

## UBE2D1, His<sub>6</sub>SUMO

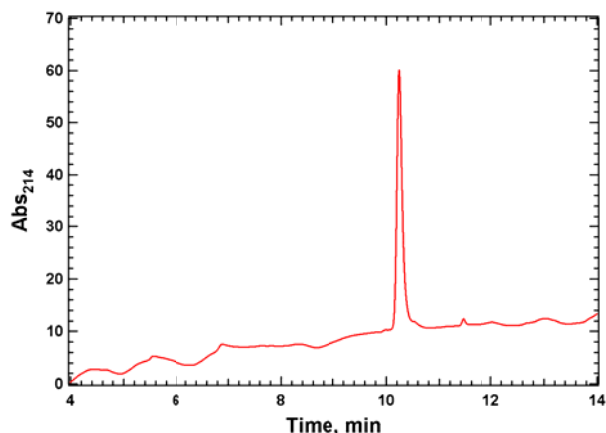
Cat. # UB210T

**Background:** UBE2D1 is a 147 amino acid (16.6kD) protein that mediates degradation of abnormal proteins or proteins with short half-lives. It is a component of a Ubiquitin ligase complex including UBE2D1, SIAH1, CACYBP/SIP, SKP1, APC and TBL1X which interacts with RNF11. It also functions in E6AP mediated ubiquitylation of p53/TP53 [1,2].

**Application:** Ubiquitin ligation reactions

### Product Information

<b>Organism</b>	Human, recombinant; Accession No. P51668
<b>Purity:</b>	≥ 95% by RP-HPLC
<b>Molecular Weight:</b>	28,621 Da
<b>Tag</b>	His <sub>6</sub> -SUMO
<b>Physical State:</b>	Liquid, 50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
<b>Quantity:</b>	20 or 75 µL of a 40 µM solution (0.8 and 3 nmoles, respectively)
<b>Solubility:</b>	>3 mg/mL
<b>Storage:</b>	-80° C. Avoid repeated freeze/thaw cycles



RP-HPLC

### References:

1. Matsuzawa, S. and Reed, J.C. Siah-1, SIP, and Ebi collaborate in a novel pathway for  $\beta$ -catenin degradation linked to p53 responses. *Mol. Cell* **7**,915-926 (2001).
2. Subramaniam, V., Li, H., Wong, M.J., Kitching, R., Attisano, L., Wrana, J., Zubovits, J., Burger, A.M., and Seth, A.K. The RIMG-H2 protein RNF11 is overexpressed in breast cancer and is a target of Smurf2 E3 ligase. *Br. J Cancer* **89**,1538-1544 (2003).

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