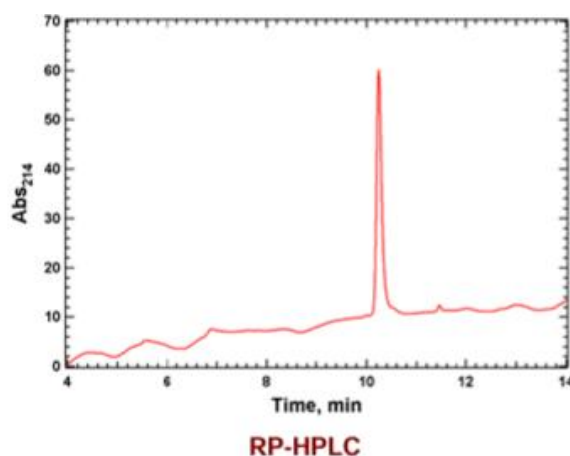


Background	UBE2D1 is a widely expressed E2 ubiquitin-conjugating enzyme that mediates the ubiquitination and subsequent degradation of abnormal or short-lived proteins by working in concert with E1 activating and various E3 ligase enzymes. It is one of four members of the UBE2D family, which includes UBE2D2, UBE2D3, and UBE2D4, and is involved in ubiquitinating diverse substrates across multiple cellular processes. UBE2D1 functions alongside several E3 ligase complexes, including those containing SIAH1, CACYBP/SIP, SKP1, APC, and TBL1X, and collaborates with the E3 ligase E6AP to mediate ubiquitination of p53/TP53. Additionally, UBE2D1 plays a role in regulating angiogenesis by controlling VEGFR2 levels, with hypoxia-responsive angiogenesis in skeletal muscle being impaired when UBE2D1 levels are elevated by TNF α stimulation. It has also been shown to interact with BRCA1/BARD1, preferentially catalyzing atypical, non-degradative K6-linked polyubiquitin chains.
Alternate Names	Ubiquitin-Conjugating Enzyme E2D 1, Ubch5, E2(17)KB1, Ubch5A, UBC4/5, Stimulator Of Fe Transport (SFT)
Application(s)	Ubiquitin ligation reactions

Product Specifications

Tag	His ₆ -SUMO
Purity	> 95% by RP-HPLC
Molecular Weight	28,261 Da (with tags)
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. P51668
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
Stability & Storage	Over 1 year at -80° C. Avoid repeated freeze/thaw cycles

Product QC



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

UBE2D1, His6-SUMO

Cat. # UB210T

References

1. Roman-Trufero M., et al., Front Cell Dev Biol, 2022;10:1058751.
2. Basic, V. T., et al., Int. J. Mol. Med.2014; 34, 228-236.
3. Nishikawa H et al., J. Biol. Chem. 2004; 279, 3916–3924.

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