

USP6 CD(529-1406) [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: TRE2, TRE17, HRP1, Ubiquitin carboxyl terminal hydrolase 6, Ubiquitin specific processing protease 6, Ubiquitin thiolesterase 6

Cat. No. 64-0045-050

Lot. No. 30155

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 specific protease 6 (USP6) is a member of the cysteine protease enzyme family and cloning of the gene was first described by Nakamura *et al.* (1992). USP6 was the first DUB to be identified as an oncogene (Oliveira and Chou, 2012). Multiple DUBs, including A20, CYLD, Cezanne, USP21 and USP15, have been shown to function as negative regulators of NFκB; however, only one DUB, USP6, induces NFκB activation (Pringle *et al.*, 2012). USP6 is translocated and overexpressed in aneurysmal bone cyst (ABC), a paediatric tumour characterized by extensive bone degradation and inflammatory recruitment (Pringle *et al.*, 2012). Recent work has shown that USP6 is part of a signalling pathway that contributes to ABC pathogenesis, raising the possibility that development of USP-specific

Continued on page 2

Physical Characteristics

Species: human

Source: insect (Sf21)

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: ~125.3 kDa

Purity: >88% by InstantBlue™ SDS-PAGE

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

Protein Sequence: Please see page 2

Quality Assurance

Purity:

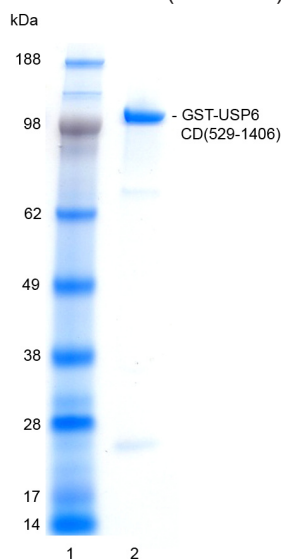
4-12% gradient SDS-PAGE

InstantBlue™ staining

Lane 1: MW markers

Lane 2: 1 µg GST-USP6

CD(529-1406)



Protein Identification:

Confirmed by mass spectrometry.

Deubiquitylase Enzyme Assay:

The activity of GST-USP6 CD(529-1406) 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-USP6 CD(529-1406) was compared confirming the deubiquitylating activity of GST-USP6 CD(529-1406).



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 2013. 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

USP6 CD(529-1406) [GST-tagged]

Deconjugating enzyme: Deubiquitylase

Alternate Names: TRE2, TRE17, HRP1, Ubiquitin carboxyl terminal hydrolase 6, Ubiquitin specific processing protease 6, Ubiquitin thiolesterase 6

Cat. No. **64-0045-050**
Lot. No. **30155**

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

inhibitors or NF-κB antagonists might be effective novel strategies for the treatment of these tumours (Ye *et al.*, 2010).

References:

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.

Nakamura T, Hillova J, Mariage-Samson R, Onno M, Huebner K, Cannizzaro LA, *et al.* (1992) A novel transcriptional unit of the tre oncogene widely expressed in human cancer cells. *Oncogene* **7**, 733-741.

Oliveira AM and Chou MM (2012) The TRE17/USP6 oncogene: a riddle wrapped in a mystery inside an enigma. *Front Biosci (Schol Ed)* **4**, 321-334.

Pringle LM, Young R, Quick L, Riquelme DN, Oliveira AM, May MJ, *et al.* (2012) Atypical mechanism of NF-κB activation by TRE17/ubiquitin-specific protease 6 (USP6) oncogene and its requirement in tumorigenesis. *Oncogene* **31**, 3525-3535.

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

Ye Y, Pringle LM, Lau AW, Riquelme DN, Wang H, Jiang T, *et al.* (2010) TRE17/USP6 oncogene translocated in aneurysmal bone cyst induces matrix metalloproteinase production via activation of NF-κB. *Oncogene* **29**, 3619-3629.

Physical Characteristics

Continued from page 1

MSPILGYWKIKGLVQPTRLLEYLEEKYEEH
LYERDEGDKWRNKKFELGLEFPNLPYYIDGD
VKLTQSMAIRYIADKHNMLGGCPKERAEISM
LEGAULD IRYGVSRIAYSKDFETLKVDFL
SKLPEMLKMFEDRLCHKTYLNGDHVTHPD
FMLYDALDVVLYMDPMCLDAFPKLVCFK
KRIEAIPOIDKYLKSSKYIAWPLQGWQATF
GGGDHPPKSDLEVLFOGPLGSKGATGLSNL
GNTCFMNSSIQCVSNTQPLTQYFISGRH
LYELNRTNPIGMKGHMAKCYGDLVQELWS
GTQKSVAPLKLRRTI AKYAPKFDGFGQQD
SQELLAFLLDGLHEDLNRVHEKPYVELKDS
DGRPDWEVAEAEAWDNHLRRNRSIIVDLFH
GQLRSQVKCKTCGHSVRFDPFNFLSLPLP
MDSYMDLEITVIKLDGTTTPVRYGLRL
NMDEKYTGLKKQLRDL CGLNSEQIL
LAEVHDSNIKNFPQDNQKVQLSVSGFL
CAFEIPVPSSPISASSPTQIDFSSSPSTNGM
FTLTTNGDLPKPIFIPNGMPNTVVP CGTEKN
FTNGMVNGHMPSLPDSPTGYIIAVHRKM
MRTELYFLSPQENRPSLFGMPLIVPCT
VHTRKKDLYDAVWIQVSWLARPLPPQEASI
HAQDRDNCMGYQYPTLRRVVQKDGNSCAW
CPQYRFRGCKIDCGEDRAFIGNAYI
AVDWHPTALHLRYQTSQERVVDKHESVEQS
RAQAEPINLDSCLRAFTSEELGESE
MYYCSKCKTHCLATKKLDLWRLPPFLII
HLKRFQFVNDQWIKSQKIVRFLRESFDP
SAFLVPRDPALCQHKPLTPQGDLSKPRI
LAREVKKVDAQSSAGKEDMLLSKSPSSLSANI
SSSPKGPSSSRKSGTSCPSSKNSSPNS
SPRTLGRSKGRLRPLQIGSKNKPSSSK
KNLDASKENGAGQICELADALSRGHM
RGGSQPELVTPQDHEVALANGFLYE
HEACNGCGDGY SNGQLGNHSEEDSTD
DQREDTHIKPIYNLYAISCHSGILSGGHY
ITYAKPNCKWYCYNDSSCEELHPDEIDTD
SAYILFYEQQGIDYAQFLPKIDGKKMADTSS
DEDESDYEKYSMLQ

Tag (**bold text**): N-terminal GST
Protease cleavage site: PreScission™ (LEVLFO▼GP)
USP6 CD(529-1406) (regular text): Start **bold italics**
(amino acid residues 529-1406)
Accession number: NP_004496



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 2013. 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