

UCH-L3 (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. UCH-L3 is a member of the lower molecular weight group of UCHs involved in the hydrolysis of small C-terminal derivatives of ubiquitin that form non-specifically during the process of protein ubiquitinylation.

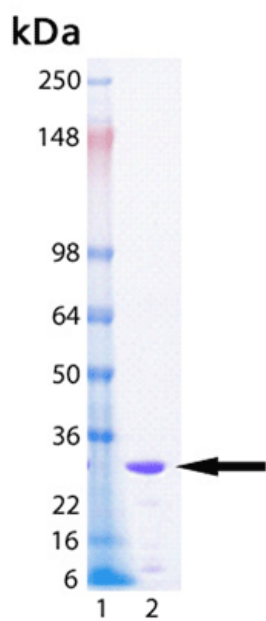
Ordering Information

[Order Online »](#)

BML-UW9745-0050	50µg
------------------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis. Lane 1: MW Marker; Lane 2: 1.0 μ g of purified UCH-L3 (human), (recombinant) (His-tag).

Handling & Storage

Use/Stability	As supplied, the enzyme is stable on ice for several hours. Activity is stable to at least 6 freeze/thaw cycles. (snap freezing in a dry/ice ethanol bath or liquid nitrogen).
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ubiquitin C-terminal hydrolase L3
Application Notes	Useful tool to study UCH-L3 kinetics and regulation, screen for potential inhibitors and activators.
Formulation	Liquid. In 50mM TRIS, pH 7.6, containing 1mM dithiothreitol (DTT).
GenBank ID	NM_006002
MW	26.2 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
Specific Activity	>1000 pmol/min/μg. Determined at 25°C with 1μM ubiquitin-AMC (Prod. No. BML-SE211) as substrate and UCH-L3 at 20pM (0.58ng/ml; see protocol).
UniProt ID	P15374

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)