

AMSH (Associated molecule with the SH3 domain of STAM)

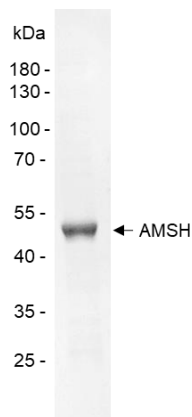
Cat. # DB301

Background	AMSH (Associated Molecule with the SH3-domain of STAM, STAM Binding Protein) is a JAMM domain-containing protein that functions as a deubiquitinase <i>in vitro</i> . AMSH mediates the cleavage of K63-linked ubiquitin chains. This cleavage is enhanced in the presence of its binding partner STAM. AMSH plays a role in the regulation of the endosomal sorting of the epidermal growth factor receptor (EGFR).
Alternate names	STAMBP, MGC126516, MGC126518, STAM-binding protein

Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	49 kDa
Quantity	25 µg
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

Product QC



SDS-Page Analysis of purified AMSH. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.

References

- 1) McCullough, J., et al., J Cell Biol, 2004. 166(4):487-92.
- 2) Tian, S., et al., Autophagy, 2021. 17(6):1367-1378.

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