

# UCH-L1 (human), (recombinant) (His- tag)

Ubiquitin C-terminal hydrolases (UCHs) are a family of cysteine hydrolases that catalyze the hydrolysis of amides, esters and thioesters of the C-terminus of ubiquitin. Mammalian neuronal cells abundantly express a deubiquitylating enzyme, ubiquitin carboxy-terminal hydrolase 1 (UCH-L1). Mutations in UCH-L1 are linked to Parkinson's disease as well as gracile axonal dystrophy (gad) in mice. In contrast to the universally expressed UCH-L3 isozyme, UCH-L1 is expressed exclusively in neurons and testis/ovary. It has been shown that UCH-L1 associates and co-localizes with monoubiquitin and elongates ubiquitin half-life and the suggestion is made that UCH-L1, with avidity and affinity for ubiquitin, ensures ubiquitin stability within neurons.

## Ordering Information

[Order Online »](#)

BML-UW9740-0050	50µg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      Ubiquitin C-terminal hydrolase L1, PGP9.5

Formulation      Liquid. In 50 mM TRIS, pH 7.6, containing 1 mM dithiothreitol (DTT).

MW      24.8 kDa

Source      Produced in *E. coli*.

Specific Activity      >100 pmol/min/μg. Determined at 25°C with 1 μM ubiquitin-AMC (BML-SE211) as substrate and UCH-L1 (see protocol).

UniProt ID      P09936



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)