLEFPNLPYYIDG
VKLTQSMAIIRYIADKHNMLGGCPKER
AEISMLEGAVLDIRYGVSRAYSKDFETLKVD
FLSKLPEMLKMFEDRLCHKTYLNGDHTHP
DFMLYDALDVVLYMDPMCLDAFPKLVCFK
KRIEAIPIQIDKYLKSSKYIAWPLQGWQATFG
GGDHPKSDLEVLFFQGPLGSMSSDRQRSD
DESPSTSSGSSDADQRDPAAPEPEEQEERKP
SATQQKKNTKLSSKTTAKLSTSAKRIQKE
LAEITLDPNCSAGPKGDNIYWRSTILGPPGS
VYEGGVFLDITFSSDYFPKPPKVTFRTRIYHCNIN
SQGVICLDILKDNWSPALTISKVLLSICSLTDCN
PADPLVGSATQYLTRAEHDRIARQWTKRYAT

Tag (**bold text**): N-terminal glutathione-S-transferase (GST)
Protease cleavage site: PreScission™ (LEVLFOVGP)
UBE2E3 (regular text): Start **bold italics** (amino acid residues 1-207)
Accession number: NP_006348.1

Quality Assurance

Purity:

4-12% gradient SDS-PAGE
InstantBlue™ staining
lane 1: MW markers
lane 2: 1 µg GST-UBE2E3



Protein Identification:

Confirmed by mass spectrometry.

E2-Ubiquitin Thioester Loading Assay:

The activity of GST-UBE2E3 was validated by loading E1 UBE1 activated ubiquitin onto the active cysteine of the GST-UBE2E3 E2 enzyme via a transthiolation reaction. Incubation of the UBE1 and GST-UBE2E3 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/GST-UBE2E3 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