

JosD2 (Josephin domain containing 2)

Cat. # DB404

Background: Josephin Domain containing protein 2 (JosD2) is a member of the Machado-Josephin Domain family of DUBs.¹ When transfected into HEK293, COS-7, or SH-SY5Y cells, JosD2 is diffusely localized in the cytoplasm, in contrast to JosD1, which localizes specifically to the plasma membrane.² JosD2 has deubiquitylating activity in vitro.¹

Alternate names: FLJ29018, Josephin-2, SBB154

Product Information

Molecular Weight:	22kDa
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) Yao, T. et al. *Proteasome recruitment and activation of the Uch37 deubiquitinating enzyme by Adrm1*. Nature Cell Biology, 2006. 8: p.994-1002.
- 2) Seki, T. et al. *JosD1, a Membrane-targeted Deubiquitinating Enzyme, Is Activated by Ubiquitination and Regulates Membrane Dynamics, Cell Motility, and Endocytosis*. Journal of Biological Chemistry, 2013. 288: p.17145-17155.

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