

AMSH (Associated molecule with the SH3 domain of STAM)

Cat. # DB301

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.

Alternate names: STAMBP, MGC126516, MGC126518, STAM-binding protein

Product Information

Molecular Weight:	49kDa
Quantity:	25µg
Physical State:	Liquid
Source:	Human Recombinant
Tag:	His6
Activity:	This enzyme cleaves K63-linked ubiquitin chains <i>in vitro</i> .
Storage:	-80° C. Avoid repeated freeze/thaw cycles

References

- 1) McCullough, J., M.J. Clague, and S. Urbe, *AMSH is an endosome-associated ubiquitin isopeptidase*. J Cell Biol, 2004. **166**(4): p. 487-92.
- 2) 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.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved