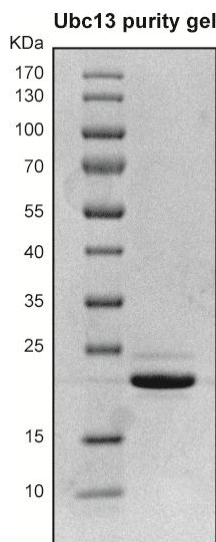


Background	Ubc13 is an E2 (ubiquitin conjugating enzyme that is involved in the conjugation of ubiquitin to target substrates along with E1 and E3 enzymes. Ubc13 acts with the UEV protein Mms2 forming a heterodimeric complex that assembles K63-linked polyubiquitin chains utilized in non-degradative ubiquitin signaling pathways (DNA repair, signal transduction). The Ubc13/Mms2 complex is unique in that it can form free K63-linked polyubiquitin chains in solution in the absence of an E3. Studies have shown Ubc13 as a potential therapeutic target for intervention in various diseases, such as different cancers, chronic inflammation, and viral infections.
Alternate Names	Ubiquitin-Conjugating Enzyme E2N, E2 Ubiquitin-Conjugating Enzyme N, UbcH-Ben, HEL-S-71
Application(s)	Ubiquitin ligation reactions

Product Specifications

Tag	His ₆
Purity	≥ 95% by SDS-PAGE
Molecular Weight	17 kDa
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	<i>S. cerevisiae</i>
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	50 mM HEPES, pH7.5, 150 mM NaCl, 10% glycerol
Activity	A typical enzyme concentration of 0.1-5 mM is used for in vitro conjugation, depending on the experimental conditions
Stability & Storage	Over 1 year at -80°C. Avoid repeated freeze/thaw cycles

Product QC



References

1. Hodge CD., et al., *Oncotarget*. 2016;7(39):64471-64504.
2. Eddins, M.J., et al., *Nat Struct Mol Biol*. 2006; **13**, 915-20.

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