

UCHL3 Antibody (Ubiquitin C-terminal hydrolase L3)

Cat. # AB101

Background

UCHL3 has

52% amino acid identity with UCHL1 and is uniformly expressed in all tissues, including brain. The activity of UCHL3 is more than 200-fold higher than UCH-L1 when a fluorogenic ubiquitin substrate is used².

Target Alternate Names: Ubiquitin carboxyl-terminal hydrolase isozyme L3, Ubiquitin thioesterase L3, UCH-L3

Target Molecular Weight: 25kDa

Product Information

Description: Chicken, polyclonal antibody to UCHL3

Species Cross Reactivity: Human

Source: Chicken

Applications: WB

Recommended Antibody Dilutions:

Western Blotting: Robust detection of 100ng of recombinant protein was possible when antibody was used at a final concentration of 5µg/mL

Storage/Purification

Polyclonal antibodies are produced by immunizing chickens with recombinant full-length protein. Antibodies are purified from egg yolks.

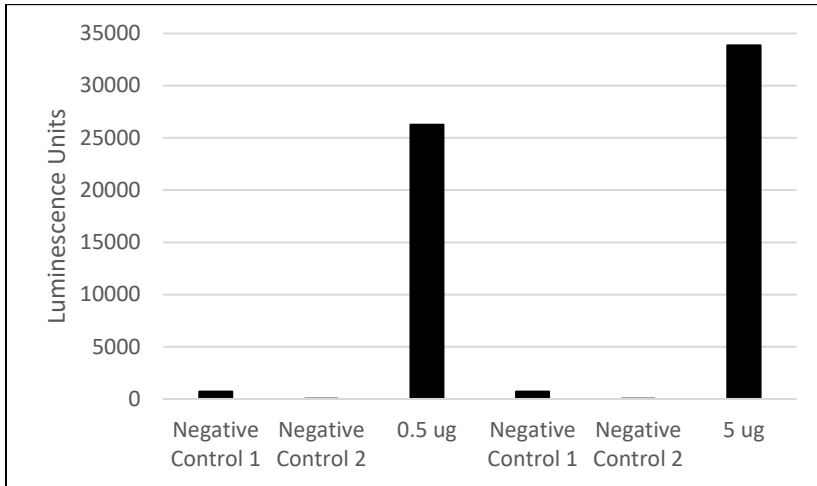
Storage: Supplied in phosphate buffered saline containing 10% glycerol.

Store at -20°C.

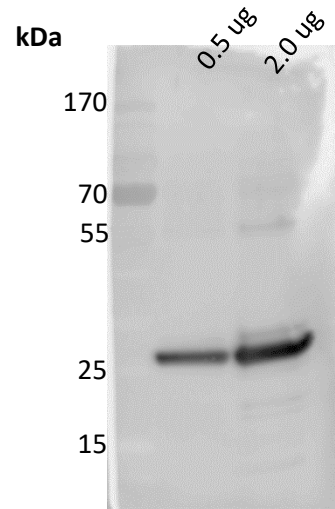
Do not aliquot the antibody.

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-L3 by AB101 using ELISA and Western Blot



0.5 or 5 ug of UCH-L3 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. UCH-L3 was detected by 0.71 ug/mL of AB101 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000). Negative Control 1: No AB101 Negative Control 2: no detection reagent.



Indicated amounts of UCH-L3 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 3.57 ug/mL of AB101. 2^o Antibody: α -Chicken HRP (1:5000).

References

- 1) Larsen C. et al. *Substrate specificity of deubiquitinating enzymes: ubiquitin C-terminal hydrolases*. Biochemistry 1998. 37: p3358-3368.
- 2) Mayer, A. N. and Wilkinson, K. D. *Detection, resolution, and nomenclature of multiple ubiquitin carboxyl-terminal esterases from bovine calf thymus*. Biochemistry 1989. 28: p166-172.

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