

UBE2G2

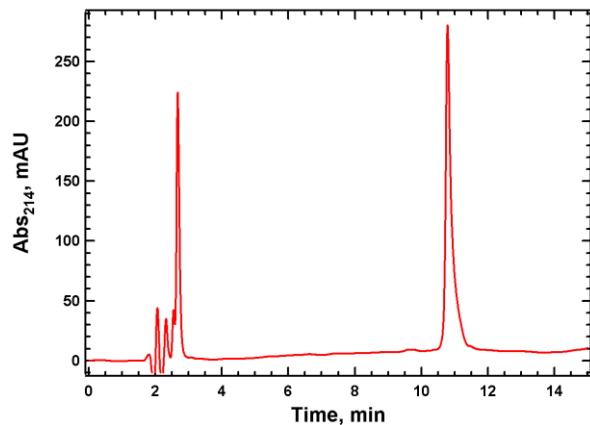
Cat. # UB227

Background: Human homolog of yeast Ubc7 involved in endoplasmic reticulum associated degradation (ERAD). Structural studies indicate that it is a single domain protein with a loop region likely involved in interaction with the RING domain to impart binding specificity or stability. It is a 165 amino acid protein expressed ubiquitously, but highly expressed in muscle.

Application: Ubiquitin ligation reactions

Product Information

Organism	Human, recombinant; Accession No. P60604
Purity:	≥ 95% by
Molecular Weight:	18565.3 Da by MS (calculated 18566.4)
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



RP-HPLC

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