

## UHL5 (Ubiquitin C-terminal hydrolase L5)

Cat. # DB102

---

**Background:** UHL5 (UCH37) is the only ubiquitin carboxy-terminal hydrolase (UCH)-family protease, that is associated with mammalian proteasomes. It is responsible for the ubiquitin isopeptidase activity in the PA700 (19S) proteasome regulatory complex<sup>1</sup>.

**Alternate names:** AD-019, CGI-70, Ubiquitin carboxyl-terminal hydrolase isozyme L5, Ubiquitin C-terminal hydrolase UCH37, Ubiquitin thioesterase L5, UCH37, UCH-L5

---

### Product Information

<b>Molecular Weight:</b>	49kDa( with tags)
<b>Quantity:</b>	25µg
<b>Physical State:</b>	50mM Tris, pH8.0, 150mM NaCl, 10% glycerol
<b>Source:</b>	Human Recombinant
<b>Tag:</b>	His6-SUMO
<b>Activity:</b>	This enzyme is active in the Ub-CHOP assay or with ubiquitin-aml (Cat. # SI260)
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles

---

### References

- 1) Yao, T. et al. *Proteasome recruitment and activation of the Uch37 deubiquitinating enzyme by Adrm1*. Nature Cell Biology, 2006. 8: p.994 – 1002.

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