

Ataxin3 (Ataxin 3)

Cat. # DB401

Background: Ataxin 3 is the disease protein in spinocerebellar ataxia type 3. Ataxin 3 contains a josephin domain and can cleave poly-ubiquitin chains *in vitro*¹. The proteolytic activity of Ataxin 3 regulates the amount of both total cellular ubiquitylation and ataxin 3 protein^{2,3}.

Alternate names: AT3, Ataxin-3, ATX3, JOS, Machado-Joseph disease protein 1, MJD, MJD1, SCA3, Spinocerebellar ataxia type 3 protein

Product Information

Molecular Weight:	41kDa
Quantity:	25 µg
Physical State:	Liquid
Source:	Human Recombinant
Tag:	His6
Activity:	This enzyme cleaves K48-linked ubiquitin chains in vitro.
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

- 1) Burnett, B. et al. *The polyglutamine neurodegenerative protein ataxin-3 binds polyubiquitylated proteins and has ubiquitin protease activity.* Hum Mol Genet. 2003. **12**(23): p3195-205.
- 2) Todi, SV et al. *Cellular turnover of the polyglutamine disease protein ataxin-3 is regulated by its catalytic activity.* J Biol Chem. 2007. **282**(40): p29348-58.
- 3) Schmitt, I et al. *Inactivation of the mouse Atxn3 (ataxin-3) gene increases protein ubiquitination.* Biochem Biophys Res Commun. 2007. **362**(3): p734-9.

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