

# USP2 (isoform 4) [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: 41 kDa ubiquitin specific protease, UBP41

Cat. No. 64-0014-050

Lot. No. 30013

Quantity: 50 µg

Storage: -70°C



FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS Page 1 of 2

## Background

Deconjugating enzymes (DCEs) are proteases that process ubiquitin or ubiquitin-like gene products, reverse the modification of proteins by a single ubiquitin or ubiquitin-like protein (UBL) and remodel polyubiquitin (or poly-UBL) chains on target proteins (Reyes-Turcu *et al.*, 2009). The deubiquitylating – or deubiquitinating – enzymes (DUBs) represent the largest family of DCEs and regulate ubiquitin dependent signaling pathways. The activities of the DUBs include the generation of free ubiquitin from precursor molecules, the recycling of ubiquitin following substrate degradation to maintain cellular ubiquitin homeostasis and the removal of ubiquitin or ubiquitin-like proteins (UBL) modifications through chain editing to rescue proteins from proteasomal degradation or to influence cell signalling events (Komander *et al.*, 2009). There are two main classes of DUB; cysteine proteases and metalloproteases. Ubiquitin specific protease 2 (USP2) is a member of the cysteine protease enzyme family and cloning of the human gene was first described by Baek *et al.* (1997). USP2 influences the nuclear translocation of NF-κB transcription factors in response to inflammation, injury and other environmental changes. In resting cells, NF-κB is bound to the inhibitory protein IκBα, and therefore maintained in an inactive state in the cytoplasm. The proteasomal degradation of IκBα is prevented by USP2 deubiquitylase activity. This regulates the expression of genes that are involved in immune responses, cell proliferation,

Continued on page 2

## Physical Characteristics

**Species:** human

**Source:** *E. coli*

**Quantity:** 50 µg

**Concentration:** 0.5 mg/ml

**Formulation:** 50 mM HEPES pH 7.5, 150 mM sodium chloride, 2 mM dithiothreitol, 10% glycerol

**Molecular Weight:** ~73 kDa

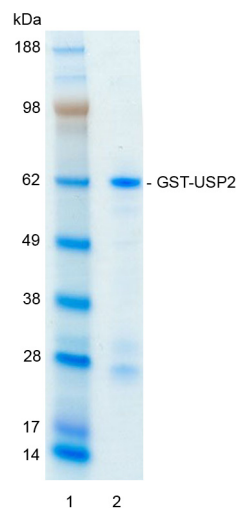
**Purity:** >68% by InstantBlue™ SDS-PAGE

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

**Protein Sequence:** Please see page 2

## Quality Assurance

**Purity:**  
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
lane 1: MW markers  
lane 2: 1 µg GST-USP2



**Protein Identification:**  
Confirmed by mass spectrometry.

### Deubiquitylase Enzyme Assay:

The activity of GST-USP2 was validated by determining the increase in fluorescence measured as a result of the enzyme catalysed cleavage of the fluorogenic substrate Ubiquitin-Rhodamine110-Glycine generating Ubiquitin and Rhodamine110-Glycine. Incubation of the substrate in the presence or absence of GST-USP2 was compared confirming the deubiquitylating activity of GST-USP2.



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 2012. 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

# USP2 (isoform 4) [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: 41 kDa ubiquitin specific protease, UBP41

Cat. No. 64-0014-050

Lot. No. 30013

Quantity: 50 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

differentiation and apoptosis (Metzig *et al.*, 2011). Deregulation of NF-κB activity is involved in the pathology of many diseases including chronic inflammation and cancer. USP2 expression itself is frequently downregulated in breast carcinomas (Metzig *et al.*, 2011). USP2 deubiquitylase activity induces cell death via the apoptosis inducing factor (AIF). AIF is a mitochondrial oxidoreductase that becomes truncated when the cell is under stress and the truncated AIF (tAIF) becomes translocated into the nucleus, and induces caspase-independent cell death. USP2 deubiquitylates and stabilizes tAIF, thus promoting AIF-mediated cell death. In contrast, the E3 ligase CHIP ubiquitylates and destabilizes tAIF, thus preventing cell death (Oh *et al.*, 2011).

## References:

Baek SH, Choi KS, Yoo YJ, Cho JM, Baker RT, Tanaka K, Chung CH (1997) Molecular cloning of a novel ubiquitin-specific protease, UBP41, with isopeptidase activity in chick skeletal muscle. *J Biol Chem* 272, 25560-25565.

Komander D, Clague MJ, Urbe S (2009) Breaking the chains: structure and function of the deubiquitinases. *Nat Rev Mol Cell Biol* 10, 550-563.

Metzig M, Nickles D, Falschlehner C, Lehmann-Koch J, Straub BK, Roth W, Boutros M (2011) An RNAi screen identifies USP2 as a factor required for TNF-alpha-induced NF-kappaB signaling. *Int J Cancer* 129, 607-618.

Oh KH, Yang SW, Park JM, Seol JH, Iemura S, Natsume T, Murata S, Tanaka K, Jeon YJ, Chung CH (2011) Control of AIF-mediated cell death by antagonistic functions of CHIP ubiquitin E3 ligase and USP2 deubiquitinating enzyme. *Cell Death Differ* 18, 1326-1336.

Reyes-Turcu FE, Ventii KH, Wilkinson KD (2009) Regulation and cellular roles of ubiquitin-specific deubiquitinating enzymes. *Ann Rev Biochem* 78, 363-397.

## Physical Characteristics

Continued from page 1

### Protein Sequence:

**MSPILGYWKIKGLVQPTRLLEYLEEKYEELH  
LYERDEGDKWRNKKFELGLEFPNLPYYIDGD  
VKLTQSMAIIRYIADKHNMLGGCPKERAEISM  
LEGAVLDIRYGVSR IAYS KDFETLKVDFL  
SKLPEMLKMFEDRLCHKTYLNGDHVTHPD  
FMLYDALDVVLYMDPMCLDAFPKLVCFK  
KR IEAIPQIDKYLKSSKYIAWPLQGWQATFG  
GGDHPPKSDLEVL FQGPLGSPNSRVD MRTSYT  
VTLPEPPAAPFPALAKELRPRSPLSPSLLL  
STFVGLLLNKAKNSKSAQGLAGLRNLGNTCF  
MNSILQCLSNTR ELRDYCLQRLYMRDLHHGS  
NAHTALVEEFAKLIQTIWTSSPNDVVSPSEFK  
TQIQRYAPRFVGYNQDQAEFLRFLLDGLH  
NEVNRVTLRPKSNPENLDHLPDDEKGRQM  
WRKYLEREDSRIGDLFVGLKSSLTCTDC  
GYCSTVDFDPFDLSLPIAKRGYPEVTLMDC  
MRLFTKEDVLDGDEKPTCCRCRGRKRCIK  
KFSIQRFPKILVHLKRFSESRI RTSKLT  
TFVNFPLRDLDLREFASENTNHA VYNYLAVS  
NHS GTTMGGHYTAYCRSPGTGEWHTFNDSS  
VTPMSSSQVRTSDAYLLFYELASPPSRM**

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

Protease cleavage site: PreScission™ (LEVL FQ▼GP)

USP2 (regular text): Start **bold italics** (amino acid residues 1-396)

Accession number: NP\_741994

UniProt number: O75604-4



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 2012. 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