

**Background** UCHL3 (Ubiquitin C-terminal hydrolase L3) is a de-ubiquitinating enzyme and a component of the ubiquitin-proteasome system. UCHL3 has 52% amino acid identity with UCHL1 and is uniformly expressed in all tissues, including brain. UCHL3 catalyzes the hydrolysis of amides, esters and thioesters of C-terminus of ubiquitin. UCHL3 plays an important role in cell cycle regulation, DNA repair, and apoptosis via deubiquitination and deneddylation activities. UCHL3 dysregulation has been associated with neurodegenerative diseases.

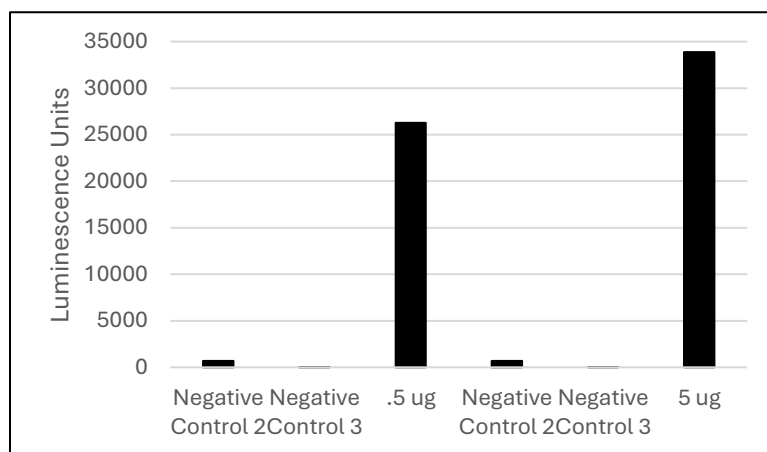
**Alternate names** Ubiquitin carboxyl-terminal hydrolase isozyme L3, Ubiquitin thioesterase L3, UCH-L3

**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.

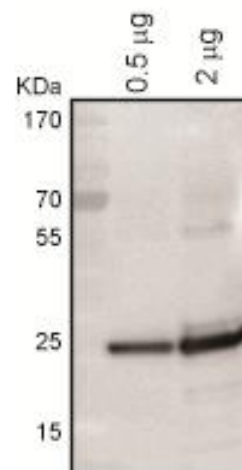
### Product Specifications

<b>Purity</b>	≥ 90% estimated by SDS-PAGE
<b>Molecular Weight</b>	25 kDa
<b>Quantity</b>	250 µg
<b>Physical State</b>	Liquid. Supplied in phosphate buffered saline pH 7.4
<b>Source &amp; Clonality</b>	Chicken. Polyclonal antibody to UCHL3.
<b>Species Reactivity</b>	Chicken, Human
<b>Immunogen</b>	Full-length human UCHL3
<b>Isotype</b>	IgY
<b>Stability &amp; Storage</b>	Over 1 year at -80° C. Avoid repeated freeze/thaw cycles

### Product QC



**Analysis of UCH-L3 using ELISA plate.** UCH-L3 of 5 µg was detected by 0.71 µg/mL of AB101 using traditional ELISA detection reagent. 2° Antibody: α-Chicken HRP (1:5000). Negative Control 1: No AB101; Negative Control



**SDS-Page Analysis of purified UCH-L3.** Indicated amounts of protein were loaded on SDS-PAGE gel followed by Western Blot using 5 µg/mL of UCH-L3. 2° Antibody: α-Chicken HRP (1:5000).

### References

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2025 LifeSensors, Inc. All Rights Reserved

1. Hirayama K., et al., Bioorg Med Chem. 2007 ;15(21) :6810-8.
2. Lei H., et al., Biochem Biophys Res Commun. 2024; 19; 734:150626.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2025 LifeSensors, Inc. All Rights Reserved