

# [C<sup>85</sup>A]Ubch5b (human), (recombinant) (His- tag)

A number of E2s in *Saccharomyces cerevisiae* and their homologues have been identified. One such family of E2s is the UBC4/5, characterised as essential for the degradation of short-lived, regulatory and abnormal proteins. Protein levels of *S. cerevisiae* UbC4/5 are up-regulated in response to stress, and their loss results in severe effects on cellular functions.

A human gene product that is 79% identical to *S. cerevisiae* UBC4/5 in amino-acid sequence was identified as Ubch5a. In addition, two other human members of this highly conserved E2 class were also cloned and designated as Ubch5b and Ubch5c, having 88% and 89% identity to Ubch5a, respectively. Members of the Ubch5a/b/c are the most active class of E2s in cell extracts. The importance of this enzyme class is underscored by the critical role of UBC4/5 in *S. cerevisiae*. Ubch5a stimulates the conjugation of ubiquitin to the tumour-repressor p53 in the presence of E6-AP and E6. Moreover, Ubch5 family is implicated in *c-fos* recognition, the modulation of which is controlled by the ubiquitin-proteasome pathway. Ubch5b and Ubch5c are associated with the signal-induced conjugation and subsequent degradation of IκBα in the presence of the SCF complexes. Ubch5c also catalyses the ubiquitination leading to the processing of p105 precursor to form p50, a subunit of the heterodimeric transcription factor NF-κB. The range and diversity of substrates and E3s with which this class of E2 enzymes interact, suggest their complex roles in cellular functions require to be studied further.

## Ordering Information

[Order Online »](#)

BML-UW9065-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## 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** Negative control for ubiquitin conjugating enzyme Ubch5b (Prod. No. BML-UW9060). The [C85A] mutation completely abolishes the ability of the enzyme to form a thiol ester.

**Biological Activity** The His<sub>6</sub>-tagged fusion proteins of Ubch5a, Ubch5b and Ubch5c all charge and support ubiquitinylation *in vitro*. Unlike their GST-tagged counterparts, the His<sub>6</sub>-tagged Ubch5 family members all appear to form thiol ester conjugates with ubiquitin at a similar rate under similar conditions. The [C<sup>85</sup>A] mutation completely abolishes the ability of the enzyme to form a thiol ester and thereby acts as an excellent negative control.

**Formulation** Liquid. In 20mM TRIS-HCl, pH7.5, containing 0.5mM DTT.

**MW** ~18kDa

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

**UniProt ID** P62837



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)