

AMSH Antibody (Associated molecule with the SH3 domain of STAM)

Cat. # AB301

Background

AMSH (Associated Molecule with the SH3-domain of STAM) is a JAMM domain-containing protein that regulates receptor endosomal sorting of the epidermal growth factor receptor (EGFR)¹. Recombinant AMSH functions as a deubiquitylase in vitro and ablation of AMSH activity by incubation of cells with AMSH siRNA enhances the degradation of EGFR¹. AMSH mediated cleavage of K63 linked ubiquitin chains is enhanced in the presence of its binding partner STAM².

Target Alternate Names: STAMBP, MGC126516, MGC126518, STAM-binding protein

Target Molecular Weight: 49kDa

Product Information

Description: Chicken, polyclonal antibody to AMSH

Species Cross Reactivity: Human

Source: Chicken

Applications: WB

Recommended Antibody Dilutions:

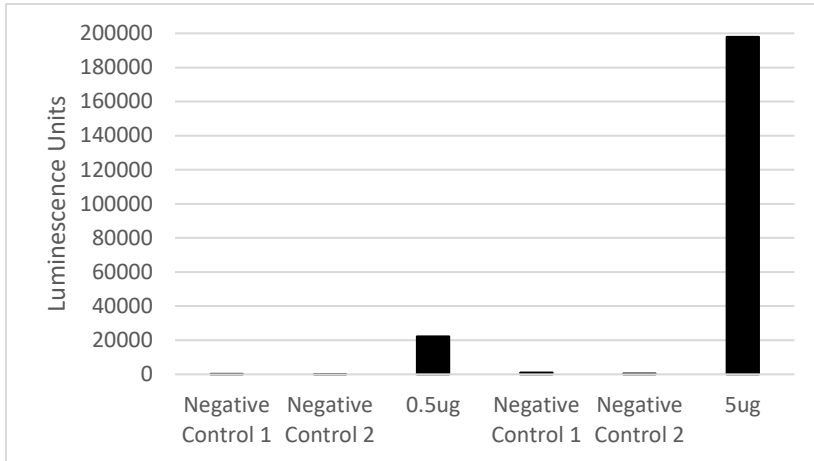
Western Blotting: Robust detection of 500ng of recombinant protein was possible when antibody was used at a final concentration of 6.88µg/mL

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

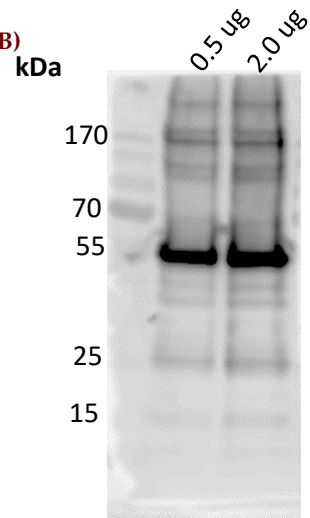
Detection of AMSH by AB301 using ELISA (A) and Western Blot (B)

(A)



0.5 or 5 ug of AMSH was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. AMSH was detected by 6.88 ug/mL of AB301 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000). Negative Control 1: No AB301; Negative Control 2: no detection reagent. 2^o Antibody: α -Chicken HRP (1:5000).

(B)



Indicated amounts of AMSH were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 6.88 ug/mL of AB304.

References

- 1) McCullough, J., et al., *Activation of the endosome-associated ubiquitin isopeptidase AMSH by STAM, a component of the multivesicular body-sorting machinery*. *Curr Biol*, 2006. 16(2): p. 160-5.
- 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.

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