

USP8core (Ubiquitin Specific Protease 8 core domain)

Cat. # DB503

Background: USP8 is associated with endosomal trafficking, it is not clear that this isopeptidase is involved in the commitment to lysosomal sorting¹. Additional research has shown that USP8 plays a critical role in the stability of receptor tyrosine kinases². USP8 has been demonstrated to regulate the stability of Nrdp1, a E3 ubiquitin ligase³.

Alternate names: Deubiquitinating enzyme 8, hUBPy, HumORF8, KIAA0055, MGC129718, Ubiquitin carboxyl-terminal hydrolase 8, Ubiquitin-specific processing protease 8, Ubiquitin thioesterase 8, UBPY

Product Information

Molecular Weight:	46kDa (With tags)
Quantity:	25µg
Physical State:	Liquid
Source:	Human Recombinant
Tag:	His6
Activity:	This enzyme is active in the Ub-CHOP assay.
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) Niendorf, S. *Essential role of ubiquitin-specific protease 8 for receptor tyrosine kinase stability and endocytic trafficking in vivo*. Mol Cell Biol. 2007. 27(13): p5029-39.
- 3) Wu, X. et al. *Stabilization of the E3 ubiquitin ligase Nrdp1 by the deubiquitinating enzyme USP8*. Mol Cell Biol. 2004. 24(17):7748-57.

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