

USP4 Antibody (Ubiquitin-specific-processing protease 4)

Cat. # AB511

Background

USP4 (UNP) has been shown cleave the ubiquitin precursor UBICEP52 and the isopeptide (ϵ -peptide linked) ubiquitin conjugate ubiquitin-(ϵ N)-lysine *in vitro*¹. USP4 interacts with the tumor suppressors Retinoblastoma protein (Rb), p107 and p130 *in vitro* and in cells and ectopic expression of USP4 is oncogenic^{2,3}.

Alternate Names: Deubiquitinating enzyme 4, Ubiquitin carboxyl-terminal hydrolase 4, Ubiquitin-specific-processing protease 4, Ubiquitin thioesterase 4, Ubiquitous nuclear protein homolog, UNP, Unph, UNPH

Molecular Weight: 120kDa

Specificity/Applications

Species Cross Reactivity: Human

Source: Chicken, IgY

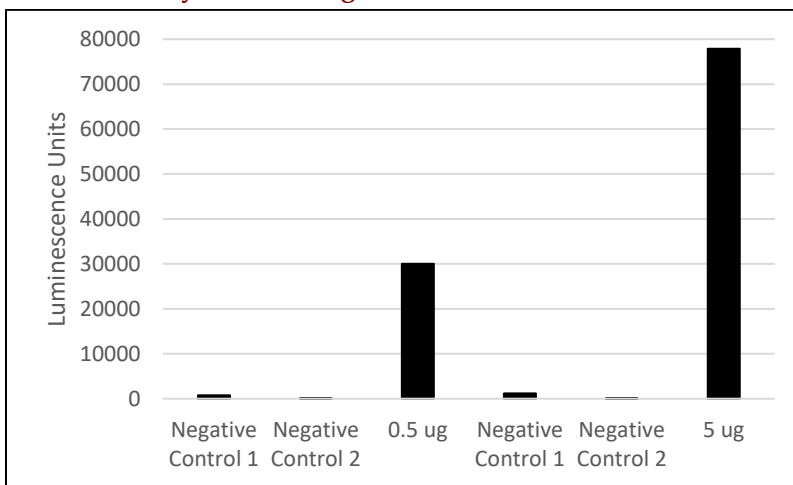
Applications: WB

Recommended Antibody Dilutions:

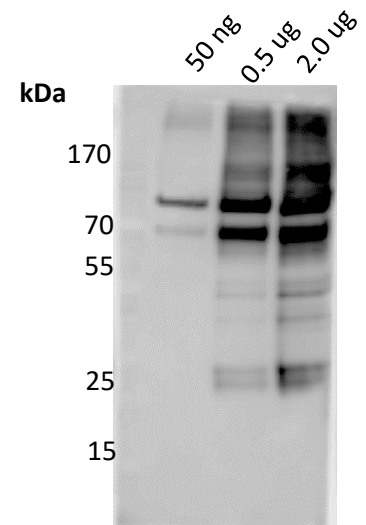
Western Blotting: Robust detection of 50ng of recombinant protein was possible when antibody was used at a final concentration of 8.43 μ g/mL

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

Detection of USP4 by AB511 using ELISA and Western Blot



0.5 or 5 ug of USP4 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. USP4 was detected by 1.69 μ g/mL of AB511 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000).
 Negative Control 1: No AB511 Negative Control 2: no detection reagent.



Indicated amounts of USP4 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 8.43 μ g/mL of AB511.
 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) Blanchette, P., et al., Association of UNP, a ubiquitin-specific protease, with the pocket proteins pRb, p107 and p130. *Oncogene*, 2001. 20(39): p. 5533-7.
- 2) Layfield, R., et al., Chemically synthesized ubiquitin extension proteins detect distinct catalytic capacities of deubiquitinating enzymes. *Anal Biochem*, 1999. 274(1): p. 40-9.
- 3) Gupta, K., M. Chevrette, and D.A. Gray, The Unp proto-oncogene encodes a nuclear protein. *Oncogene*, 1994. 9(6): p. 1729-31.

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