

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*¹

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

Product Information

Molecular Weight:	41kDa
Quantity:	25µg
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

References

- 1) N. Tzvetkov and P. Breuer. *Josephin domain-containing proteins from a variety of species are active de-ubiquitination enzymes*. Biol. Chem. 2007. 388: p973-978.

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