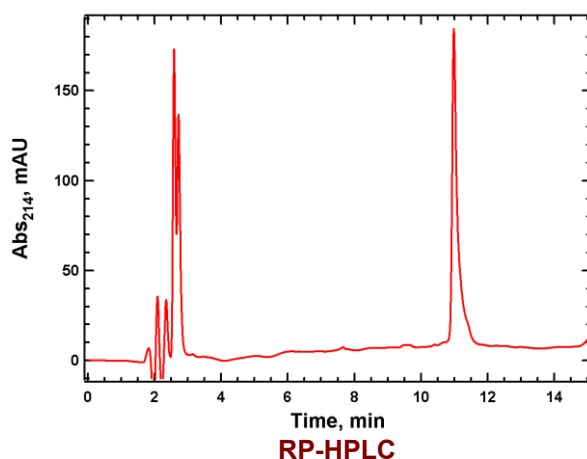


<b>Background</b>	UBE2D2 is an E2 (ubiquitin conjugating enzyme) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E3 enzymes. For the initial mono ubiquitination of substrates, UBE2D2 is involved as donor and UBE2C and UBE2S are involved in subsequent K-11 linked chain extension. UBE2D enzymes have been implicated in number of key pathways, such as inflammatory and cancer pathways, including the ubiquitylation of p53 (involving the E3 ligase Hdm2) and IκBα (involving the E3 ligase SCF). In addition, UBE2D enzymes function in transcriptional control as donors for ubiquitination of histone tail by the Polycomb protein Ring1B and DNA methylation regulator UHRF1.
<b>Alternate Names</b>	Ubch5b, E2-17K2
<b>Application(s)</b>	Ubiquitin ligation reactions

### Product Specifications

<b>Tag</b>	His <sub>6</sub>
<b>Purity</b>	> 95% by RP-HPLC
<b>Molecular Weight</b>	17,729.6 Da by MS (calculated 17,729.3)
<b>Quantity</b>	20 µL or 75 µL of a 40 µM solution (0.8 and 3 nmoles, respectively)
<b>Species</b>	Human, recombinant; Accession No. P62837
<b>Expression System</b>	<i>E. Coli</i>
<b>Physical State</b>	Liquid
<b>Buffer</b>	50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
<b>Solubility</b>	> 3 mg/mL
<b>Stability &amp; Storage</b>	Over 1 year at -80° C. Avoid repeated freeze/thaw cycles

### Product QC



### References

1. Roman-Trufero M., et al., Front Cell Dev Biol, 2022;10:1058751.
2. DaRosa, P. A., et al., Mol. Cell, 2018; 72, 753–7.
3. Yang, J., et al., Cancer Res, 2021; 81, 898–909.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2025 LifeSensors, Inc. All Rights Reserved