

UBE2L3

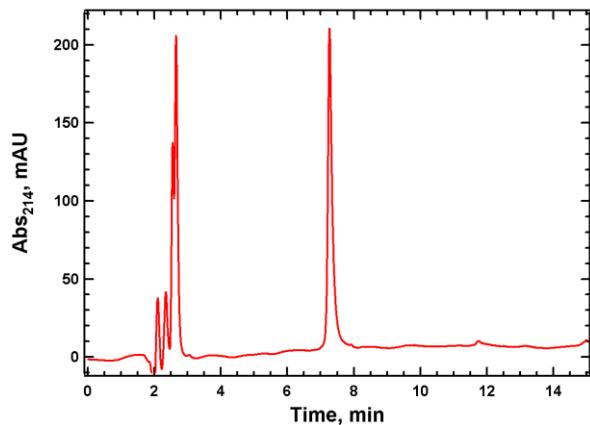
Cat. # UB202

Background: UBE2L3 (UbcH7) is an E2 (ubiquitin conjugating enzyme) ligase involved in the conjugation of ubiquitin to target substrates, along with E1 and E3 enzymes. UbcH7 has been shown to work with HECT domain E3's (E6-AP), in addition to RING containing E3's. UbcH7 is homologous to the UbcH5 family.

Application: Ubiquitin ligation reactions

Product Information

Organism	Human, recombinant; Accession No. P68036
Purity:	≥ 95% by RP-HPLC
Molecular Weight:	17861.9 Da by MS (calculated 17861.6)
Tag	none
Physical State:	Liquid, 50 mM Tris, pH 7.5; 150 mM NaCl; 10 mM DTT; 10% glycerol
Quantity:	20 µL or 75 µL of a 40 µM solution (0.8 and 3 nmoles, respectively)
Solubility:	>3 mg/mL
Storage:	-80° C. Avoid repeated freeze/thaw cycles



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