

# UbcH13/Mms2 (human), (recombinant) (His- tag)

A heterodimer composed of the catalytically active ubiquitin-conjugating enzyme UbcH13 and its catalytically inactive paralog hMms2, which lacks an active site cysteine residue.

The UbcH13/Mms2 complex catalyzes synthesis of Lys-linked, ubiquitin chains, which serve as non-proteolytic tags involved in error-free post-replicative DNA repair and NF-κB signal transduction.

Citations: 1

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-UW9565-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Due to similarities in molecular weight (18.4kDa and 16.4kDa) it is difficult to separate the two components on SDS-PAGE under standard conditions. UW9565 has been characterised in a thiolester assay [in the presence (A) and absence (B) of ATP] using biotinylated ubiquitin (product code UW8705) as shown opposite. Major bands observed are for biotinylated ubiquitin and the biotinyl-ubiquitin thiol ester linked UbchH13 as marked. Please note that, due to the efficiency of ubiquitin transfer under the conditions used, it is not possible to observe the presence of the biotinyl-ubiquitin thiol ester linked E1.

## Handling & Storage

**Handling** Avoid 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.

**Formulation** Liquid. In 50mM HEPES, pH 8.0, 50mM sodium chloride, 10% glycerol, 1mM DTT.

**MW** ~18.4kDa

**Purity** ≥95% (SDS-PAGE)

**Source** Produced in *E. coli*.

**UniProt ID** P61088 (UbcH13), Q15819 (Mms2)

**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)