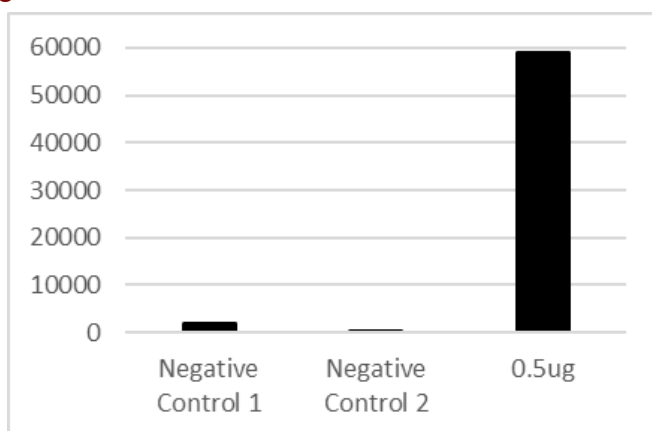


Background	UCHL1 (PGP9.5) is a 25 kDa protein specifically expressed in neurons and cells of the diffuse neuroendocrine system, as well as in their tumors. It constitutes more than 1% of the total brain protein but is almost absent from other tissues. Dysfunction of UCHL1 has been implicated in Parkinson's disease and lung cancer.
Alternate names	Neuron cytoplasmic protein 9.5, PARK5, PGP9.5, PGP 9.5, Ubiquitin carboxyl-terminal hydrolase isozyme L1, Ubiquitin thioesterase L1, UCH-L1
Application(s)	ELISA, Western Blot. Robust detection of 100 ng of recombinant protein was possible when antibody was used at a final concentration of ~5 µg/ml.

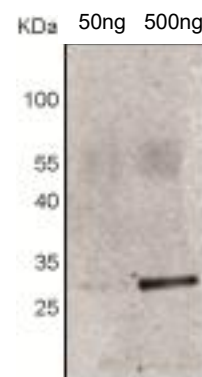
Product Specifications

Purity	>95%
Quantity	50 µg, 25 µg
Molecular Weight	25 KDa
Physical State	Supplied in phosphate buffered saline with 10% glycerol
Source & Clonality	Chicken. Polyclonal antibody to UCHL1.
Species Reactivity	Chicken
Immunogen	Full-length human UCHL1
Isotype	IgY
Stability & Storage	Over 1 year at -20° C. Avoid repeated freeze/thaw cycles. Aliquoting is not recommended

Product QC



Analysis of UCH-L1 using ELISA plate. UCH-L1 of 0.5µg was detected by 7.86 µg/mL of AB101 using traditional ELISA detection reagent. 2° Antibody: α-Chicken HRP (1:5000). Negative Control 1: No AB104; Negative Control 2: no detection reagent. AB104.



SDS-Page Analysis of purified UCH-L1. Indicated amounts of UCHL1 were loaded on SDS-PAGE gel followed by Western Blot using 7.86 µg/mL of UCH-L1. 2° Antibody: α-Chicken HRP (1:5000).

References

1. Wang X., et al. Ubiquitin C-terminal hydrolase-L1: A new cancer marker and therapeutic target with dual effects *Oncol Lett.* 25(3):123 (2023).
2. Liu Y, et al. The UCH-L1 gene encodes two opposing enzymatic activities that affect alpha-synuclein degradation and Parkinson's disease susceptibility. *Cell* 111: 209–218 (2002).
3. Larsen C. et al. Substrate specificity of deubiquitinating enzymes: ubiquitin C-terminal hydrolases. *Biochemistry* 37:3358–3368 (1998).
4. 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).
5. Wilkinson KD et al. The neuron-specific protein PGP 9.5 is a ubiquitin carboxyl-terminal hydrolase. *Science* 246:670-673 (1989).