

UBA7 [6His-tagged]

E1 Activating Enzyme

Alternate Name: UBE1L

Cat. No. 61-0007-010
Lot. No. 30080

Quantity: 10 µg
Storage: -70°C



FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS

CERTIFICATE OF ANALYSIS Page 1 of 2

Background

The enzymes of the ISGylation pathway play a pivotal role in the immune response. Three classes of enzymes are involved in the process of ISGylation an activating enzyme (E1), conjugating enzymes (E2s) and protein ligases (E3s). UBA7 is a member of the E1 activating enzyme family and cloning of the human gene was first described by Kok *et al.* (1993). The UBA7 gene has been mapped to chromosome 3p21 by high resolution fluorescence in situ hybridization (Carritt *et al.* 1992). UBA7 and UBE2L6 have been reported to function as the E1 and E2 enzymes respectively for ISG conjugation forming a thioester intermediate through Cys-131 of UBE2L6 (Takeuchi *et al.* 2005). UBA7 (Ube1L) knockout (KO) mice are deficient in ISGylation and are fertile with no obvious phenotype. (Kim *et al.* 2006). The expression of UBA7 was found to be reduced in many lung cancer cell lines and although originally thought to be a tumor suppressor gene candidate recent studies have demonstrated that UBA7 does not suppress the development of lung adenoma or thymic lymphoma in a K-rasLA2 cancer model (Kok *et al.* 1993).

References:

Carritt B, Kok K, van den Berg A, Osinga J, Pilz A, et al. (1992) A gene from human chromosome region 3p21 with reduced expression in small cell lung cancer. *Cancer Res* **52**, 1536-1541.

Kim KI, Yan M, Malakhova O, Luo JK, Shen MF, et al. (2006) Ube1L and protein ISGylation are not essential for alpha/beta interferon signaling. *Mol Cell Biol* **26**, 472-479.

Kok K, Hofstra R, Pilz A, van den Berg A, Terpstra P, et al. (1993) A gene in the chromosomal region 3p21 with greatly reduced expression in lung cancer is similar to the gene for ubiquitin-activating enzyme. *Proc Natl Acad Sci USA* **90**, 6071-6075.

Continued on page 2

Physical Characteristics

Species: human

Protein Sequences: Please see page 2

Source: Insect sf21

Quantity: 10 µg

Concentration: 0.5 mg/ml

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

Molecular Weight: 116.53 kDa

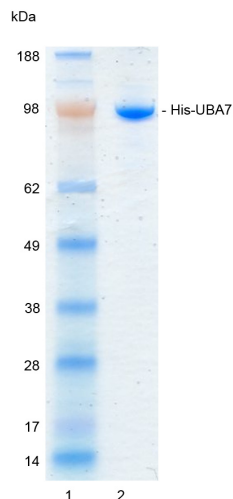
Purity: >98% by InstantBlue™ SDS-PAGE

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-UBA7



Protein Identification:

Confirmed by mass spectrometry.

E1 Thioester ISG15 Loading Assay:

The activity of His-UBA7 was validated by loading ISG15 onto the active cysteine of His-UBA7. Incubation of the His-UBA7 enzyme in the presence of ISG15 and ATP at 30°C was compared at two time points, T₀ and T₁₀ minutes. Sensitivity of the ISG15/His-UBA7 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 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

UBA7 [6His-tagged]

E1 Activating Enzyme

Alternate Name: UBE1L

Cat. No. 61-0007-010

Lot. No. 30080

Quantity: 10 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

Takeuchi T, Iwahara S, Saeki Y, Sasajima H, Yokosawa H (2005) Link between the ubiquitin conjugation system and the ISG15 conjugation system: ISG15 conjugation to the UbcH6 ubiquitin E2 enzyme. *J Biochem* 138, 711-719.

Physical Characteristics

Continued from page 1

Protein Sequence:

MSY YHHHHHDYDIP TTE NLY FOGAMGS
GIQRPTSTSSLVAAA **MDALD**ASKLL
DEELYSRQLYVLGSPAMQRIQGARVLVS
GLQGLGAEVAKNLVLMGVGSLTLHDPHPTCWS
DLAAQFLLSEQDLERSRAEASQELLAQLN
RAVQVVVHTGDITEDLLLDLDFQVVVLTAAKLE
EQLKVGTLCHKHGVCFLAADTRGLVGLFCD
FGEDFTVQDPTEAEP LTA A I Q H I S Q G
SPGILTLRKGANTHYFRDGDLVTFSGIEGM
VELNDCDPRSIHVREDGSLEIGDTTTFSTRYL
RGGAI TEVKRPKTVRHKSLDTALLQPHV
VAQSSQEVHHAHCLHQAF CALHKFQHLHGR
PPQPWDPVDAETVVGLARDLEPLKRTEEE
PLEEPLDEALVRTVALSSAGVLSPMVAML
GAVAAQEV LKAISRKFMP LDQWLYFDALD
CLPEDGELLSPEDCALRGSRYDGGIAVF
GAGFQEKLRROHYLLV GAGAIGCELLKV
FALVGLGAGNSGGLTVVDMDH IERSNLSRQ
FLFRSQDVGRPKAEVAAAAARGLNPDQVI
PLTYPLDPTTEHIYGDNFFSRVDGVAALDS
FQARRYVAARCTHYLKPLLEAGTSGTWGSAT
VFM PHVTEAYRAPASAAASEDAPYPVCT
VRYFPSTA EHTLQWARHEFEELFRLSAET
INHHQQAHTSLADMDEPQTLTLLKPVLGVL
RVRPQNWQDCVAWALGHWKL CFHYGIKQLL
RHFPNKVLEDGTPFWSGPKQCPQPLEFDT
NQDTHLLYVLAANLYAQM HGLPGSQD
WTALRELLKLLPQDPQMAPIFASNLELA
SASAEFGPEQQKELNKALEVWSVGPPLKPLM
FEKDDSNFHVDFVVAASLRCONYGI
PVNRAQSKRIVGQIIPAIATTTAAVAGLL
GLELYKVVSGPRPRSAFRHSYLHLAENY
LIRYMPFAPAIQTFHHLKWT SWDR LKV
PAGQPRTL ESSL AHLQE QHGLRVRILLHG
SALLYAAGWSPEKQAQHLPLRVTEL VQQLT
GQAPAPGQRV LVLELSCEGDEDDTAFPPLHYEL

Tag (bold text): N-terminal His

Protease cleavage site: TEV (**ENLYF**▼**QG**)

UBA7 (regular text): Start **bold italics** (amino acid residues 1-1012)

Accession number: NP_003326.2



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