

UCHL5 Antibody (Ubiquitin C-terminal hydrolase L5)

Cat. # AB102

Background

UCHL5 (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 complex¹.

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

Target Molecular Weight: 49kDa (With tags)

Product Information

Description: Chicken, polyclonal antibody to UCHL5

Species Cross Reactivity: Human

Source: Chicken

Applications: WB

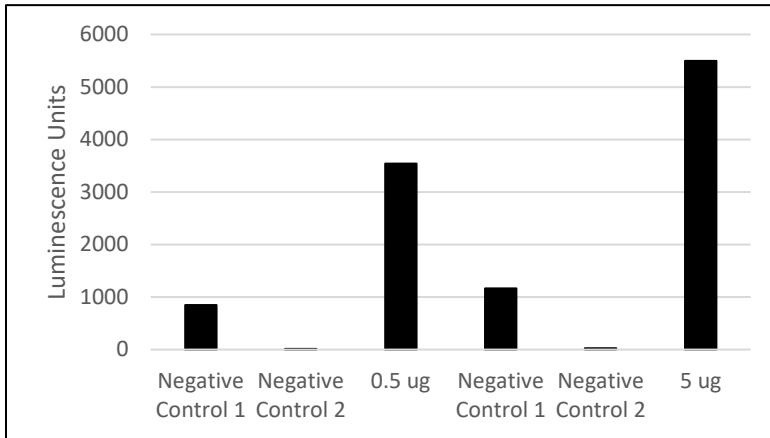
Recommended Antibody Dilutions:

Western Blotting: Robust detection of 500ng of recombinant protein was possible when antibody was used at a final concentration of 10.14µg/mL

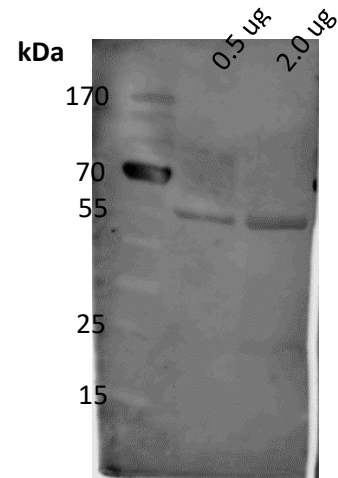
Storage: Store at -20°C; Avoid repeated freeze thaw. Supplied in phosphate buffered saline pH 7.4.

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

Detection of UCH-L5 by AB102 using ELISA and Western Blot



0.5 or 5 ug of UCH-L5 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. UCH-L5 was detected by 2.03 ug/mL of AB102 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000). Negative Control 1: No AB102 Negative Control 2: no detection reagent.



Indicated amounts of UCH-L5 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 10.14 ug/mL of AB102. 2^o Antibody: α -Chicken HRP (1:5000).

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 © 2007 LifeSensors, Inc. All Rights Reserved