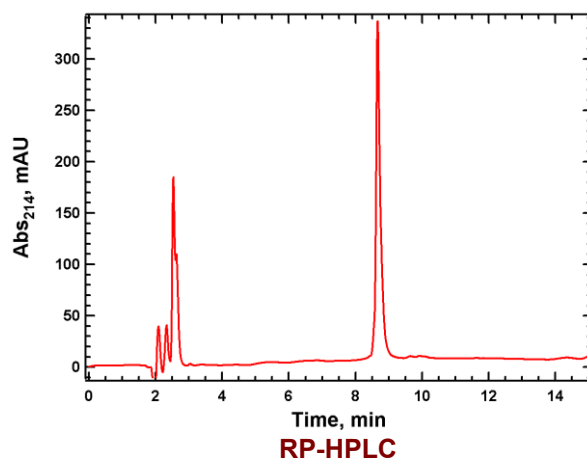


Background	UBE2S (also known as E2-24K or E2-EPF) is an E2 ubiquitin-conjugating enzyme that mediates the transfer of ubiquitin to target substrates in coordination with E1 and E3 enzymes. It has been identified as an E2 enzyme for the anaphase-promoting complex/cyclosome (APC/C), where it promotes the elongation of K11-linked polyubiquitin chains on APC/C substrates that have been initially ubiquitinated by either UBE2C or UBE2D1. This chain elongation enhances the degradation of APC/C substrates by the proteasome. Additionally, UBE2S targets von Hippel–Lindau (VHL) for proteasomal degradation, leading to the stabilization of HIF-1 α .
Alternate Names	Ubiquitin-Conjugating Enzyme E2 S, Ubiquitin-Conjugating Enzyme E2-24 KD (E2-24K), E2-EPF, E2-EPF5, EPF5
Application(s)	Ubiquitin ligation reactions

Product Specifications

Tag	None
Purity	$\geq 95\%$ by RP-HPLC
Molecular Weight	23,845.7 Da by MS (calculated 23,845.3)
Quantity	20 μ l or 75 μ l of a 40 μ M solution (0.8 and 3 nmoles, respectively)
Species	Human, recombinant; Accession No. Q16763
Expression System	<i>E. Coli</i>
Physical State	Liquid
Buffer	20 mM Tris, pH 7.4, 150 mM NaCl, 10% glycerol, 10 mM DTT
Solubility	> 3 mg/ml
Stability & Storage	1 year at -80°C. Avoid repeated freeze/thaw cycles

Product QC



References

1. Zhang M., et al., Heliyon. 2024;10(2):e24465.
2. Xiao K, Peng S, Lu J, et al. Cell Death Dis. 2023;14(7):408.
3. Wu, T. et al. Proc Natl Acad Sci U S A. 2010; 107, 1355-60.
4. Garnett, M.J. et al. Nat Cell Biol. 2009; 11, 1363-9.

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