

# Ataxin-3L [6His-tagged]

## Deconjugating enzyme: Deubiquitylase

Alternate Name: Machado-Joseph disease protein 1-like

Cat. No. **64-0034-050**  
Lot. No. **30077**

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. Ataxin-3L is a cysteine protease and a member of the Machado-Joseph Domain (MJD) enzyme family. Cloning of the human gene was first described by Gerhard *et al.* (2004). Machado-Joseph disease (MJD), the most common form of spinocerebellar ataxia worldwide, is a progressive and ultimately fatal neurodegenerative disorder caused by polyQ expansion in ataxin-3, a conserved and ubiquitous protein known to bind polyubiquitin chains and to function as a deubiquitylating enzyme. Ataxin-3 has been linked to protein homeostasis

Continued on page 2

### Physical Characteristics

**Species:** human

**Source:** *E. coli*

**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:** ~43 kDa

**Purity:** >92% by InstantBlue™ SDS-PAGE

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

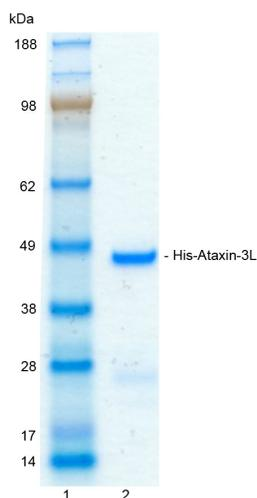
### Protein Sequence:

**MGSSHHHHHSSGLEVLFGPGSMDFIF**  
HEKQEGFLCAQHCNNLLQGEYFSPVELA  
SIAHQLDDEERMMAEGGVTSSEYLAFLQPS  
ENDDTGFFSIQVISNALKFWGLEIIHFNNPEY  
QKLGIDPINERSFICNYKQHWFTIRKFGKHWFN  
LNSLLAGPELISDTCLANFLARLQQQAYSVFV  
VKGDLPDCEADQLLQIISVEEMDTPKLNK  
KLVKQKEHRVYKTVLEKVSEESDESSTSDQDEED  
FQRALELSRQETNREDEHLRSTIELSMQGS  
SNTSQDLPKTSCVTPASEQPKKIKEDY  
FEKHQQEQKQQQQSDLPGHSSYLHERPTTSS  
RAIESDLSDDISEGTVQAAVDITLEIMRKNLKI  
GK

Tag (**bold text**): N-terminal His  
Protease cleavage site: PreScission™ (LEVLFG▼GP)  
Ataxin-3L (regular text): Start **bold italics** (amino acid residues 1-355)  
Accession number: NP\_001129467

### Quality Assurance

**Purity:**  
4-12% gradient SDS-PAGE  
InstantBlue™ staining  
Lane 1: MW markers  
Lane 2: 1 µg His-Ataxin-3L



### Protein Identification:

Confirmed by mass spectrometry.

### Deubiquitylase Enzyme Assay:

The activity of His-Ataxin-3L 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-Ataxin-3L was compared confirming the deubiquitylating activity of His-Ataxin-3L.



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

# Ataxin-3L [6His-tagged]

## Deconjugating enzyme: Deubiquitylase

Alternate Name: Machado-Joseph disease protein 1-like

Cat. No. 64-0034-050

Lot. No. 30077

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

maintenance, transcription, cytoskeleton regulation and myogenesis (Matos *et al.*, 2011). Ataxin-3L shares 85% sequence identity with ataxin-3 although it has recently been shown that the Josephin domain of ataxin-3L demonstrates substantially higher deubiquitylating activity than the ataxin-3 Josephin domain (Weeks *et al.*, 2011).

### References:

Gerhard DS, Wagner L, Feingold EA, Shenmen CM, Grouse LH, Schuler G, *et al.* (2004) The status, quality, and expansion of the NIH full-length cDNA project: the Mammalian Gene Collection (MGC). *Genome Res* **14**, 2121-2127.

Komander D, Clague MJ, Urbe S (2009) Breaking the chains: structure and function of the deubiquitinases. *Nat Rev Mol Cell Biol* **10**, 550-563.

Matos CA, de Macedo-Ribeiro S, Carvalho AL (2011) Polyglutamine diseases: the special case of ataxin-3 and Machado-Joseph disease. *Prog Neurobiol* **95**, 26-48.

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

Weeks SD, Grasty KC, Hernandez-Cuevas L, Loll PJ (2011) Crystal structure of a Josephin-ubiquitin complex: evolutionary restraints on ataxin-3 deubiquitinating activity. *J Biol Chem* **286**, 4555-4565.



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