

## USP7 (Ubiquitin Specific Protease 7)

Cat. # DB502

- 
- Background:** USP7 (HAUSP) was originally identified by its binding activity to a herpes viral protein, ICP0<sup>1</sup>. USP7 plays a key role in regulating the ubiquitylation of the RING-finger E3 ligase Mdm2 (and its human homolog Hdm2) by deubiquitylating and stabilizing Mdm2<sup>2,3</sup>. Additional targets of USP7 have been identified including the forkhead transcription factor, FOXO4<sup>4</sup>. Full length USP7 cleaves Ub-PLA<sub>2</sub> with high efficiency<sup>5</sup>.
- Alternate names:** Deubiquitinating enzyme 7, HAUSP, Herpesvirus-associated ubiquitin-specific protease, TEF1, Ubiquitin carboxyl-terminal hydrolase 7, Ubiquitin-specific-processing protease 7, Ubiquitin thioesterase 7
- 

### Product Information

<b>Molecular Weight:</b>	135kDa
<b>Quantity:</b>	25µg
<b>Physical State:</b>	Liquid
<b>Source:</b>	Human Recombinant
<b>Tag:</b>	His6
<b>Activity:</b>	This enzyme is active in the Ub-CHOP assay.
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

---

### References

- 1) Meredith, M., A. Orr, and R. Everett, *Herpes simplex virus type 1 immediate-early protein Vmw110 binds strongly and specifically to a 135-kDa cellular protein*. Virology, 1994. **200**(2): p. 457-69.
- 2) Li, M., et al., *A dynamic role of HAUSP in the p53-Mdm2 pathway*. Mol Cell, 2004. **13**(6): p. 879-86.
- 3) Cummins, J.M., et al., *Tumour suppression: disruption of HAUSP gene stabilizes p53*. Nature, 2004. **428**(6982): p. 1 p following 486.
- 4) van der Horst, A., et al., *FOXO4 transcriptional activity is regulated by monoubiquitination and USP7/HAUSP*. Nat Cell Biol, 2006. **8**(10): p. 1064-73.
- 5) Nicholson, B., et al., *Characterization of ubiquitin and ubiquitin-like-protein isopeptidase activities*. Protein Sci, 2008.

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