

UbcH1 (human), (recombinant) (His- tag)

Ubiquitylation 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.

Among ubiquitin-conjugating enzymes, the mammalian ubiquitin conjugating enzyme UbcH1, also known as E2-25k, is unique in its ability to catalyse the synthesis *in vitro* of unanchored Lys⁴⁸-linked poly-ubiquitin chains from mono- or poly-ubiquitin, E1, and ATP. The properties of truncated and chimeric forms of E2-25k suggest that the polyubiquitin chain synthesis activity of this E2 depends on specific interactions between its conserved 150-residue core domain and its unique 50-residue tail domain. Similarly E2-25k can catalyse the cyclisation of longer poly-ubiquitin chains, including tetra- and penta-ubiquitin. Although cyclic tri-ubiquitin resists hydrolysis by the PA700 or isopeptidase T deubiquitylating enzymes, it can be disassembled to ubiquitin monomers by isopeptidase(s) in a red blood cell extract. Thus, if cyclic poly-ubiquitin forms *in vivo*, it will not accumulate as a dead-end product.

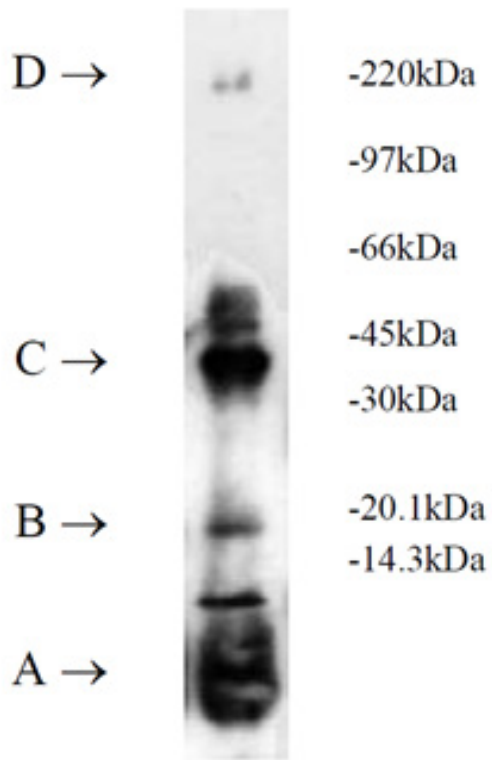
Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-UW9020-0100	100µg
-----------------	-------



BML-UW9020 has been characterised 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 Ubch1; D – biotinyl-ubiquitin 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.

Biological Activity Recombinant UbchH1 charges and supports ubiquitinylation *in vitro*.

Formulation Liquid. In 20mM TRIS-HCl, pH 7.5, containing 0.5mM DTT.

MW ~25kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified using Ni²⁺ chelation chromatography with Ni²⁺ in limiting concentration.

Source Produced in *E. coli*.

UniProt ID P61086

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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