

UCH-L1 (Ubiquitin C-terminal hydrolase L1)

Cat. # DB104

Background

UCHL1, or Ubiquitin C-Terminal Hydrolase L1, is a deubiquitinating enzyme that plays a vital role in protein homeostasis within cells. UCHL1 is highly expressed in neurons and is often used as a marker for neuronal cells. It is involved in various cellular processes, including ubiquitin recycling, proteostasis, and axonal transport. Mutations or dysregulation of UCHL1 have been linked to neurodegenerative disorders like Parkinson's and Alzheimer's diseases, highlighting its importance in maintaining neuronal health and proteostasis.

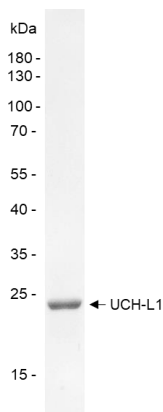
Alternate names

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

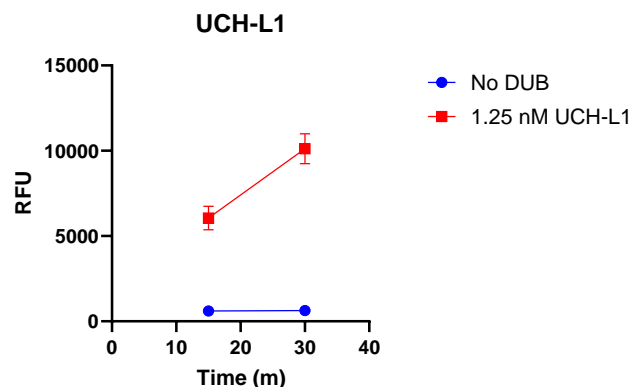
Product Information

Purity	≥ 95% by SDS-PAGE
Molecular Weight	25 kDa
Quantity	25 µg
Physical State	Liquid
Source	Human Recombinant
Tag	His6
Storage	-80°C. Avoid repeated freeze/thaw cycles

Product QC



SDS-Page Analysis of purified UCH-L1. Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



Activity Assay of UCH-L1. 1.25 nM UCH-L1 was tested in a Ub-Rh110 assay showing robust DUB activity.

References

1. Sharma, A., et al., Int J Mol Sci., 2020. 21(11):3910.
2. Wang, K., et al., Expert Opin Ther Targets, 2017. 21(6):627-638.

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