

USP8 Antibody (Ubiquitin Specific Protease 8)

Cat. # AB504

Background

USP8 is associated with endosomal trafficking, it is not clear that this isopeptidase is involved in the commitment to lysosomal sorting¹. Additional research has shown that USP8 plays a critical role in the stability of receptor tyrosine kinases². USP8 has been demonstrated to regulate the stability of Nrdp1, a E3 ubiquitin ligase³.

Target Alternate Names:	Deubiquitinating enzyme 8, hUBPy, HumORF8, KIAA0055, MGC129718, Ubiquitin carboxyl-terminal hydrolase 8, Ubiquitin-specific processing protease 8, Ubiquitin thioesterase 8, UBP
Target Molecular Weight:	127kDa (With tags)

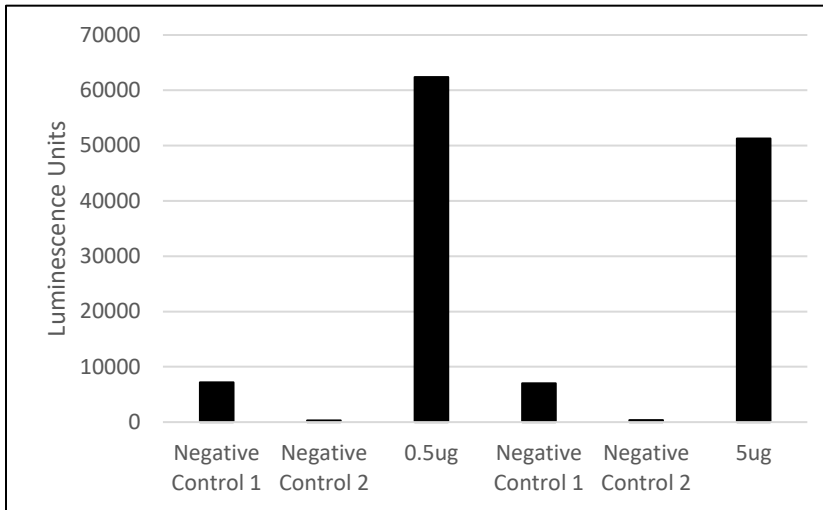
Product Information

Description:	Chicken, polyclonal antibody to USP8
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 8.28µg/mL
Storage:	Store at -20°C; Avoid repeated freeze thaw. Supplied in phosphate buffered saline pH 7.4.

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

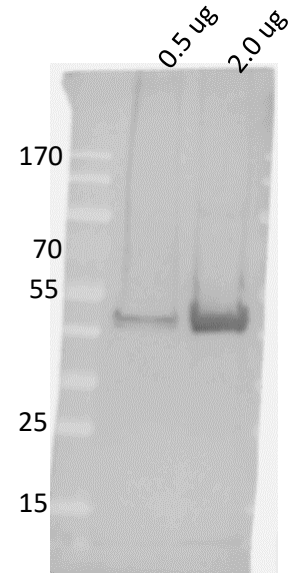
Detection of USP8 by AB504 using ELISA (A) and Western Blot (B)

(A)



0.5 or 5 ug of USP8 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. USP8 was detected by 8.28 ug/mL of AB504 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000). Negative Control 1: No USP8 antigen; Negative Control 2: no detection reagent. 2^o Antibody: α -Chicken HRP (1:5000).

(B)



Indicated amounts of USP8 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 8.28 ug/mL of AB504.

References

- 1) Niendort, S. *Essential role of ubiquitin-specific 8 for receptor tyrosine kinase stability and endocytic trafficking in vivo*. Mol Cell Biol. 2007. 27(13): p5029-39.
- 2) McCullough, J., M.J. Clague, and S. Urbe, *AMSH is an endosome-associated ubiquitin isopeptidase*. J Cell Biol, 2004. 166(4): p. 487-92.
- 3) Wu, X. et al. *Stabilization of the E3 ubiquitin ligase Nrdp1 by the deubiquitinating enzyme USP8*. Mol Cell Biol. 2004. 24(17):7748-57.

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