

UCH-L1 Antibody (Ubiquitin thioesterase L1)

Cat. # AB104

Background: UCHL1 (PGP9.5) is a 25 kDa protein; it is highly specific to neurons and to cells of the diffuse neuroendocrine system and their tumors^{1,2}. It comprises >1% of total brain protein but is almost absent from other tissues. It has been implicated both in Parkinson's disease and in lung cancer^{3,4}.

Target Alternate names: Neuron cytoplasmic protein 9.5, PARK5, PGP9.5, PGP 9.5, Ubiquitin carboxyl-terminal hydrolase isozyme L1, Ubiquitin thioesterase L1, UCH-L1

Product Information

Description: Chicken, polyclonal antibody to UCHL1

Species: Chicken

Isotype: IgY

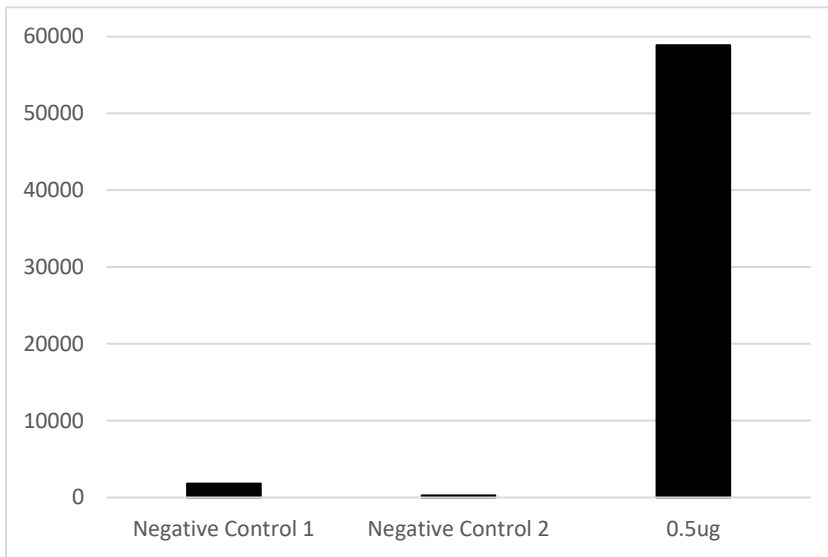
Quantity: 50; 250 µg

Application: WB, 5 µg/mL will detect 100 ng of recombinant UCHL-1

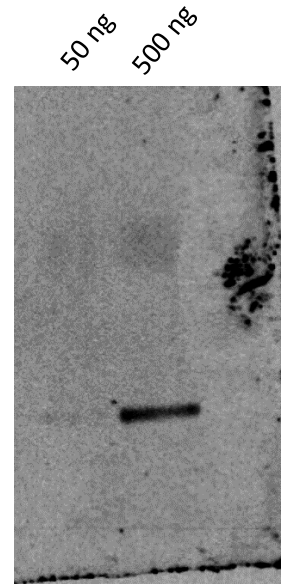
Storage: -20° C in PBS with 10% glycerol. **N.B.** aliquotting is not recommended

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved

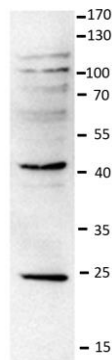
Detection of UCH-L1 by AB104 using ELISA and Western Blot



0.5ug of UCHL1 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. UCHL1 was detected by 7.86ug/mL of AB104 using traditional ELISA detection reagent. 2^o Antibody: α -Chicken HRP (1:5000).
 Negative Control 1: No AB104 Negative Control 2: no detection reagent.
 2^o Antibody: α -Chicken HRP (1:5000).



Indicated amounts of UCHL1 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 7.86 ug/mL of AB104.



Western blot analysis with crude HEK293 cell lysate

References

- 1) Liu Y, et al. The UCH-L1 gene encodes two opposing enzymatic activities that affect alphasynuclein degradation and Parkinson's disease susceptibility. *Cell* **111**: 209–218 (2002).
- 2) Larsen C. et al. Substrate specificity of deubiquitinating enzymes: ubiquitin C-terminal hydrolases. *Biochemistry* **37**:3358–3368 (1998).
- 3) Day et al. The structure of the human gene encoding protein gene product 9.5 (PGP9.5), a neuron-specific ubiquitin C-terminal hydrolase. *Biochem J.* **268(2)**: p521-4 (1990).
- 4) KD Wilkinson et al. The neuron-specific protein PGP 9.5 is a ubiquitin carboxyl-terminal hydrolase. *Science* **246**:670-673 (1989).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2009 LifeSensors, Inc. All Rights Reserved