

USP5 [6His-tagged]

Deubiquitylating Enzyme

Alternate Names: Isopeptidase T, ISOT

Cat. No. 64-0002-050

Lot. No. 1735

Quantity: 50 µg

Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 2

Background

The Deubiquitylating enzymes (DUBs) 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 protein (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 processing protease 5 (USP5) is a member of the cysteine protease enzyme family and cloning of the human gene was first described by Wilkinson *et al.* (1995). USP5 protein contains an N-terminal zinc finger ubiquitin-binding domain (ZNF-UBP), a ubiquitin-specific processing protease (UBP) domain containing the active-site cys and his boxes, and two ubiquitin-associated domains (UBA1 and UBA2) (Reyes-Turcu *et al.*, 2006). Crystal structures of the ZNF-UBP domain have revealed a deep binding pocket where the C-terminal diglycine motif of ubiquitin is inserted explaining the specificity of USP for an unmodified C-terminus on the proximal subunit of polyubiquitin (Reyes-Turcu *et al.*, 2006). Recently USP5 has been identified in association with the 26S proteasome alongside other proteins

Continued on page 2

Physical Characteristics

Species: human

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: ~98 kDa

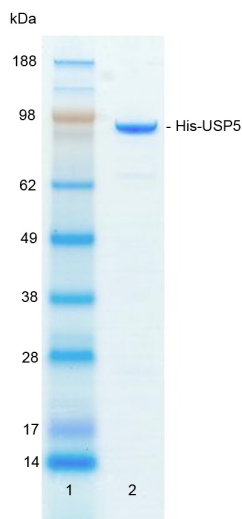
Purity: >90% by InstantBlue™ SDS-PAGE

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 His-USP5



Protein Identification:

Confirmed by mass spectrometry.

Deubiquitylating Enzyme Assay:

The activity of His-USP5 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-USP5 was compared confirming the deubiquitylating activity of His-USP5.



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

USP5 [6His-tagged] Deubiquitylating Enzyme

Alternate Names: Isopeptidase T, ISOT

Cat. No. **64-0002-050**
Lot. No. **1735**

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

known in complex as the UBL interactome (Besche *et al.*, 2009).

References:

Besche H, Haas W, Gygi S, Goldberg A (2009) Isolation of mammalian 26S proteasomes and p97/VCP complexes using the ubiquitin-like domain from HHR23B reveals novel proteasome-associated proteins. *Biochemistry* **48**, 2538-49.

Komander D, Clague MJ, Urbe S (2009) Breaking the chains: structure and function of the deubiquitinases. *Nat Rev Mol Cell Biol* **10**, 550-63.

Reyes-Turcu FE, Horton JR, Mullally JE, Heroux A, Cheng X, Wilkinson KD (2006) The ubiquitin binding domain ZnF UBP recognizes the C-terminal diglycine motif of unanchored ubiquitin. *Cell* **124**, 1197-208.

Wilkinson KD (1995) Roles of ubiquitylation in proteolysis and cellular regulation. *Annu Rev Nutr* **15**, 161-89.

Physical Characteristics

Continued from page 1

Protein Sequence:

MGS SHHHHS SGLEVLFQGPSPNS
PLGSM AELSEEALLSVLPTIRVPK
AGDRVHKDECAFSFDTPSEGGLYIC
MNTFLGFGKQYVERHFNKTGQRVYLHL
RRTRRPKEEDPATGTGDPPrKkPTR
LAIGVEGGFDLSEKFELEDVKIVILP
DYLEIARDGLGGLPDIVRDRVTSAVEALL
SADSASRKQEVQAWDGEVRQVSKHAFS
LKQLDNPARI PPGWKCSKCDMRENL
WLNLT DGSILCGRRYFDGSGGNHAVE
HYRETGYPLAVKLGITPDGADVSYDED
DMVLDPSLAEHL SHFGIDMLKMOKTKD
TMTELEIDMNQRIGEWELIQESGVPLK
PLFGPGYTGIRNLGNSCYLNSVVQVLF
SIPDFQRKYVDKLEKIFQNAPTDPTQD
FSTQVAKLGHGLLSGEYSKVPESGDG
ERVPEQKEVQDGIAPRMFKALIGKGHPEF
STNRQQAQEFFLHLINMVERNCRSSNP
NEVFRFLVEEKIKCLATEKVKYTORVDY
IMQLPVPMDAALNKEELLEEEK
KRQAE E EKMA LPELVRAQVPFSS
CLEAYGAPEQVDDFWSTALQAKS
VAVKTRRFASFDPDYLVIQIKKFT
FGLDWVPK KLDVSIEMPEELD IS
QLRGTGLQPGE EELPDIAPPLVTPDEP
KGS LGFYGN EDEDSFCSPHFSSPTSP
MLDESVI IQLVEMGFPM DACRKAVYYT
GNSGAEAMNWVM SHMDDPDFANPLIL
PGSSGPGSTSAADPPPEDCVTTIVSMGF
SRDQALKALRATNNSLERAVDWIFSHID
DLDAEAAMDISEGRSAADSI SESVPVGP
KVRDGP GKYQLFAFISHMGTSTMCGHYV
CHIKKEGRWVIYNDQKVCASEKPPKDLGY
IYFYQRVAS

Tag (**bold text**): N-terminal His
Protease cleavage site: PreScission™ (**LEVL**FQ▼**GP**)
USP5 (regular text): Start **bold italics** (amino acid residues 1-858)
Accession number: P45974



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