

## USP51core (Ubiquitin Specific Protease 51 catalytic domain)

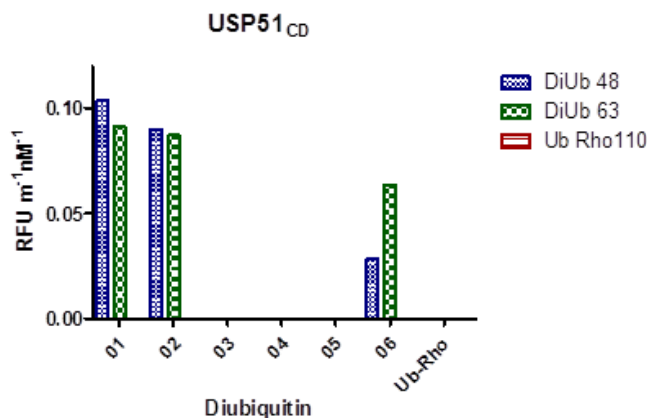
Cat. # DB507

**Background:** USP51, a member of the USP family of cysteine proteases, contains a UBP-type Zn-finger and is expressed in the prostate, brain lung, aorta, and kidney [1].

**Alternate names:** Deubiquitinating enzyme 51, Ubiquitin carboxyl-terminal hydrolase 51, Ubiquitin-specific-processing protease 51, Ubiquitin thioesterase 51

### Product Information

<b>Accession No.</b>	Q70EK9
<b>Molecular Weight:</b>	56 kDa (With tags)
<b>Quantity:</b>	25 µg
<b>Physical State:</b>	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
<b>Source:</b>	Human Recombinant
<b>Tag:</b>	His <sub>6</sub> -SUMO
<b>Suggested Substrate:</b>	UB-CHOP reporter
<b>Storage:</b>	-80°C. Avoid repeated freeze/thaw cycles



Assay Conditions: 50nM USP51, 200nM substrate, 50mM HEPEs pH 7.5, 150mM NaCl, 5mM DTT, 0.05% CHAPs

### References

1. Quesada, V., Diaz-Perales, A., Gutierrez-Fernandez, A., Garabaya, C., Cal, S., and Lopez-Otin, C., Cloning and enzymatic analysis of 22 novel human ubiquitin-specific proteases. *Biochem Biophys Res Commun.* **314**:54-62 (2004).

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