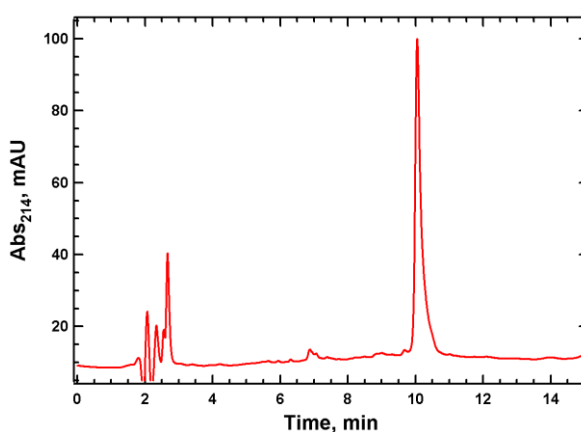


Background	UBE2G2, the human homolog of yeast Ubc7, is involved in the ubiquitination of misfolded ER proteins, targeting them for proteasomal degradation. Structural studies indicate that this single-domain protein possesses a loop region likely involved in interactions with RING domain-containing E3 ligases, which may confer binding specificity and stability. UBE2G2 is ubiquitously expressed, with particularly high levels in muscle tissues. UBE2G2 also plays a key role in sterol-regulated degradation pathways, including the ERAD-mediated turnover of HMG-CoA reductase.
Alternate Names	Ubiquitin-Conjugating Enzyme E2G2 (UBC7 Homolog, Yeast), Ubiquitin Carrier Protein G2, Ubiquitin Conjugating Enzyme 7
Application(s)	Ubiquitin ligation reactions

Product Specifications

Tag	His ₆ -SUMO
Purity	≥ 95% by RP-HPLC
Molecular Weight	30,585.1 Da by MS (calculated 30,585.8)
Quantity	20 µl or 75 µl of a 40 µM solution (0.8 and 3 nmoles, respectively)
Species	Human, recombinant; Accession No. P60604
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
Solubility	> 3 mg/mL
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC**RP-HPLC****References**

1. Hosseini SM, et al. Clin Chim Acta. 2019; 498:126-134.
2. Kikkert M, et al. J Biol Chem. 2004;279(9):3525–3534.

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