

JosD1 Antibody (Josephin domain containing 1)

Cat. # AB403

Background

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

Alternate Names: dJ508I15.2, Josephin-1, KIAA0063

Molecular Weight: 22kDa

Specificity/Applications

Species Cross Reactivity: Human

Source: Chicken

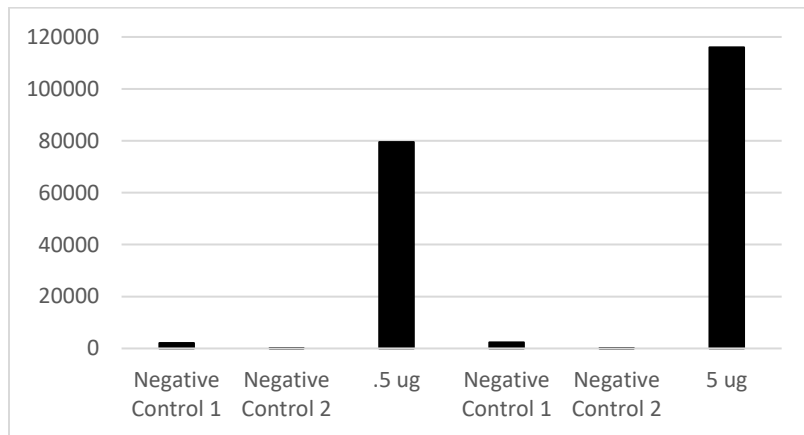
Applications: WB

Recommended Antibody Dilutions:

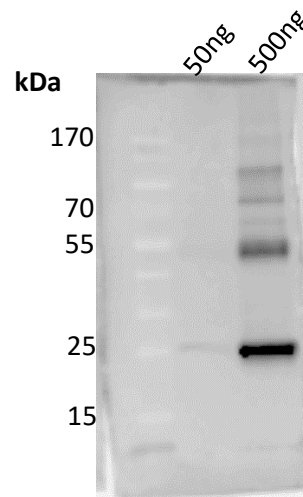
Western Blotting: Robust detection of 50ng of recombinant protein was possible when antibody was used at a final concentration of 12.57µg/mL

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

Detection of JosD1 by AB403 using ELISA and Western Blot



0.5 or 5 ug of JosD1 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. JosD1 was detected by 2.51 µg/mL of AB403 using traditional ELISA detection reagent. 2^o Antibody: α-Chicken HRP (1:5000).
 Negative Control 1: No AB403 Negative Control 2: no detection reagent.



Indicated amounts of JosD1 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 12.57 µg/mL of AB403. 2^o Antibody: α-Chicken HRP (1:5000).

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 © 2007 LifeSensors, Inc. All Rights Reserved