

ATXN3-Like (Ataxin 3-like)

Cat. # DB402

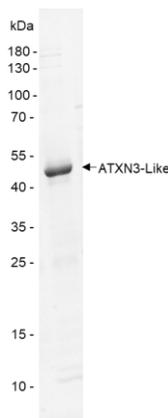
Background Ataxin 3-like is highly homologous to ataxin 3 but currently no homolog has been found in an organism other than humans. Ataxin 3-like contains a josephin domain and can cleave Ub-AMC *in vitro*.

Alternate names Machado-Joseph disease protein 1-like, MJDL, ATXN3L

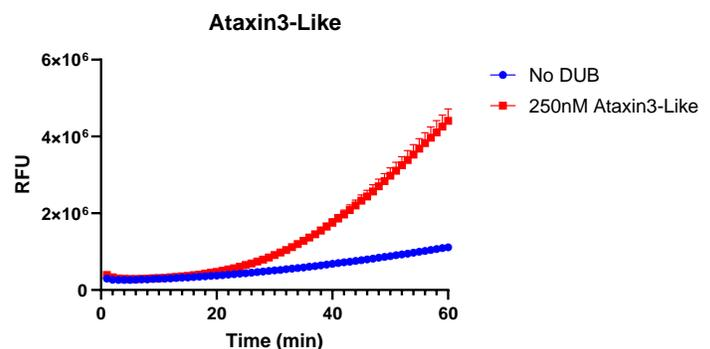
Product Information

Purity	≥ 85% by SDS-PAGE
Molecular Weight	41 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Activity	This enzyme is active in the Ub-CHOP assay.
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified ATXN3-Like. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of ATXN3-Like. Ataxin3-like Protein was tested in a Ub-CHOP assay showing robust DUB activity.

References

1. Ge, F., et al., *Oncotarget*, 2015. 6(25):21369-78.
2. Lee, DH., et al., *Int J Mol Sci.*, 2020. 21(5):1755.

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