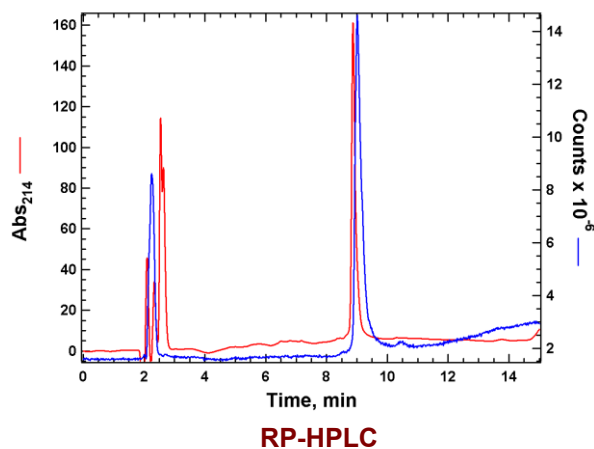


Background	UBE2Z is a ubiquitin-conjugating enzyme (E2) that appears to be a specific substrate for the Uba6 ubiquitin-activating enzyme (E1). Although the specific cellular functions of UBE2Z remain poorly characterized, it has been implicated in proapoptotic pathways. The protein exists in multiple splice variants, including a full-length isoform of 354 amino acids and a shorter isoform (isoform 2) of 181 amino acids. LifeSensors currently offers the 181-amino acid isoform.
Alternate Names	Uba6-Specific E2 Conjugating Enzyme 1, Ubiquitin-Conjugating Enzyme E2 Z, Ubiquitin Carrier Protein Z, FLJ13855, HOYS7, Use1
Application(s)	Ubiquitin ligation reactions

Product Specifications

Tag	None
Purity	≥ 95% by RP-HPLC
Molecular Weight	28,075.4 Da by MS (calculated 28,075.1)
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	Human, recombinant; Accession No. Q9H832
Expression System	<i>E. coli</i>
Physical State	Liquid.
Buffer	25 mM Tris, pH 7.4; 150 mM NaCl; 10 mM DTT; 10% glycerol
Solubility	> 3 mg/ml
Stability & Storage	1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC



References

1. Cheng, Q., et al., (2025). The emerging role and mechanism of E2/E3 hybrid enzyme UBE2O in human diseases. *Biomedicines*, 13(5), 1082.
2. Lee, D., et al., (2024). UBA6 inhibition accelerates lysosomal TRPML1 depletion and exosomal secretion in lung cancer cells. *International Journal of Molecular Sciences*, 25(5), 2843.
3. Jin, J., et al., (2007). Dual E1 activation systems for ubiquitin differentially regulate E2 enzyme charging. *Nature*, 447(7147), 1135–1138.
4. Park, K. M., et al., (2007). Identification of novel regulators of apoptosis using a high-throughput cell-based screen. *Molecular Cells*, 23(2), 170–174.

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