

## AMSH (Associated molecule with the SH3 domain of STAM) Catalytic domain

Cat. # DB302

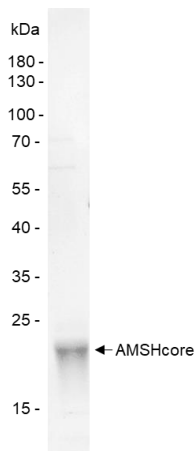
**Background** AMSH (Associated Molecule with the SH3-domain of STAM) is a JAMM domain-containing protein that regulates receptor endosomal sorting of the epidermal growth factor receptor (EGFR)<sup>1</sup>. Recombinant AMSH functions as a deubiquitylase *in vitro* and ablation of AMSH activity by incubation of cells with AMSH siRNA enhances the degradation of EGFR<sup>1</sup>. AMSH mediated cleavage of K63 linked ubiquitin chains is enhanced in the presence of its binding partner STAM<sup>2</sup>.

**Alternate names** AMSHcore

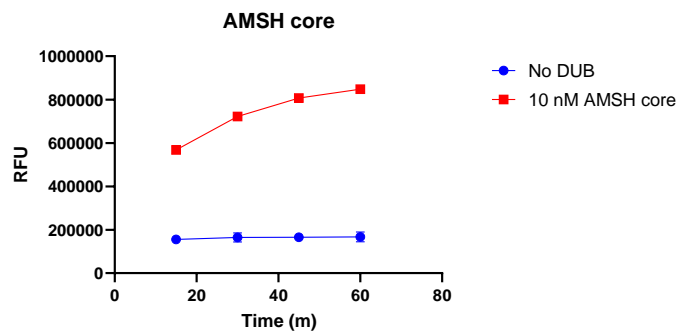
### Product Information

<b>Molecular Weight</b>	20.5 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid
<b>Source</b>	Human, recombinant, residues 250-424
<b>Tag</b>	His <sub>6</sub>
<b>Activity</b>	This enzyme cleaves K63-linked ubiquitin chains <i>in vitro</i> .
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



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



**Activity Assay of AMSH core.** 10 nM AMSHcore was tested in a Ub-Rh110 assay showing robust DUB activity.

### References

- 1) Davies, C., et al., J Mol Biol., 2011. 413(2):416-29.
- 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