

Otub1 (Otubain 1)

Cat. # DB201

Background: Otubain 1 is a member of the OTU class of ubiquitin isopeptidases. It was originally identified as a protein capable of binding ubiquitin aldehyde¹. Researchers have shown that it is able to cleave K48-linked ubiquitin chains *in vitro*¹. Otubain 1 has been shown to be involved in T cell function and is a substrate for *Yersinia* protein kinase^{2,3}.

Alternate Names: Deubiquitinating enzyme OTUB1, FLJ20113, FLJ40710, HSPC263, MGC111158, OTB1, OTU1, Otubain 1, OTU domain-containing ubiquitin aldehyde-binding protein 1, Ubiquitin-specific-processing protease OTUB1, Ubiquitin thioesterase protein OTUB1

Product Information:

Molecular Weight:	32kDa
Quantity:	25µg
Physical State:	Liquid
Purity:	>95%
Source:	Human Recombinant
Tag:	His6
Activity:	This enzyme cleaves K48- and K63- linked ubiquitin chains <i>in vitro</i>
Storage:	Store at -80°C, avoiding repeated freeze-thaw cycles.

References

- 1) Juris, SJ et al. *Identification of otubain 1 as a novel substrate for the Yersinia protein kinase using chemical genetics and mass spectrometry*. FEBS Lett. 2006. **580**(1): p179-83.
- 2) Soares, L et al. *Two isoforms of otubain 1 regulate T cell anergy via GRAIL*. Nat Immunol. 2004. **5**(1): p45-54.
- 3) Balakirev, MY et al. *Otubains: a new family of cysteine proteases in the ubiquitin pathway*. EMBO Rep. 2003. **4**(5): p517-22.

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