

Otub2 Antibody (Otubain 2)

Cat. # AB203

Background

Otubain 1 and 2 were originally isolated from HeLa cells by affinity purification with Ubalddehyde¹. Crystal structure determination for Otub2 revealed a unique organization of the enzymes active site². Recently, Otubains 1 and 2 were shown to negatively regulate virusinduced IFN induction and the antiviral response via deubiquitylation of TRAF3 and -6³.

Alternate Names: C14orf137, Deubiquitinating enzyme OTUB2, FLJ21916, MGC3102, OTB2, OTU2, Otubain 2, OTU domain-containing ubiquitin aldehyde-binding protein 2, Ubiquitin-specific-processing protease OTUB2, Ubiquitin thioesterase protein OTUB2

Molecular Weight: 22kDa

Specificity/Applications

Species Cross Reactivity: Human

Source: Chicken

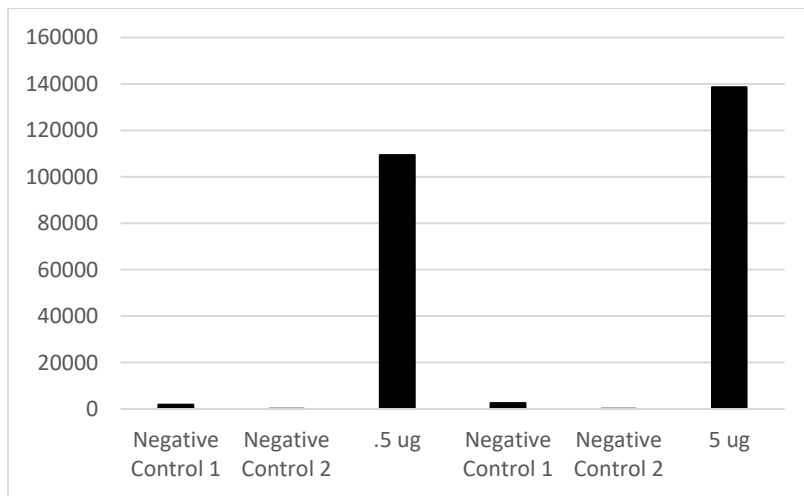
Applications: WB

Recommended Antibody Dilutions:

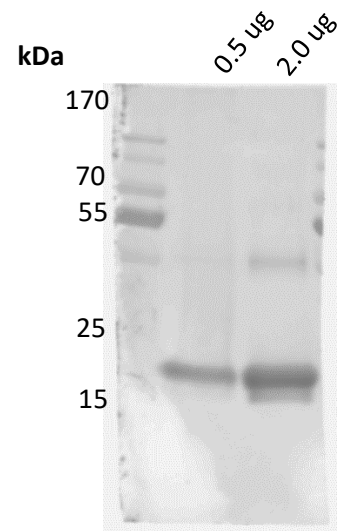
Western Blotting: Robust detection of 0.5ug of recombinant protein was possible when antibody was used at a final concentration of 9.71 µg/mL

Storage: Supplied in phosphate buffered saline. Store at -20°C or below. Avoid repeated freeze/thaw cycles.

Detection of Otub2 by AB203 using ELISA and Western Blot



0.5 or 5 ug of Otub2 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. Otub2 was detected by 1.94 µg/mL of AB203 using traditional ELISA detection reagent. 2^o Antibody: α-Chicken HRP (1:5000).
 Negative Control 1: No AB203 Negative Control 2: no detection reagent.



Indicated amounts of Otub2 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 9.71 µg/mL of AB203. 2^o Antibody: α-Chicken HRP (1:5000).

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. Balakirev, M.Y., Tcherniuk, S.O., Jaquinod, M., and Chroboczek, J., *Otubains: a new family of cysteine proteases in the ubiquitin pathway. EMBO Rep.* 4, 517-22 (2003).
2. Nanao, M.H., Tcherniuk, S.O., Chroboczek, J., Dideberg, O., Dessen, A., and Balakirev, M.Y., *Crystal structure of human otubain 2. EMBO Rep.* 5, 783-8 (2004).
3. Li, S., Zheng, H., Mao, A.P., Zhong, B., Li, Y., Liu, Y., Gao, Y., Ran, Y., Tien, P., and Shu, H.B., *Regulation of virus-triggered signaling by OTUB1- and OTUB2-mediated deubiquitination of TRAF3 and TRAF6. J Biol Chem.* 285, 4291-7 (2010).

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