

# UBE2G1 (Ubc7) [GST-tagged]

## E2 – Ubiquitin Conjugating Enzyme

Alternate Names: E217K, UBC7, UBE2G

Cat. No. **62-0027-100**  
Lot. No. **1394**

Quantity: 100 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 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). UBE2G1 is a member of the E2 conjugating enzyme family and cloning of the human gene was first described by Watanabe *et al.* (1996). UBE2G1 shares 74% sequence identity with UBC7 from *C. elegans* and a high degree of homology with UBC7 from other species. Expression of UBE2G1 and a helix-loop-helix transcription factor and member of the MYC/MAX superfamily (ROX/MNT) is decreased in medullablastoma tumours. Haploinsufficiency of the human 17p13.3 region is associated with 35% to 50% of medullablastomas, indicating the presence of one or more tumour suppressor genes which have not yet been identified (Cvekl *et al.*, 2004).

### References:

Cvekl A, Jr., Zavadii J, Birshstein BK, Grotzer MA, Cvekl A (2004) Analysis of transcripts from 17p13.3 in medulloblastoma suggests ROX/MNT as a potential tumour suppressor gene. *Eur J Cancer* **40**, 2525-32.

Watanabe TK, Kawai A, Fujiwara T, Maekawa H, Hirai Y, Nakamura Y, Takahashi E (1996) Molecular cloning of UBE2G, encoding a human skeletal muscle-specific ubiquitin-conjugating enzyme homologous to UBC7 of *C. elegans*. *Cytogenet Cell Genet* **74**, 146-8.

### 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:** ~46 kDa

**Purity:** >98% by InstantBlue™ SDS-PAGE

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

### Protein Sequence:

**MSPILGYWKIKGLVQPTRLLEYLEEKYEEH  
LYERDEGDKWRNKKFELGLEFPNLPYYIDGD  
VKLTQSMAIRYIADKHNMLGGCPKERAEISM  
LEGAVLDIRYGVSR IAYS KDFETLKVDFL  
SKLPEMLKMFEDRLCHKTYLNGDHSVTHPD  
FMLYDALDVVL YMDP MCLDAFPKLVCFK  
KR IEAIPQIDKYLKSSKYIAWPLQGWQATFG  
GGDHPKSDLEVLVFOGPLGSMTELOSALLLR  
RQLAELNKNPVEGFSAGLIDDNDLYRWEVLI  
IGPPDTLYEGGVFKAHLTFPKDYPLRPPKM  
KFITEIWHPNVDKNGDVCISILHEPGEDKY  
GYEKPEERWLPHTVETIMISVISMLADPNGD  
SPANVDAAKEWREDRNGEFKRKVARCVRK  
SQETA FE**

Tag (**bold text**): N-terminal GST

Protease cleavage site: PreScission™ (LEVLVFOG↓GP)

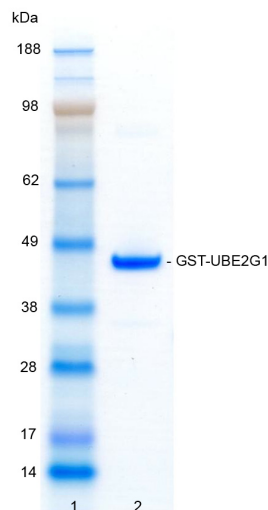
UBE2G1 (regular text): Start **bold italics** (amino acid residues 1-170)

Accession number: NP\_003333

### Quality Assurance

#### Purity:

4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 1 µg GST-UBE2G1



#### Protein Identification:

Confirmed by mass spectrometry.

#### E2-Ubiquitin Thioester Loading Assay:

The activity of GST-UBE2G1 was validated by loading E1 UBE1 activated ubiquitin onto the active cysteine of the GST-UBE2G1 E2 enzyme via a transthioylation reaction. Incubation of the UBE1 and GST-UBE2G1 enzymes in the presence of ubiquitin and ATP at 30°C was compared at two time points, T<sub>0</sub> and T<sub>10</sub> minutes. The sensitivity of this ubiquitin/GST-UBE2G1 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