

## USP28 (Ubiquitin Specific Protease 28)

Cat. # DB525

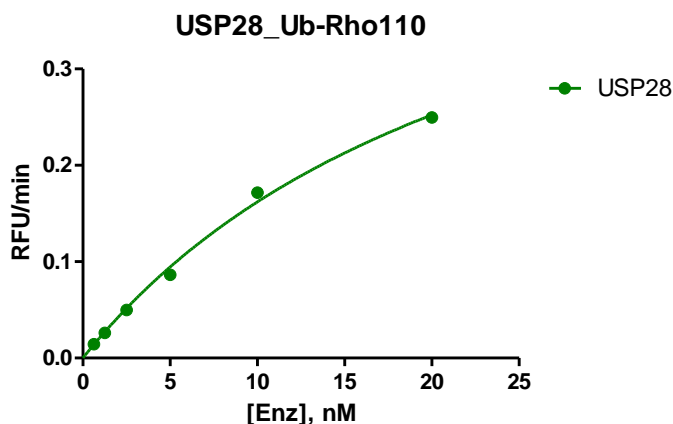
**Background:** USP28 has been shown to be required for myc stability in tumor cells<sup>1</sup>. USP28 is also highly expressed in several cancers, including colon and breast<sup>2</sup>.

**Alternate names:** Deubiquitinating enzyme 28, KIAA1515, Ubiquitin carboxyl-terminal hydrolase 28, Ubiquitin-specific-processing protease 28, Ubiquitin thioesterase 28

### Product Information

<b>Molecular Weight:</b>	123kDa
<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 with ubiquitin rhodamine 110 (Cat. # SI230) and in the Ub-CHOP assay (Cat. # PR1001).
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

### Sample Data



Initial rates of increasing concentrations of USP28 activity were assessed with ubiquitin-rhodamine 110 (Cat. # SI230) at 100nM. Assays were performed in 20mM Tris-HCl, pH 7.5, 150mM NaCl, 2mM DTT, 0.02% Tween20 at 30°C.

### References

- 1) Popov, N., et al., *The ubiquitin-specific protease USP28 is required for MYC stability*. Nat Cell Bio., 2007 **9**(7): p. 765-74.
- 2) Popov, N., et al., *Fbw7 and Usp28 regulate myc protein stability in response to DNA damage*. Cell Cycle, 2007. **6**(19): p. 2327-31.

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