

UBE2S

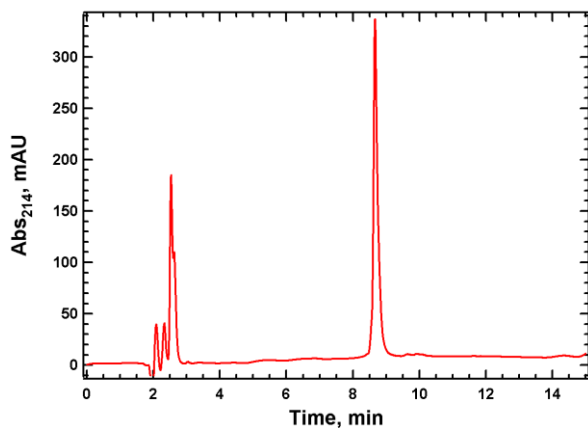
Cat. # UB221

Background: UBE2S (E2-24K, E2-EPF) is an E2 (ubiquitin conjugating enzyme) that is involved in the conjugation of ubiquitin to target substrates along with E1 and E3 enzymes. It has been identified as an E2 for the anaphase-promoting complex (APC/C) promoting the elongation of K11-linked polyubiquitin chains on APC/C substrates which have been initiated by either the E2 UBE2C or UBE2D1. This enhances degradation of APC/C substrates by the proteasome. UBE2S also targets von Hippel-Lindau (VHL) for proteasome degradation, resulting in the stabilization of HIF-1 α .

Application: Ubiquitin ligation reactions

Product Information

| | |
|--------------------------|---|
| Organism | Human, recombinant; Accession No. Q16763 |
| Purity: | $\geq 95\%$ by RP-HPLC |
| Molecular Weight: | 23,845.7 Da by MS (calculated 23,845.3) |
| Tag | none |
| Physical State: | Liquid, 20 mM Tris, pH 7.4, 150 mM NaCl, 10% glycerol, 10 mM DTT |
| Quantity: | 20 or 75 μ L of a 40 μ M solution (0.8 or 3 nmoles, respectively) |
| Solubility: | >3 mg/mL |
| Storage: | -80 $^{\circ}$ C. Avoid repeated freeze/thaw cycles |



RP-HPLC

References

1. Wu, T. et al. UBE2S drives elongation of K11-linked ubiquitin chains by the anaphase-promoting complex. Proc Natl Acad Sci U S A 107, 1355-60.
2. Williamson, A. et al. Identification of a physiological E2 module for the human anaphase-promoting complex. Proc Natl Acad Sci U S A 106, 18213-8 (2009).
3. Garnett, M.J. et al. UBE2S elongates ubiquitin chains on APC/C substrates to promote mitotic exit. Nat Cell Biol 11, 1363-9 (2009).

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved