

# USP4 [6His-tagged]

## Deconjugating enzyme: Deubiquitylase

**Alternate Names:** Deubiquitinating enzyme 4, Ubiquitin carboxyl-terminal hydrolase 4, Ubiquitin-specific-processing protease 4, Ubiquitin thioesterase 4

**Cat. No.** 64-0001-050  
**Lot. No.** 30031

**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 signalling 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 carboxyl-terminal hydrolase 4 (Ubiquitin Specific Protease 4; USP4) is a member of the cysteine protease enzyme family and cloning of the human gene was first described by Gupta *et al.* (1993). In 1995, USP4 was identified as a proto-oncogene related to USP6, showing a consistently elevated gene expression level in small cell tumours and lung adenocarcinomas suggesting that it may have a possible causative role in neoplasia (Gray *et al.*, 1995). USP4 has been implicated in a number of other processes, including protein quality control in the endoplasmic

Continued on page 2

### Physical Characteristics

**Species:** human

**Protein Sequence:** Please see page 2

**Source:** *E. coli* expression

**Quantity:** 50 µg

**Concentration:** 0.5 mg/ml

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

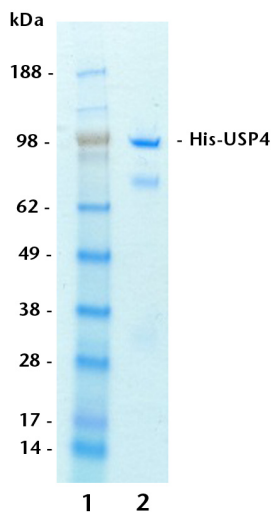
**Molecular Weight:** ~111 kDa

**Purity:** >76% by InstantBlue™ SDS-PAGE

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

### Quality Assurance

**Purity:**  
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 1 µg His-USP4



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

**Deubiquitylating Enzyme Assay:**  
The activity of His-USP4 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 His-USP4 was compared confirming the deubiquitylating activity of His-USP4.



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](mailto: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

# USP4 [6His-tagged]

## Deconjugating enzyme: Deubiquitylase

**Alternate Name:** Deubiquitinating enzyme 4, Ubiquitin carboxyl-terminal hydrolase 4, Ubiquitin-specific-processing protease 4, Ubiquitin thioesterase 4

**Cat. No.** 64-0001-050  
**Lot. No.** 30031

**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

reticulum and p53 and Wnt signalling. USP4 has also been reported to inhibit the kinase TAK1 that is ubiquitylated by the AKT regulator TRAF6 (Uras *et al.*, 2012).

#### References:

Gray DA, Inazawa J, Gupta K, Wong A, Ueda R and Takahashi T (1995) Elevated expression of Unph, a proto-oncogene at 3p21.3, in human lung tumors. *Oncogene* **10**, 2179-2183.

Gupta K, Copeland NG, Gilbert DJ, Jenkins NA and Gray DA (1993) Unp, a mouse gene related to the tre oncogene. *Oncogene* **8**, 2307-2310.

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

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

Uras IZ, List T and Nijman SM (2012) Ubiquitin-specific protease 4 inhibits mono-ubiquitination of the master growth factor signaling kinase PDK1. *PLoS one* **7**, e31003.

### Physical Characteristics

Continued from page 1

#### Protein Sequence:

**MGSSHHHHHSSGLEVLFQGP**GSMAEGGGCR  
ERPDAETQKSELGPLMRTTLQGAQWYLID  
SRWFKQWKKYVGFDSWDMYVGEHNLFP  
PIDNSGLFSDPESQTLKEHLIDELDYVLP  
TEAWNKLNLNWCVEGQQPIVRKVVEH  
GLFVKHCKVEVYLLLELKLKENSPTNVLSCHF  
SKADTIATIEKEMRKLFNIPAEERTRL  
WNKYMSNTYEQLSKLDNTVQDAGLYQGQV  
LVIEPQNE DGTWPRQTLQSKSSTAPSRN  
FTTSPKSSASPYSSVSASLIANGDSTSTCG  
MHSSGVSRRGGSGFSASYNCOEPPSSHIQ  
PGLCGLGNLGNTECFMNSALQCLSNAPLTDY  
FLKDEYEAENRDNPLGMKGEIAEAYAELIKQ  
MWSGRDAHVAPRMFKTQVGRFAPQFSGYQ  
QQDSQELAFLLDGLHEDLNRVKKPYLELKD  
ANGRPDAVVAKEAWENHRLRNDSDVIVD  
TFHGLFKSTLVCECAKVSVTFDFPCYLTL  
PLPLKDRVMEVFLVPADPHCRPTQYRVTV  
PLMGAVSDLCEALSRLSGIAAENMVVAD  
VYNHRFHKIFQMDGLNHIMPRDIFVYEV  
CSTSVDGSECVTLPVYFRERKSRPSSST  
SSASALYGPQLLSVPHKHLTLESLYQAVCD  
RISRYVKQPLPDEFSGSPLEPGACNGSRNS  
CEGEDEEMEHQEEGKEQLSETEGSGEDE  
PGNDPSETTQKKIKGQPCPKRLFTFSLVN  
SYGTADINSLAADGKLLKLNRSSTLAMDW  
DSETRRLYDEQSEAYEKHVSMLQPQKKK  
TTVALRDCIELFTTMETLGEHDPWYCPNCK  
KHQOATKFDLWLSLPKILVVHLKRFSYNRY  
WRDKLDTVVEFPPIRGLNMFVNCNLSAR  
PYVVDLIAVSNHYGAMGVGHYTAYAKN  
KLNKQWYFDDSNVSLASEDQIVTKAAYV  
LFYQRRDDEFYKTPSLSSSGSSDGGTRP  
SSSQQGFGDDEACSMDTN

Tag (**bold text**): N-terminal His  
Protease cleavage site: PreScission™ (LEVLFQ▼GP)  
USP4 (regular text): Start **bold italics** (amino acid residues 1-963)  
Accession number: NP\_003354



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