

## Yod1 (YOD1 OUT deubiquitinating enzyme 1 homolog)

Cat. # DB202

---

**Background:** Yod1 is the mammalian homolog of the *S. cerevisiae* protein Otu1. Otu1 has been demonstrated to bind Cdc48p and deubiquitylate poly-ubiquitin chains<sup>1</sup>.

**Alternate Names:** DKFZp451J1719, OTUD2, PRO0907

---

### Product Information:

<b>Molecular Weight:</b>	38kDa
<b>Quantity:</b>	25µg
<b>Physical State:</b>	Liquid
<b>Purity:</b>	>90%
<b>Source:</b>	Human Recombinant
<b>Tag:</b>	His6
<b>Activity:</b>	This enzyme cleaves Ub-AMC
<b>Storage:</b>	Store at -80°C, avoiding repeated freeze-thaw cycles.

---

### References

- 1) Rumpf S and Jentsch S. *Functional division of substrate processing cofactors of the ubiquitin-selective Cdc48 chaperone*. Mol Cell. 2006. **21**(2): p261-9.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2007 LifeSensors, Inc. All Rights Reserved