

UbcH7 (human), (recombinant) (His- tag)

Ubiquitinylation of proteins constitutes an important cellular mechanism for targeting short-lived proteins for degradation by the 26S proteasome.

Three classes of enzymes are involved in the conjugation of ubiquitin to proteins. E1, the ubiquitin activating enzyme, activates ubiquitin through the ATP-dependent formation of a high-energy thiol ester bond between the carboxyl terminus of ubiquitin and the active-site cysteine within E1. This E1-activated ubiquitin is transferred to a cysteine residue of an E2, or ubiquitin-conjugating enzyme (UbC). E2 enzymes, either by themselves or in conjunction with E3 enzymes (ubiquitin ligases), then transfer ubiquitin to target proteins forming stable isopeptide bonds resulting in multi-ubiquitin chain formation. It is the diverse combinations of E2-E3 complexes which are thought to define substrate specificity.

UbcH7 is a class I enzyme which functions in the stress response and the control of transcription factors¹. The enzyme is ubiquitously expressed with high levels of expression seen in adult muscle². It has been demonstrated to participate in the ubiquitinylation of p53, c-Fos and NF-κB. UbcH7 is one of two E2s (UbcH5 being the other) with which HECT domain proteins interact with UbcH7 being able to efficiently substitute for UbcH5 in E6-AP-dependent ubiquitinylation.

Citations: 1

[View Online »](#)

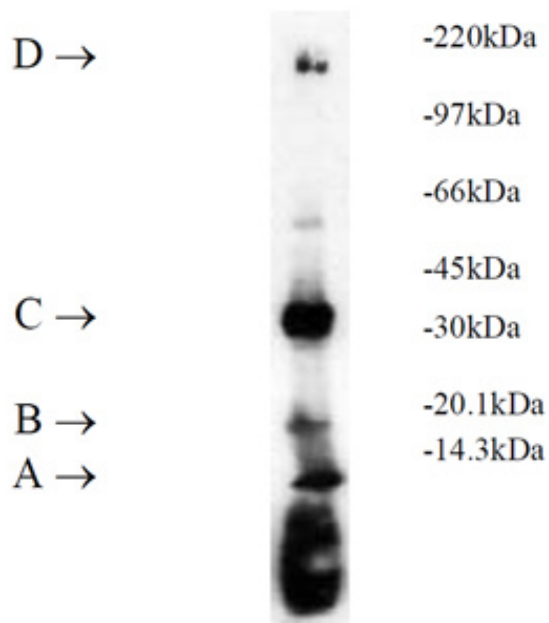
Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| BML-UW9080-0100 | 100µg |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



BML-UW9080 has been characterized in a thiolester assay using biotinylated ubiquitin as shown opposite. Labels are as follows: A – biotinylated ubiquitin; B – probably biotinylated di-ubiquitin; C – biotinyl-ubiquitin thiol ester linked Ubch7; D – biotinylubiquitin thiol ester linked E1.

Handling & Storage

Use/Stability Enzyme is stable to multiple freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

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

Biological Activity Recombinant Ubch7 charges and supports ubiquitinylation *in vitro*.

Formulation Liquid. In 50 mM HEPES, pH 7.5, containing 150 mM NaCl and 1 mM DTT.

MW ~22kDa

Purity ≥95% (SDS-PAGE)

Source Produced in *E. coli*.

UniProt ID P68036



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)