



6405 Mira Mesa Blvd, Ste 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

## ***AMSH, His-tag Recombinant***

Catalog #: 81080  
Lot#: 230329

**Description:** Human AMSH recombinant protein, full length, with His-tag, expressed in E. coli.

**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). 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. AMSH mediated cleavage of K63 linked ubiquitin chains is enhanced in the presence of its binding partner STAM.

**Amt:** 25 µg

**Concentration:** 1.0 mg/ml

**MW:** 49 kDa

**Storage:** Stable for ≥1 year at –80°C.

**Applications:** This enzyme cleaves K63-linked ubiquitin chains in vitro. Suggested Substrate: DiUbiquitin K63

### **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.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)