

USP14 (human), (recombinant) (untagged)

USP14 is a ubiquitin-specific protease. Little is known about the biological roles of substrates of USP14, but depletion of its homolog in mice cause defects in synaptic transmission.

Citations: 1

[View Online »](#)

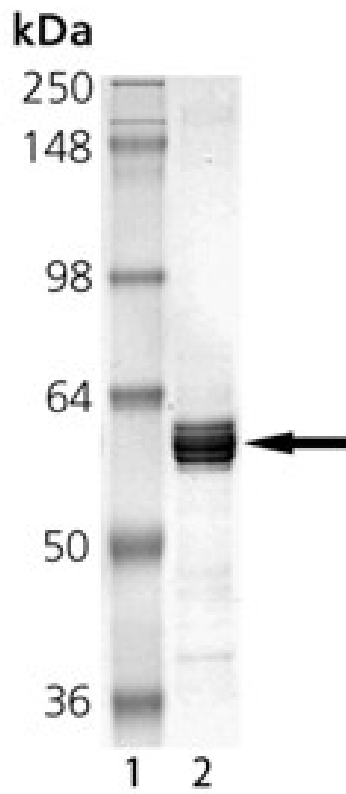
Ordering Information

[Order Online »](#)

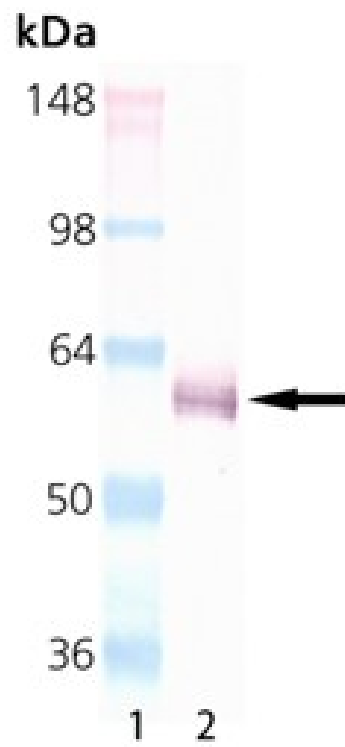
BML-UW9840-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 2.0 μ g of purified USP-14 (human), (Recombinant) protein.



Western Blot analysis: Lane 1: MW marker, Lane 2: 100ng of purified USP-14 (human), (Recombinant) protein probed with USP-14 mAb.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ubiquitin-specific protease 14
Formulation	Liquid. In 10mM TRIS containing 50mM NaCl and 1mM DTT, pH 7.4.
MW	56.1 kDa
Purity	≥90% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	P54578

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-
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)