

JosD2 Antibody (Josephin domain containing 2)

Cat. # AB404

Background

Josephin Domain containing protein 2 has deubiquitylating activity *in vitro*¹

Alternate Names: FLJ29018, Josephin-2, SBB154

Molecular Weight: 22kDa

Specificity/Applications

Species Cross Reactivity: Human

Source: Chicken

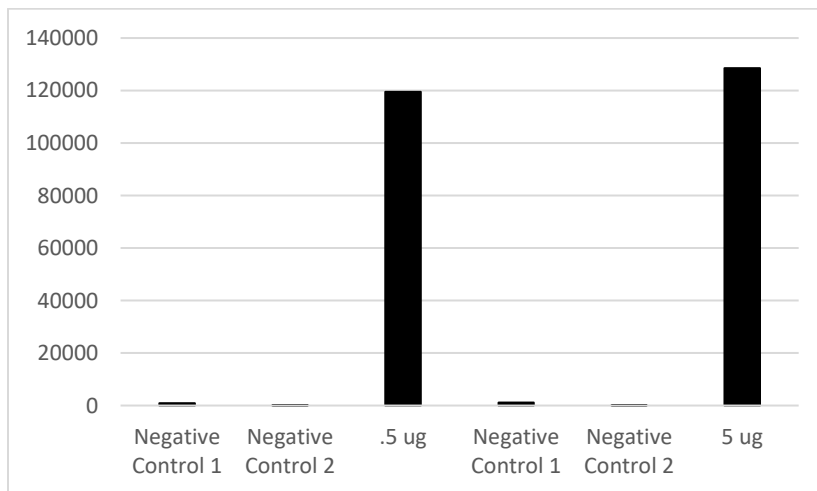
Applications: WB

Recommended Antibody Dilutions:

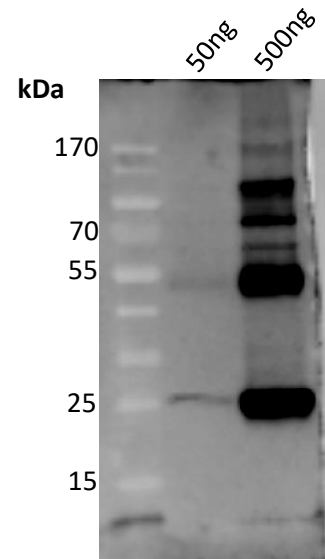
Western Blotting: Robust detection of 500ng of recombinant protein was possible when antibody was used at a final concentration of 8.57 µg/mL

Storage: Supplied in phosphate buffered saline. Store at -20°C or below. Avoid repeated freeze/thaw cycles.

Detection of JosD2 by AB404 using ELISA and Western Blot



0.5 or 5 ug of JosD2 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. JosD2 was detected by 1.71 µg/mL of AB404 using traditional ELISA detection reagent. 2^o Antibody: α-Chicken HRP (1:5000). Negative Control 1: No AB404 Negative Control 2: no detection reagent.



Indicated amounts of JosD2 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 8.57 µg/mL of AB404 2^o Antibody: α-Chicken HRP (1:5000).

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved

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.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved