

# Otub1 Antibody (Otubain 1)

Cat. # AB201

## Background

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

**Target 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.

**Target Molecular Weight:** 32kDa

## Product Information

**Description:** Chicken, polyclonal antibody to Otub1

**Species Cross Reactivity:** Human

**Source:** Chicken

**Applications:** WB

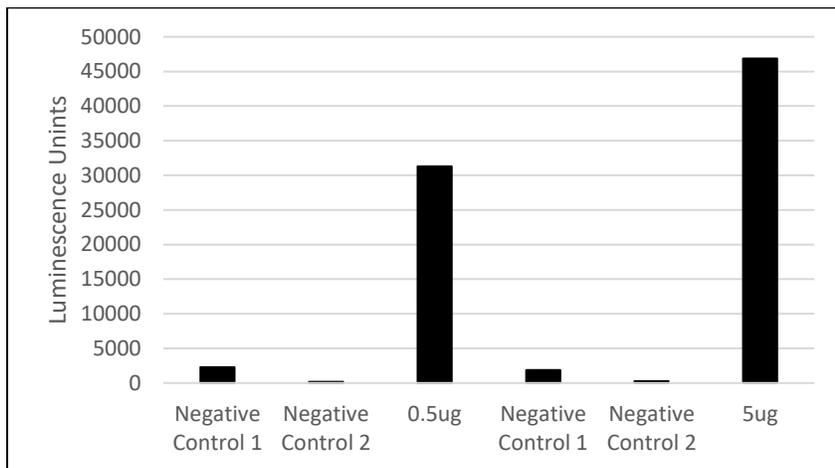
**Recommended Antibody Dilutions:**

Western Blotting: Robust detection of 50ng of recombinant protein was possible when antibody was used at a final concentration of 4.57µg/mL

**Storage:** Store at -20°C; Avoid repeated freeze thaw. Supplied in phosphate buffered saline pH 7.4.

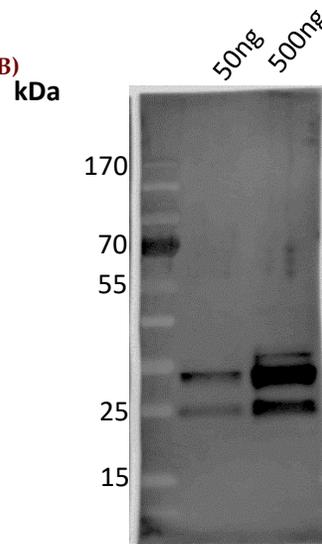
## Detection of Otub1 by AB201 using ELISA (A) and Western Blot (B)

(A)



0.5 or 5 ug of Otub1 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. Otub1 was detected by 4.57 µg/mL of AB201 using traditional ELISA detection reagent. 2<sup>o</sup> Antibody: α-Chicken HRP (1:5000). Negative Control 1: No AB201; Negative Control 2: no detection reagent. 2<sup>o</sup> Antibody: α-Chicken HRP (1:5000).

(B)



Indicated amounts of Otub1 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 4.57 µg/mL of AB201.

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

## 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