

UbcH8 (human), (recombinant) (His- tag)

UbcH8 is a class I ubiquitin-conjugating (E2) enzyme highly similar in structure to UbcH7. UbcH8 has been demonstrated to act in conjunction with Parkin, a ring-finger-containing E3 ubiquitin protein ligase implicated in the pathogenesis of autosomal recessive familial Parkinson's Disease, and with Staring, an ring-finger E3 ligase responsible for ubiquitinylation of syntaxin 1. UbcH8 is a primary target gene of interferon- α and γ , and recent studies have indicated that UbcH8 is the E2 enzyme for the ubiquitin-like protein ISG15 (also known as UCRP). These studies suggest overlap between the ISG15 and ubiquitin conjugation pathways, and have important implications for understanding the regulation and function of ISG15 in the IFN α/β response.

Citations: 2

[View Online »](#)

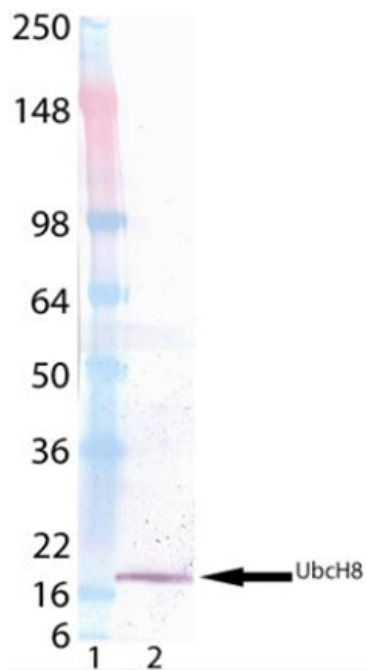
Ordering Information

[Order Online »](#)

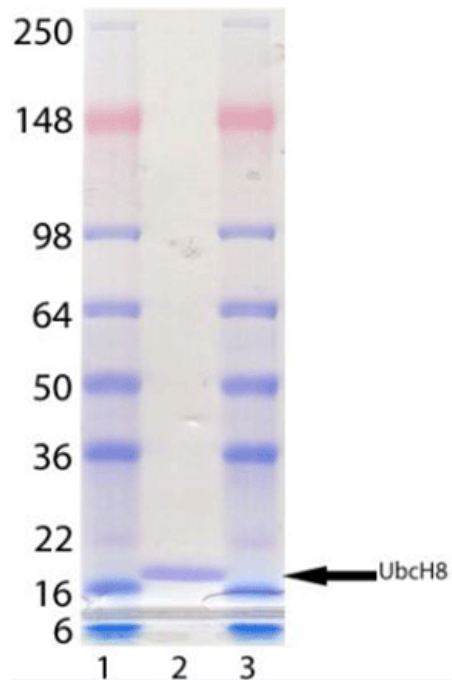
BML-UW9135-0100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of BML-UW9135: Lane 1: MW Marker, Lane 2: 100ng of UbchH8 (human), (recombinant) (His-tag) probed with Anti-6X His-tag® antibody.



SDS-PAGE analysis of Prod. No. BML-UW9135: Lane 1: MW marker, Lane 2: 2µg UbchH8 (human), (recombinant) (His-tag), Lane 3: MW marker.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Biological Activity

This enzyme supports ubiquitinylation *in vitro*.

Formulation

Liquid. In 20mM TRIS-HCl, pH 8.0, containing 0.5mM DTT.

MW

~18 kDa

Purity

≥95% (SDS-PAGE)

Source

Produced in *E. coli*.

Technical Info / Product Notes

Typical enzyme concentration to support conjugation *in vitro* is 100nM to 1µM depending upon conditions.

UniProt ID

O14933

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)