

Ataxin3 (Ataxin 3)

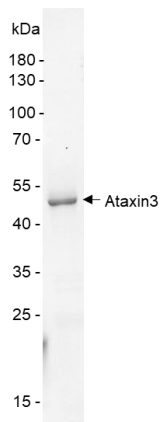
Cat. # DB401

Background	Ataxin-3 (MJD1, SCA3) is a josephin domain-containing protein that functions as a deubiquitinase <i>in vitro</i> . It mediates the cleavage of long polyubiquitin chains while it has weak or no activity against chains of 4 or less ubiquitins. Ataxin-3 is the disease protein in spinocerebellar ataxia type 3 and plays a role in protein homeostasis maintenance, transcription, cytoskeleton regulation, myogenesis and degradation of misfolded chaperone substrates.
Alternate names	AT3, Ataxin-3, ATX3, JOS, Machado-Joseph disease protein 1, MJD, MJD1, SCA3, Spinocerebellar ataxia type 3 protein

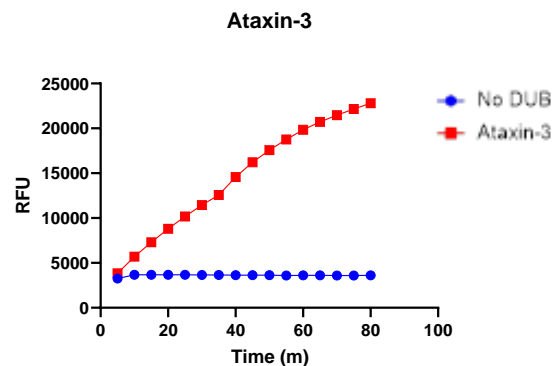
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	41 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Activity	This enzyme cleaves K48-linked ubiquitin chains <i>in vitro</i> .
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified Ataxin3. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of Ataxin-3. Ataxin-3 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

1. Carvalho, A., et al., *Adv Exp Med Biol.*, 2018. 1049:275-288.
2. Sowa, A., et al., *Front Immunol.*, 2022. 13:870966.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved