

## UCHL1 (Ubiquitin C-terminal hydrolase L1)

Cat. # DB104

**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<sup>1,2</sup>. 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<sup>3,4</sup>.

**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

<b>Molecular Weight:</b>	25kDa
<b>Quantity:</b>	25µg
<b>Physical State:</b>	Liquid
<b>Source:</b>	Human Recombinant
<b>Tag:</b>	His6
<b>Storage:</b>	-80°C. Avoid repeated freeze/thaw cycles

### References

- 1) KD Wilkinson et al. *The neuron-specific protein PGP 9.5 is a ubiquitin carboxyl-terminal hydrolase.* Science 1989. 246: p670-673.
- 2) 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. 1990. 268(2): p521-4.
- 3) Liu Y, et al. *The UCH-L1 gene encodes two opposing enzymatic activities that affect alphasynuclein degradation and Parkinson's disease susceptibility.* Cell 2002. 111: p209–218
- 4) Larsen C. et al. *Substrate specificity of deubiquitinating enzymes: ubiquitin C-terminal hydrolases.* Biochemistry 1998. 37: p3358–3368.

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