

# TRIAD1 [6His-tagged]

## E3 Ligase

Alternate Names: Triad domain-containing protein 1; *Drosophila Ariadne* homolog 2; ARIH2

Cat. No. 63-0029-025

Lot. No. 30026

Quantity: 25 µg

Storage: -70°C

FOR RESEARCH USE ONLY

NOT FOR USE IN HUMANS



CERTIFICATE OF ANALYSIS Page 1 of 1

## Background

The enzymes of the ubiquitylation pathway play a pivotal role in a number of cellular processes including the regulated and targeted proteasome-dependent degradation of substrate proteins. Three classes of enzymes are involved in the process of ubiquitylation; activating enzymes (E1s), conjugating enzymes (E2s) and protein ligases (E3s). Triad Domain-Containing Protein 1 (TRIAD1) is a member of the E3 protein ligase family and cloning of the human gene was first described by van der Reijden *et al.* (1999). TRIAD1 contains a TRIAD motif containing two RING domains which flank a conserved cysteine-rich (C6HC) domain designated DRIL (double RING finger-linked domain) (Martelijn *et al.*, 2005). TRIAD is thought to be involved in protein translation, interacting with Ubch7 (UBE2L3) to polyubiquitylate eIF4E2 targeting it for proteasomal degradation (Tan *et al.*, 2003). More recently these proteins have been referred to as Ring in-between Ring E3 ligases (RBRs) that function like RING-HECT hybrids regulating processes such as translation and immune signaling (Wenzel *et al.*, 2011).

## References:

Martelijn JA, van Ernst L, Erpelinck-Verschueren CA, Nikoloski G, Menke A, de Witte T, Lowenberg B, Jansen JH, van der Reijden BA (2005) The E3 ubiquitin-protein ligase Triad1 inhibits clonogenic growth of primary myeloid progenitor cells. *Blood* 106, 4114-23.

Tan NG, Ardley HC, Scott GB, Rose SA, Markham AF, Robinson PA (2003) Human homologue of ariadne promotes the ubiquitylation of translation initiation factor 4E homologous protein, 4EHP. *FEBS Lett* 554, 501-4.

van der Reijden BA, Erpelinck-Verschueren CA, Lowenberg B, Jansen JH (1999) TRIADs: a new class of proteins with a novel cysteine-rich signature. *Protein Sci* 8, 1557-61.

Wenzel DM, Lissounov A, Brzovic PS, Klevit RE (2011) UBCH7 reactivity profile reveals parkin and HHAR1 to be RING/HECT hybrids. *Nature* 474, 105-8.

## Physical Characteristics

Species: human

Source: Sf21 insect cell-baculovirus expression

Quantity: 25 µg

Concentration: 0.5 mg/ml

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

Molecular Weight: ~61.2 kDa

Purity: >98% by InstantBlue™ SDS-PAGE

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

## Protein Sequence:

**MSYYHHHHHDYDIP****TTENLYFQ**GAMGSM  
VDMNSQGS DSDNEEDYDPNCEEEEEEEED  
DPGDIEDYVGVASDVEQQGADAFDPEEYQFTCL  
TYKSEGALNEHMTSLASVLKVSHSVAKLILVNFH  
WQVSEILDYKSNQAQLLVEARVQPNPSKHVPTSH  
PHCAVCMQFVRKENLLSLACQHQFRCSCWEQHC  
LVKDGVGVGVSQMAQDCPLRTPEDFVFPPLPNEEL  
REKYRRLFRDYVESHYQLQLCPGADCPMVI  
RVQE PRARRVQCNRCNEVFCFKCRQMYHAPTDCATIRK  
WLTCKADDSETANYISAHTKDCPKNCICIEKNG  
GCNHMQCSKCKHDFCWMCLGDWKTGSEYYECS  
RYKENPDIVNQSQAQAREALKKYLFFERWENHNK  
SLQLEAQT YQRIHEKIQERVMNNLGTWIDWQYLO  
NAAKLLAKCRYTQTYTPYAYYMEGPRKKLFY  
QQAQLEAEIENLSWKVERADSYDRGDLNQMH  
IEQRRTLLKDFHDT

Tag (bold text): N-terminal His

Protease cleavage site: TEV (ENLYF↓QG)

TRIAD1 (regular text): Start **bold italics** (amino acid residues 1-493)

Accession number: NP\_006312

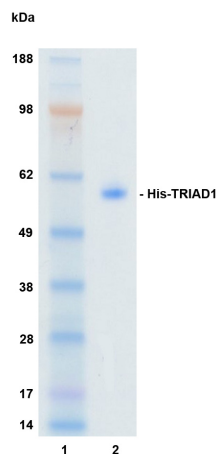
## Quality Assurance

### Protein Identification:

Confirmed by mass spectrometry.

### Purity:

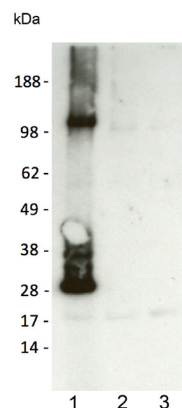
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 1 µg His-TRIAD1



### E3 ligase assay:

The ubiquitin conjugating activity of His-TRIAD1 was validated through its ability to catalyse the generation of polyubiquitin chains in the presence of the E1 activating enzyme His-UBE1, the E2 conjugating enzyme His-UBE2D1 (Ubch5a) (several E2s were tested, data generated with this E2 is provided by way of example) and ubiquitin. Incubation of His-

TRIAD1 for 120 minutes at 37°C in the presence of ubiquitin, His-UBE1, His-UBE2D1 and ATP (Lane 1) was compared alongside two control reactions with either ATP (Lane 2) or His-TRIAD1 (Lane 3) excluded from the reaction. Ubiquitin conjugates were identified by Western blotting using an anti-ubiquitin conjugate antibody and these were observed only in the presence of both ATP and His-TRIAD1.



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: v1.0.1