

ATG7 [6His-tagged]

E1 Activating Enzyme

Alternate Name: APG7, GSA7

Cat. No. 61-0008-050

Lot. No. 30081

Quantity: 50 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 autophagy pathway play a pivotal role in the degradation of cytoplasmic constituents and organelles. Structures known as autophagosomes sequester portions of the cytoplasm which are degraded by the lysosome and recycled back into the cell (Kuma *et al.* 2004). Three classes of enzymes are involved in autophagy; E1-like activating enzymes, E2-like conjugating enzymes and E3-like ligases. Ubiquitin-like proteins (ubl) Autophagy 12 (ATG12) and ATG8 are activated by ATG7 the E1-like activating enzyme. ATG12 and ATG8 are transferred to two E2-like conjugating enzymes ATG10 and ATG3 respectively. Cloning of the human ATG7 gene was first described by Yuan *et al.* (1999) and ATG7 shares 38% sequence identity to its yeast homologue Apg7. ATG7 can also activate ATG8 mammalian homologues, GABARAP and GATE-16. It forms a homodimer via the C-terminal region that is important for enzyme substrate interaction and E1-E2 complex formation (Komatsu *et al.* 2005). ATG7 Δ FAP a mutant form of the ATG7 E1-like activating enzyme which lacks the (Phe-Asp-Pro) FAP motif has been shown to be unable to form an E2 substrate intermediate with ATG3 and the ubl Microtubule-Associated Protein 1, Light Chain 3 (MAPLC3) (Tanida *et al.* 2012). ATG7 has been shown to be involved in a novel pathway in which the inhibition of caspase-8 results in autophagic death induced by receptor-interacting protein (RIP), Jun amino-terminal kinase, and beclin-1 (Yu *et al.* 2004).

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: 81.34 kDa

Purity: >95% by InstantBlue™ SDS-PAGE

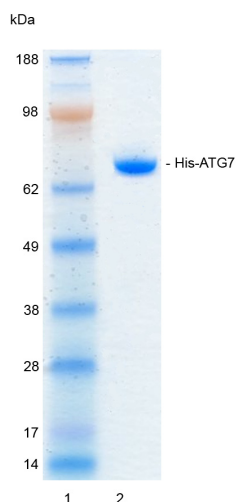
Stability/Storage: 12 months at -70°C; aliquot as required

Protein Sequences: Please see page 2

Quality Assurance

Purity:

4-12% gradient SDS-PAGE
InstantBlue™ staining
Lane 1: MW markers
Lane 2: 1 µg His-ATG7



Protein Identification:

Confirmed by mass spectrometry.

E1 Thioester GABARAP Loading Assay:

The activity of His-ATG7 was validated by loading GABARAP onto the active cysteine of His-ATG7. Incubation of the His-ATG7 enzyme in the presence of GABARAP and ATP at 30°C was compared at two time points, T₀ and T₁₀ minutes. Sensitivity of the GABARAP/His-ATG7 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

ATG7 [6His-tagged]

E1 Activating Enzyme

Alternate Name: APG7, GSA7

Cat. No. 61-0008-050

Lot. No. 30081

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

References:

Komatsu M, Waguri S, Ueno T, Iwata J, Murata S, *et al.* (2005) Impairment of starvation-induced and constitutive autophagy in Atg7-deficient mice. *J Cell Biol* **169**, 425-434.

Kuma A, Hatano M, Matsui M, Yamamoto A, Nakaya H, *et al.* (2004) The role of autophagy during the early neonatal starvation period. *Nature* **432**, 1032-1036.

Tanida I, Yamasaki M, Komatsu M, Ueno T (2012) The FAP motif within human ATG7, an autophagy-related E1-like enzyme, is essential for the E2-substrate reaction of LC3 lipidation. *Autophagy* **8**, 88-97.

Yu L, Alva A, Su H, Dutt P, Freundt E, *et al.* (2004) Regulation of an ATG7-beclin 1 program of autophagic cell death by caspase-8. *Science* **304**, 1500-1502.

Yuan W, Stromhaug PE, Dunn WA, Jr. (1999) Glucose-induced autophagy of peroxisomes in *Pichia pastoris* requires a unique E1-like protein. *Mol Biol Cell* **10**, 1353-1366.

Physical Characteristics

Continued from page 1

Protein Sequence:

MSY YHHHHHHDYDIPTT**ENLYF**QGAMGS
MAAATGDPGLSKLQFAPFSSALDVGFWHEL
TQKKLNEYRLDEAPKDIKGYYYNGDSAGL
PARLTLEFSAFDMSAPTPARCCPAIGTLYNT
NTLESFKTADKLLLEQAANEIWESI KSG
TALNPVLLNKFLLLTFADLKKYHFYWF
CYPALCLPESLPLIQGPVGLDQRFSLKQIEA
LECA YDNLCQTEGV T ALPYFLIKYDENM
VLVSL LKHYSDF FQGQRTKITIGVYDPCN
LAQYPGWPLRNFLVLAHRWSSSFQSVEVVC
FRDRTMQGARDVAHSIIFEVKLPEMAFSPDCP
KAVGWEKNQKGGMGRMVNLSECMDPKRLAES
SVDLNLKLMCWRLVPTLDL DKVVSVKCLLL
GAGTLGCNVAR TLMGWGVRHITFVDNAKISYS
NPVRQPLYEFEDCLGGGKPKALAAADRLQKIF
PGVNARGFNMSIPMPGHPVNFSSVTLEQARRD
VEQLEQLIESHDVVFL LMDTRESRWLP AVI
AASKRKLVINAAALGFDTFVVMRHGLKPKQ
GAGDLCPNHPVASADLLGSSLFANIPGYKLG
CYFCNDVVAPGDSTRDR TLDQQCTVSRPGLA
VIAGALAVELMVSVLQHPEGGYAIASSSDDRM
NEPPTSLGLVPHQIRGFLSRFDNVL PVSLAFD
KCTACSSKVL DQYEREGFNFLAKVFNSSHS
FLEDLTGLTLLHQETQAAEIWDMSDDETI

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

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

ATG7 (regular text): Start **bold italics** (amino acid residues 1-706)

Accession number: NP_006386



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